

AssayMAP Bravo Platform

Labware Reference Guide



Notices

Manual Part Number

D0012628 Revision A

October 2022

Copyright

© Agilent Technologies, Inc. 2022

No part of this manual may be reproduced in any form or by any means (including electronic storage and retrieval or translation into a foreign language) without prior agreement and written consent from Agilent Technologies, Inc. as governed by United States and international copyright laws.

Contact Information

Agilent Technologies Inc.
Automation Solutions
5301 Stevens Creek Blvd.
Santa Clara, CA 95051
USA

Web:
<https://www.agilent.com>

Contact page:
<https://www.agilent.com/en/contact-us/page>

Documentation feedback:
documentation.automation@agilent.com

Acknowledgements

Microsoft® and Windows® are either registered trademarks or trademarks of the Microsoft Corporation in the United States and other countries.

Warranty

The material contained in this document is provided "as is," and is subject to being changed, without notice, in future editions. Further, to the maximum extent permitted by applicable law, Agilent disclaims all warranties, either express or implied, with regard to this manual and any information contained herein, including but not limited to the implied warranties of merchantability and fitness for a particular purpose. Agilent shall not be liable for errors or for incidental or consequential damages in connection with the furnishing, use, or performance of this document or of any information contained herein. Should Agilent and the user have a separate written agreement with warranty terms covering the material in this document that conflict with these terms, the warranty terms in the separate agreement shall control.

Technology Licenses

The hardware and/or software described in this document are furnished under a license and may be used or copied only in accordance with the terms of such license.

Restricted Rights Legend

If software is for use in the performance of a U.S. Government prime contract or subcontract, Software is delivered and licensed as "Commercial computer software" as defined in DFAR 252.227-7014 (June 1995), or as a "commercial item" as defined in FAR 2.101(a) or as "Restricted computer software" as defined in FAR 52.227-19 (June 1987) or any equivalent agency regulation or contract clause. Use, duplication or disclosure of Software is subject to Agilent Technologies' standard commercial license terms, and non-DOD Departments and Agencies of the U.S. Government will receive no greater than Restricted Rights as defined in FAR 52.227-19(c)(1-2) (June 1987). U.S. Government users will receive no greater than Limited Rights as defined in FAR 52.227-14 (June 1987) or DFAR 252.227-7015 (b)(2) (November 1995), as applicable in any technical data.

Safety Notices

WARNING

A **WARNING** notice denotes a hazard. It calls attention to an operating procedure, practice, or the like that, if not correctly performed or adhered to, could result in personal injury or death. Do not proceed beyond a **WARNING** notice until the indicated conditions are fully understood and met.

CAUTION

A **CAUTION** notice denotes a hazard. It calls attention to an operating procedure, practice, or the like that, if not correctly performed or adhered to, could result in damage to the product or loss of important data. Do not proceed beyond a **CAUTION** notice until the indicated conditions are fully understood and met.

Contents

Preface	v
About this guide	vi
Reporting problems	viii
1. Labware options	1
Labware options for each application	2
Labware options for each utility	12
2. Labware photo reference	21
12 Column, Low Profile Reservoir, Natural PP	23
8 Row, Low Profile Reservoir, Natural PP	24
Reservoir, Seahorse 201254-100, PP, no walls, pyramid bottoms	25
Reservoir, Axygen Scientific RES-SW96-LP, 86mL pyramid bottom	26
96 AbGene PCR plate	27
96 AbGene 1127, 1mL Deep Well, Square Well, Round Bottom	28
96 AbGene U-Bottom Thermal Insert + AbGene 1127, Square Well, Round Bottom	29
96 Agilent, 2mL Square Deep Well	30
96 Bio-Rad PCR, Hard-Shell, Low-Profile, Full Skirt	31
96 Red PCR Insert + Bio-Rad PCR, Hard-Shell, Low-Profile, Full Skirt	32
96 Corning 9017, Flat Bottom, Polystyrene, Clear	33
96 Costar 3363, PP Conical Bottom	34
96 Eppendorf 30129300, PCR, Full Skirt, PolyPro	35
96 Red PCR Insert + Eppendorf 30129300, PCR, Full Skirt	36
96 Eppendorf 96-1000_U-bottom, Clear PolyPro	37
96 Eppendorf 96-500_V-bottom, Clear PolyPro	38
96 Greiner 650201, U-Bottom Standard, PolyPro	39
96 Greiner 650201_U-Bottom, Clear PolyPro	40
96 Greiner U-Bottom Thermal Insert + Greiner 650201_U-Bottom, Clear PolyPro	41
96 Greiner 650207_U-Bottom, White PolyPro	42
96 Greiner 651201_V-Bottom, Clear PolyPro	43
96 Greiner V-Bottom Thermal Insert + Greiner 651201, V-Bottom, Clear PolyPro	44
96 Greiner 652270, PCR, Full Skirt, PolyPro	45
96 Red PCR Insert + Greiner 652270, PCR, Full Skirt, PolyPro	46
96 Greiner 655101 PS Clr Rnd Well Flat Btm	47
96 Greiner 675801, Half Area, Flat Bottom, UV Star	48
96 Nunc 269620, Flat Bottom, Polystyrene	49
96 Thermo Matrix 3732, V-bottom, 0.75ml Storage Tubes	50
96 Thermo Matrix 3735, V-bottom, 500µl Storage Tubes	51
96 Thermo Matrix 3744, V-bottom, 500µl ScrewTop Storage Tubes	52
96 Waters 186005837, Clear PolyPro	53

Contents

96AM 250 uL Tip Loading Station	54
96AM Cartridge Rack and Receiver Plate	55
96AM Receiver Plate	56
96AM Cartridge Seating Station.	57
96AM Tip Wash Station.	58
96 V11 Manual Fill Reservoir	59
96 V11 LT 250 Tip Box 19477.002	60
Equilibration Holder + Reservoir, Seahorse 201254-100.	61
Evotip Holder with Evotips	62
Bravo Plate Riser.	63
Universal Lid PS Black	64

Preface

This guide provides a reference to the AssayMAP labware options that appear as selections in the Agilent Protein Sample Prep Workbench software.

This preface contains the following topics:

- [“About this guide” on page vi](#)
- [“Reporting problems” on page viii](#)

About this guide

Overview

This guide provides a reference to the labware that you can use on the AssayMAP Bravo Platform.

Software version

This guide documents labware for the following software versions:

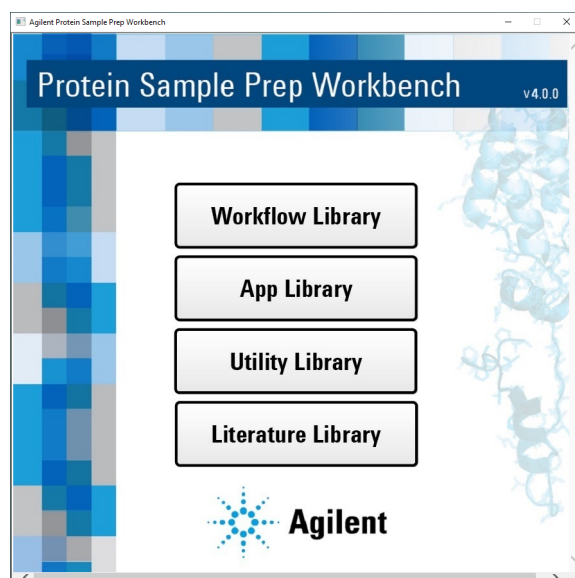
- Protein Sample Prep Workbench 4.0
- VWorks Automation Control 14.1.1

What is new in this revision

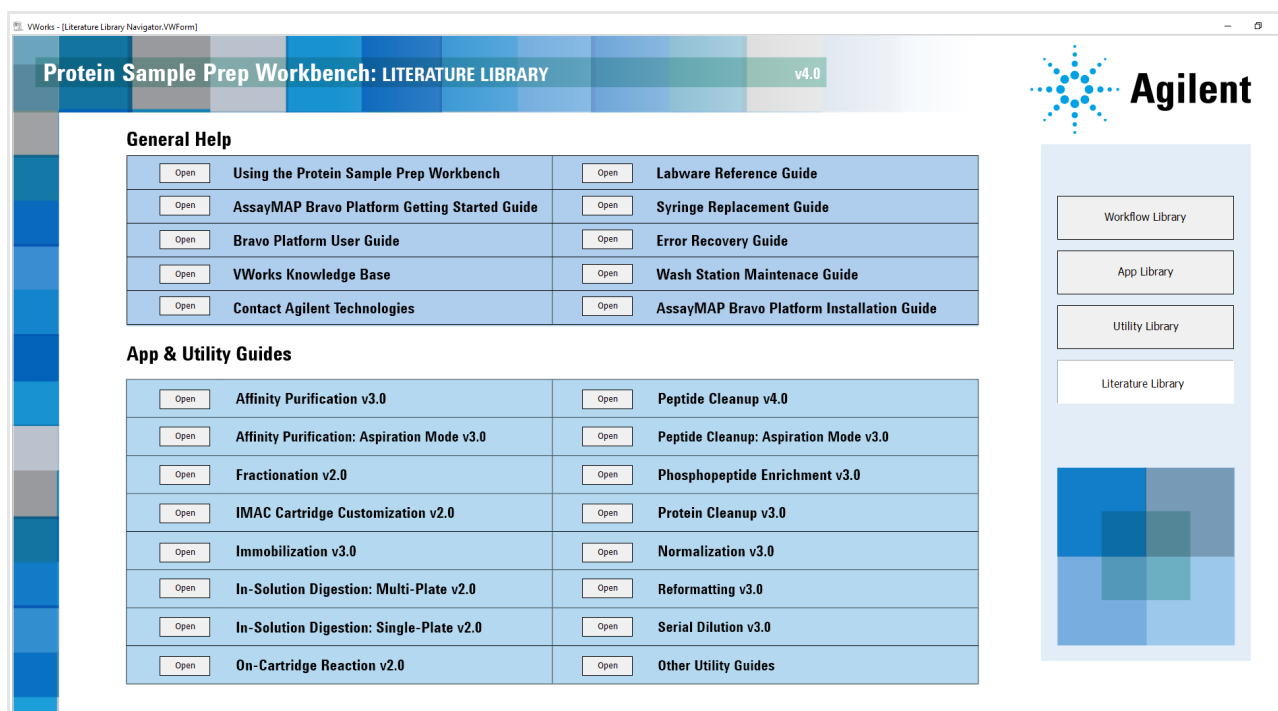
Item	See...
Removed the tables that listed default labware for the apps and utilities.	–
Added the Abgene deep well plate as an option at location 9 in the Affinity Purification app.	"Affinity Purification v3.0 labware options" on page 3
Added labware reference tables.	"Affinity Purification: Aspiration Mode v3.0 labware options" on page 3 "Peptide Cleanup: Aspiration Mode v3.0 labware options" on page 9 "Evosep - 96AM Workflow v1.0 labware options" on page 4
Added the Eppendorf PCR plate as a labware option for locations 4–9 in the Fractionation app.	"Fractionation v2.0 labware options" on page 4
Added the Agilent deep well plate as an option at locations 3 and 7 in the Peptide Cleanup, Phosphopeptide Enrichment, and Protein Cleanup apps.	"Peptide Cleanup v4.0 labware options" on page 9 "Phosphopeptide Enrichment v3.0 labware options" on page 10 "Protein Cleanup v3.0 labware options" on page 11
Updated the photo reference chapter, including <ul style="list-style-type: none">• Added the Agilent 96 Deep Well plate and 96 Channel Wash Station.• Added labware and accessories for the Agilent Evosep - 96AM, Semi-Automated, Full Plate Workflow.• Updated the URL links.	"96 AbGene PCR plate" on page 27 "96 Agilent, 2mL Square Deep Well" on page 30 "96AM Tip Wash Station" on page 58 "Equilibration Holder + Reservoir, Seahorse 201254-100" on page 61 "Evotip Holder with Evotips" on page 62

Accessing the user guides

You can find the user guides for the AssayMAP applications and utilities in the Protein Sample Prep Workbench.



The Literature Library provides links to the user guides.



The app quick start guides are available in the App Library and additional utility guides are available in the Utilities Library.

Reporting problems

If you find a problem with the AssayMAP Bravo Platform, contact Agilent Automation Solutions Technical Support. For contact information, go to <https://www.agilent.com/en/contact-us/page>.

To report problems with...	Have the following information ready
Hardware	<ul style="list-style-type: none">• Instrument serial number from the Bravo serial number label• Short description of the problem
Software	<ul style="list-style-type: none">• Instrument serial number from the Bravo serial number label• Short description of the problem• Relevant software version number (for example, automation control software, diagnostics software, and firmware)• Error message text (or screen capture of the error message dialog box)• Relevant files, such as log files
AssayMAP Cartridges	<ul style="list-style-type: none">• Cartridge type• Lot number• Short description of the problem

For instructions on how to resolve common error messages, see the [AssayMAP Bravo Platform Error Recovery Guide](#).

1 Labware options

This chapter provides a reference to the labware options in the Protein Sample Prep Workbench.

The topics are:

- [“Labware options for each application” on page 2](#)
- [“Labware options for each utility” on page 12](#)

Labware options for each application

The topics in this section are:

- [“AssayMAP Cartridges and pipette tips” on page 2](#)
- [“Affinity Purification v3.0 labware options” on page 3](#)
- [“Affinity Purification: Aspiration Mode v3.0 labware options” on page 3](#)
- [“Evosep - 96AM Workflow v1.0 labware options” on page 4](#)
- [“Fractionation v2.0 labware options” on page 4](#)
- [“IMAC Cartridge Customization v2.0 labware options” on page 4](#)
- [“Immobilization v3.0 labware options” on page 5](#)
- [“In-Solution Digestion: Multi-Plate v2.0 labware options” on page 6](#)
- [“In-Solution Digestion: Single Plate v2.0 labware options” on page 7](#)
- [“On-Cartridge Reaction v2.0 labware options” on page 8](#)
- [“Peptide Cleanup v4.0 labware options” on page 9](#)
- [“Peptide Cleanup: Aspiration Mode v3.0 labware options” on page 9](#)
- [“Phosphopeptide Enrichment v3.0 labware options” on page 10](#)
- [“Protein Cleanup v3.0 labware options” on page 11](#)

IMPORTANT

See the [“Labware photo reference” on page 21](#) for dimensionally equivalent alternatives to the part numbers listed in the following tables.

AssayMAP Cartridges and pipette tips

Most AssayMAP applications require cartridges preloaded in the 96AM Cartridge & Tip Seating Station (Agilent G5409-20025) at deck location 2. For cartridge details, see the [Agilent AssayMAP Cartridge Selection Guide](#).

The In-Solution Digestion applications and several utilities have the option of using 250- μ L pipette tips (Agilent 19477-002) in the seating station.

Affinity Purification v3.0 labware options

Labware	Manufacturer part number	Deck location options
12 Column, Low Profile Reservoir, Natural PP	Agilent 201280-100	3, 5, 6, 7, 8
8 Row, Low Profile Reservoir, Natural PP	Agilent 201282-100	3, 5, 6, 7, 8
96 ABgene 1127, 1mL Deep Well, Square Well, Round Bottom	ABgene AB-1127	3–9
96 Eppendorf 30129300, PCR, Full Skirt, PolyPro	Eppendorf 30129300	4, 7, 9
96 Greiner 652270, PCR, Full Skirt, PolyPro	Greiner 652270	4, 7, 9*
96 Greiner 650201_U-Bottom, Clear PolyPro	Greiner 650201	3–9
96 Greiner 650207_U-Bottom, White PolyPro	Greiner 650207	3–9
96 Greiner 651201_V-Bottom, Clear PolyPro	Greiner 651201	3–9
96 Costar 3363, PP Conical Bottom	Corning Costar 3363	3–9
96 Greiner 675801, Half Area, Flat Bottom, UV Star	Greiner 675801	4, 7, 9
96 V11 Manual Fill Reservoir	Agilent G5498B#049	3, 5, 6, 8
*The Greiner PCR plate is not compatible with the 25 µL cartridges at deck locations 7 and 9.		

Affinity Purification: Aspiration Mode v3.0 labware options

Labware	Manufacturer part number	Deck location options*
12 Column, Low Profile Reservoir, Natural PP	Agilent 201280-100	3–9
8 Row, Low Profile Reservoir, Natural PP	Agilent 201282-100	3–9
96 ABgene 1127, 1mL Deep Well, Square Well, Round Bottom	ABgene AB-1127	3–9
96 Greiner 652270, PCR, Full Skirt, PolyPro	Greiner 652270	3–9
96 Greiner 650201_U-Bottom, Clear PolyPro	Greiner 650201	3–9
96 Greiner 675801, Half Area, Flat Bottom, UV Star	Greiner 675801	3–9
96 V11 Manual Fill Reservoir	Agilent G5498B#049	3–9
*This application has not yet been optimized for the 25 µL cartridges.		

1 Labware options

Labware options for each application

Evosep - 96AM Workflow v1.0 labware options

Labware	Manufacturer part number	Deck location options
250-µL pipette tips	Agilent 19477-002	5, 6*
96 Abgene PCR	Fisher Scientific AB-0800	4
96 Eppendorf 30129300, PCR, Full Skirt, PolyPro	Eppendorf 30129300	4
96 Bio-Rad PCR, Hard-Shell, Low-Profile, Full Skirt	Bio-Rad HSP-9611	4
96 Greiner652270, PCR, Full Skirt, PolyPro	Greiner 652270	4
Reservoir, Seahorse 201254-100, PP, 1 Well, pyramid bottom	Agilent 201254-100	3, 9
Evotip Holder with Evotips	<i>Holder:</i> Evosep EV1142 <i>Evotips:</i> Evosep EV2001	7
Equilibration Holder + Reservoir, Seahorse 201254-100	<i>Holder:</i> Evosep EV1142 <i>Reservoir:</i> Agilent 201254-100	8

*If less than 12 full columns are used in the sample plate, the Tip Setup protocol is required to prepare the tip layout. The Tip Setup protocol uses empty pipette tip boxes at locations 8 and 9.

Fractionation v2.0 labware options

Labware	Manufacturer part number	Deck location options
96 ABgene 1127, 1mL Deep Well, Square Well, Round Bottom	ABgene AB-1127	3
96 Eppendorf 30129300, PCR, Full Skirt, PolyPro	Eppendorf 30129300	4–9
96 Greiner 650201_U-Bottom, Clear PolyPro	Greiner 650201	4–9

IMAC Cartridge Customization v2.0 labware options

Labware	Manufacturer part number	Deck location options
12 Column, Low Profile Reservoir, Natural PP	Agilent 201280-100	3–9
8 Row, Low Profile Reservoir, Natural PP	Agilent 201282-100	3–9
96 ABgene 1127, 1mL Deep Well, Square Well, Round Bottom	ABgene AB-1127	3–8
96 Eppendorf 30129300, PCR, Full Skirt, PolyPro	Eppendorf 30129300	7, 9
96 Greiner 652270, PCR, Full Skirt, PolyPro	Greiner 652270	7, 9*
96 Bio-Rad PCR, Hard-Shell, Low-Profile, Full Skirt	Bio-Rad HSP-9611	7, 9*

Labware	Manufacturer part number	Deck location options
96 Greiner 650201 U-Bottom, Clear PolyPro	Greiner 650201	3–9
96 Greiner 650207_U-Bottom, White PolyPro	Greiner 650207	3–9
96 Greiner 651201_V-Bottom, Clear PolyPro	Greiner 651201	3–9
96 Costar 3363, PP Conical Bottom	Corning Costar 3363	3–9
96 Greiner 675801, Half Area, Flat-Bottom, UV Star	Greiner 675801	7, 9
96 V11 Manual Fill Reservoir	Agilent G5498B#049	4, 5, 6, 8, 9
Reservoir, Axygen Scientific RES-SW96-LP, 86mL, pyramid bottom	Axygen Scientific RES-SW96-LP	4, 5, 6, 8, 9
Reservoir, Seahorse 201254-100, PP, no walls, pyramid bottoms	Agilent 201254-100	4, 5, 6, 8, 9
*The Greiner and BioRad PCR plates are not compatible with the 25 µL cartridges at deck location 7.		

Immobilization v3.0 labware options

Labware	Manufacturer part number	Deck location options
12 Column, Low Profile Reservoir, Natural PP	Agilent 201280-100	3, 5, 6, 7, 8
8 Row, Low Profile Reservoir, Natural PP	Agilent 201282-100	3, 5, 6, 7, 8
96 ABgene 1127, 1mL Deep Well, Square Well, Round Bottom	ABgene AB-1127	3–8
96 Eppendorf 30129300, PCR, Full Skirt, PolyPro	Eppendorf 30129300	4, 7, 9
96 Greiner 652270, PCR, Full Skirt, PolyPro	Greiner 652270	4, 7, 9*
96 Greiner 650201_U-Bottom, Clear PolyPro	Greiner 650201	3–9
96 Greiner 650207_U-Bottom, White PolyPro	Greiner 650207	3–9
96 Greiner 651201_V-Bottom, Clear PolyPro	Greiner 651201	3–9
96 Costar 3363, PP Conical Bottom	Corning Costar 3363	3–9
96 Greiner 675801, Half Area, Flat-Bottom, UV Star	Greiner 675801	4, 7, 9
96 V11 Manual Fill Reservoir	Agilent G5498B#049	3, 5, 6, 8
*The Greiner PCR plate is not compatible with the 25 µL cartridges at deck location 7.		

1 Labware options

Labware options for each application

In-Solution Digestion: Multi-Plate v2.0 labware options

Labware	Manufacturer part number	Deck location options
Reagent Setup		
96 Red PCR Insert + 96 Eppendorf 30129300, PCR, Full Skirt	Agilent insert G5498B#013 and Eppendorf 30129300	4
96 Greiner 650207_U-Bottom, White PolyPro	Greiner 650207	5, 7
250- μ L pipette tips	Agilent 19477-002	6
96 ABgene 1127, Deep Well, Square Well, Round Bottom	ABgene AB-1127	8
In-Solution Digestion		
96 Greiner 650207_U-Bottom, White PolyPro	Greiner 650207	2, 5–9
96 ABgene 1127, Deep Well, Square Well, Round Bottom	ABgene AB-1127	3
96 Red PCR Insert + 96 Eppendorf 30129300, PCR, Full Skirt	Agilent insert G5498B#013 and Eppendorf 30129300	4

In-Solution Digestion: Single Plate v2.0 labware options

Labware	Manufacturer part number	Deck location options
Lid, Universal*	Agilent 200858-100	4–9
96 Eppendorf 30129300, PCR, Full Skirt, PolyPro	Eppendorf 30129300	4–9
96 Bio-Rad PCR, Hard-Shell, Low-Profile, Full Skirt	Bio-Rad HSP-9611	4–9
96 Greiner 652270, PCR, Full Skirt, PolyPro	Greiner 652270	4–9
96 Greiner 650201_U-Bottom, Clear PolyPro	Greiner 650201	4–9
96 Greiner 650207_U-Bottom, White PolyPro	Greiner 650207	4–9
96 Greiner 651201_V-Bottom, Clear PolyPro	Greiner 651201	4–9
96 Costar 3363, PP Conical Bottom	Corning Costar 3363	4–9
96 ABgene 1127, 1mL Deep Well, Square Well, Round Bottom	ABgene AB-1127	4–9
96 Eppendorf 96-500_V-bottom, Clear PolyPro	Eppendorf 96/500	4–9
96 Eppendorf 96-1000_U-bottom, Clear PolyPro	Eppendorf 96/1000	4–9
96 Waters 186005837, Clear PolyPro	Waters 186005837	4–9
12 Column, Low Profile Reservoir, Natural PP Lidded	Agilent 201280-100	5–9
8 Row, Low Profile Reservoir, Natural PP Lidded	Agilent 201282-100	5–9
Reservoir, Seahorse 201254-100, PP, 1 Well, pyramid bottom	Agilent 201254-100	5–9
96 Red PCR Insert + Eppendorf 30129300, PCR, Full Skirt	Agilent insert (provided) and Eppendorf 30129300	4
96 Red PCR Insert + Bio-Rad PCR, Hard-Shell, Low-Profile, Full Skirt	Agilent insert (provided) and Bio-Rad HSP-9611	4
96 Red PCR Insert + Greiner 652270, PCR, Full Skirt, PolyPro	Agilent insert (provided) and Greiner 652270	4
96 Greiner U-Bottom Thermal Insert + Greiner 650201_U-Bottom, Clear PolyPro	Agilent G5498B#126 and Greiner 650201	4
96 Greiner V-Bottom Thermal Insert + Greiner 651201_V-Bottom, Clear PolyPro	Agilent G5498B#126 and Greiner 651201	4
96 AbGene U-Bottom Thermal Insert + AbGene 1127, Square Well, Round Bottom	Agilent G5498B#127 and ABgene AB-1127	4
*Lidded plates are allowed at these deck locations, not the lid by itself. If you plan to use a lidded plate, ensure that you use only the approved lid type.		

1 Labware options

Labware options for each application

On-Cartridge Reaction v2.0 labware options

Labware	Manufacturer part number	Deck location options
12 Column, Low Profile Reservoir, Natural PP	Agilent 201280-100	3–9
8 Row, Low Profile Reservoir, Natural PP	Agilent 201282-100	3–9
Reservoir, Seahorse 201254-100, PP, no walls, pyramid bottoms	Agilent 201254-100	3–9
Reservoir, Axygen Scientific RES-SW96-LP, 86mL pyramid bottom	Axygen Scientific RES-SW96-LP	3–9
96 ABgene 1127, 1mL Deep Well, Square Well, Round Bottom	ABgene AB-1127	3–9
96 Bio-Rad PCR, Hard-Shell, Low-Profile, Full Skirt	Bio-Rad HSP-9611	3–9*
96 Eppendorf 30129300, PCR, Full Skirt, PolyPro	Eppendorf 30129300	3–9
96 Greiner 652270, PCR, Full Skirt, PolyPro	Greiner 652270	3–9*
96 Greiner 650201_U-Bottom, Clear PolyPro	Greiner 650201	3–9
96 Greiner 650207_U-Bottom, White PolyPro	Greiner 650207	3–9
96 Greiner 651201_V-Bottom, Clear PolyPro	Greiner 651201	3–9
96 Costar 3363, PP Conical Bottom	Corning Costar 3363	3–9
96 Greiner 675801, Half Area, Flat-Bottom, UV Star	Greiner 675801	3–9
96 V11 Manual Fill Reservoir	Agilent G5498B#049	3–9
96 Red PCR Insert + Eppendorf 30129300, PCR, Full Skirt	Agilent insert (provided) and Eppendorf 30129300	4
96 Red PCR Insert + Bio-Rad PCR, Hard-Shell, Low-Profile, Full Skirt	Agilent insert (provided) and Bio-Rad HSP-9611	4*
96 Red PCR Insert + Greiner 652270, PCR, Full Skirt, PolyPro	Agilent insert (provided) and Greiner 652270	4*
96 Greiner U-Bottom Thermal Insert + Greiner 650201_U-Bottom, Clear PolyPro	Agilent G5498B#126 and Greiner 650201	4
96 Greiner V-Bottom Thermal Insert + Greiner 651201 V-Bottom, Clear PolyPro	Agilent G5498B#126 and Greiner 651201	4
96 AbGene U-Bottom Thermal Insert + AbGene 1127, Square Well, Round Bottom	Agilent G5498B#127 and ABgene AB-1127	4
*The Greiner and BioRad PCR plates are not compatible with the 25 µL cartridges at deck locations 4, 7, and 9.		

Peptide Cleanup v4.0 labware options

Labware	Manufacturer part number	Deck location options
12 Column, Low Profile Reservoir, Natural PP	Agilent 201280-100	3, 5–8
8 Row, Low Profile Reservoir, Natural PP	Agilent 201282-100	3, 5–8
96 Agilent, 2mL Square Deep Well labware	Agilent 204353-100	3, 7
96 ABgene 1127, 1mL Deep Well, Square Well, Round Bottom	ABgene AB-1127	3–8
96 Eppendorf 30129300, PCR, Full Skirt, PolyPro	Eppendorf 30129300	4, 7, 9
96 Greiner 652270, PCR, Full Skirt, PolyPro	Greiner 652270	4, 7, 9*
96 Greiner 650201_U-Bottom, Clear PolyPro	Greiner 650201	3–9
96 Greiner 650207_U-Bottom, White PolyPro	Greiner 650207	3–9
96 Greiner 651201_V-Bottom, Clear PolyPro	Greiner 651201	3–9
96 Costar 3363, PP Conical Bottom	Corning Costar 3363	3–9
96 Greiner 675801, Half Area, Flat-Bottom, UV Star	Greiner 675801	4, 7, 9
96 V11 Manual Fill Reservoir	Agilent G5498B#049	5, 6, 8
*The Greiner PCR plate is not compatible with the 25 µL cartridges at deck locations 7 and 9.		

Peptide Cleanup: Aspiration Mode v3.0 labware options

Labware	Manufacturer part number	Deck location options*
12 Column, Low Profile Reservoir, Natural PP	Agilent 201280-100	3, 5–8
8 Row, Low Profile Reservoir, Natural PP	Agilent 201282-100	5–8
96 ABgene 1127, 1mL Deep Well, Square Well, Round Bottom	ABgene AB-1127	3–8
96 Eppendorf 30129300, PCR, Full Skirt, PolyPro	Eppendorf 30129300	4, 7, 9
96 Greiner 652270, PCR, Full Skirt, PolyPro	Greiner 652270	4, 7, 9
96 Greiner 650201_U-Bottom, Clear PolyPro	Greiner 650201	3–9
96 Greiner 650207_U-Bottom, White PolyPro	Greiner 650207	3–9
96 Greiner 675801, Half Area, Flat-Bottom, UV Star	Greiner 675801	4, 7, 9
96 V11 Manual Fill Reservoir	Agilent G5498B#049	5–8
*This application has not yet been optimized for the 25 µL cartridges.		

1 Labware options

Labware options for each application

Phosphopeptide Enrichment v3.0 labware options

Labware	Manufacturer part number	Deck location options*
12 Column, Low Profile Reservoir, Natural PP	Agilent 201280-100	3, 5–8
8 Row, Low Profile Reservoir, Natural PP	Agilent 201282-100	3, 5–8
96 Agilent, 2mL Square Deep Well labware	Agilent 204353-100	3, 7
96 ABgene 1127, 1mL Deep Well, Square Well, Round Bottom	ABgene AB-1127	3–8
96 Eppendorf 30129300, PCR, Full Skirt, PolyPro	Eppendorf 30129300	4, 7, 9
96 Greiner 652270, PCR, Full Skirt, PolyPro	Greiner 652270	4, 7, 9*
96 Bio-Rad PCR, Hard-Shell, Low-Profile, Full Skirt	Bio-Rad HSP-9611	4, 7, 9*
96 Greiner 650201_U-Bottom, Clear PolyPro	Greiner 650201	3–9
96 Greiner 650207_U-Bottom, White PolyPro	Greiner 650207	3–9
96 Greiner 651201_V-Bottom, Clear PolyPro	Greiner 651201	3–9
96 Costar 3363, PP Conical Bottom	Corning Costar 3363	3–9
96 Greiner 675801, Half Area, Flat-Bottom, UV Star	Greiner 675801	4, 7, 9
96 V11 Manual Fill Reservoir	Agilent G5498B#049	5, 6, 8
Reservoir, Axygen Scientific RES-SW96-LP, 86mL pyramid bottom	Axygen Scientific RES-SW96-LP	5, 6, 8
Reservoir, Seahorse 201254-100, PP, no walls, pyramid bottoms	Agilent 201254-100	5, 6, 8

*The Greiner and BioRad PCR plates are not compatible with the 25 µL cartridges at deck locations 7 and 9.

Protein Cleanup v3.0 labware options

Labware	Manufacturer part number	Deck location options*
12 Column, Low Profile Reservoir, Natural PP	Agilent 201280-100	3, 5–8
8 Row, Low Profile Reservoir, Natural PP	Agilent 201282-100	3, 5–8
96 Agilent, 2mL Square Deep Well labware	Agilent 204353-100	3, 7
96 ABgene 1127, 1mL Deep Well, Square Well, Round Bottom	ABgene AB-1127	3–8
96 Eppendorf 30129300, PCR, Full Skirt, PolyPro	Eppendorf 30129300	4, 7, 9
96 Greiner 652270, PCR, Full Skirt, PolyPro	Greiner 652270	4, 7, 9*
96 Greiner 650201_U-Bottom, Clear PolyPro	Greiner 650201	3–9
96 Greiner 650207_U-Bottom, White PolyPro	Greiner 650207	3–9
96 Greiner 651201_V-Bottom, Clear PolyPro	Greiner 651201	3–9
96 Costar 3363, PP Conical Bottom	Corning Costar 3363	3–9
96 Greiner 675801, Half Area, Flat-Bottom, UV Star	Greiner 675801	4, 7, 9
96 V11 Manual Fill Reservoir	Agilent G5498B#049	5, 6, 8
*The Greiner PCR plate is not compatible with the 25 µL cartridges at deck locations 7 and 9.		

Labware options for each utility

This section includes the following tables:

- “Cartridge Transfer utility v2.0 required labware” on page 12
- “Pipette Tip Transfer utility v2.0 required labware” on page 12
- “Reagent Aliquot utility v2.0 labware options” on page 14
- “Reagent Transfer utility v3.0 labware options” on page 15
- “Reformatting utility v3.0 labware options” on page 16
- “Serial Dilution utility v3.0 labware options” on page 17
- “Single Liquid Addition utility v2.0 labware options” on page 18
- “Syringe Test utility v2.0 labware options” on page 18
- “Syringe Wash utility v3.0 labware options” on page 19
- “System Startup/Shutdown utility v3.0 labware options” on page 19

IMPORTANT

See the “[Labware photo reference](#)” on page 21 for dimensionally equivalent alternatives to the part numbers listed in the following tables.

Cartridge Transfer utility v2.0 required labware

Labware	Manufacturer part number	Deck location
96AM Cartridge & Tip Seating Station	Agilent G5409-20025	2
96AM Cartridge Rack and Receiver Plate with cartridges (all types)	See Agilent AssayMAP Cartridge Selection Guide	6

Pipette Tip Transfer utility v2.0 required labware

Labware	Manufacturer part number	Deck location
96AM Cartridge & Tip Seating Station	Agilent G5409-20025	2
96 V11 LT 250 Tip Box (250-µL pipette tips)	Agilent 19477-002	6

Normalization v3.0 labware options

Labware	Manufacturer part number	Deck location options*
96 V11 LT 250 Tip Box (250-µL pipette tips)	Agilent 19477-002	3
96AM Cartridge & Tip Seating Station	Agilent G5409-20025	5
96 ABgene 1127, 1mL Deep Well, Square Well, Round Bottom	ABgene AB-1127	2, 6
96 Eppendorf 30129300, PCR, Full Skirt, PolyPro	Eppendorf 30129300	2, 6
96 Greiner 652270, PCR, Full Skirt, PolyPro	Greiner 652270	2, 6
96 Greiner 650201_U-Bottom, Clear PolyPro	Greiner 650201	2, 6
96 Greiner 650207_U-Bottom, White PolyPro	Greiner 650207	2, 6
96 Greiner 651201_V-Bottom, Clear PolyPro	Greiner 651201	2, 6
96 Costar 3363, PP Conical Bottom	Corning Costar 3363	2, 6
96 Greiner 675801, Half Area, Flat-Bottom, UV Star	Greiner 675801	2, 6
12 Column, Low Profile Reservoir, Natural PP	Agilent 201280-100	4
8 Row, Low Profile Reservoir, Natural PP	Agilent 201282-100	4
96 V11 Manual Fill Reservoir	Agilent G5498B#049	4
Reservoir, Axygen Scientific RES-SW96-LP, 86mL pyramid bottom	Axygen Scientific RES-SW96-LP	4
*At deck locations 2 and 6 the labware must be seated on Plate Risers.		

1 Labware options

Labware options for each utility

Reagent Aliquot utility v2.0 labware options

Labware	Manufacturer part number	Deck location options
96AM Cartridge & Tip Seating Station	Agilent G5409-20025	2
96 V11 LT 250 Tip Box (250-µL pipette tips) (required)	Agilent 19477-002	6
96 Eppendorf 30129300, PCR, Full Skirt, PolyPro	Eppendorf 30129300	4, 7
96 Bio-Rad PCR, Hard-Shell, Low-Profile, Full Skirt	Bio-Rad HSP-9611	4, 7
96 Greiner 652270, PCR, Full Skirt, PolyPro	Greiner 652270	4, 7
96 ABgene 1127, 1mL Deep Well, Square Well, Round Bottom	ABgene AB-1127	4, 7
96 Eppendorf 96-500 V-bottom, Clear, PolyPro	Eppendorf 96/500	4, 7
96 Eppendorf 96-1000 U-bottom, Clear, PolyPro	Eppendorf 96/1000	4, 7
96 Waters 186005837, Clear PolyPro	Waters 186005837	4, 7
96 Greiner 650201_U-Bottom, Clear PolyPro	Greiner 650201	4, 7
96 Greiner 650207_U-Bottom, White PolyPro	Greiner 650207	4, 7
96 Greiner 651201_V-Bottom, Clear PolyPro	Greiner 651201	4, 7
96 Costar 3363, PP Conical Bottom	Corning Costar 3363	4, 7
12 Column, Low Profile Reservoir, Natural PP	Agilent 201280-100	7
8 Row, Low Profile Reservoir, Natural PP	Agilent 201282-100	7
Reservoir, Seahorse 201254-100, PP, no walls, pyramid bottoms	Agilent 201254-100	7
Reservoir, Axygen Scientific RES-SW96-LP, 86mL pyramid bottom	Axygen Scientific RES-SW96-LP	7
96 V11 Manual Fill Reservoir	Agilent G5498B#049	7

Reagent Transfer utility v3.0 labware options

Labware	Manufacturer part number	Deck location options
96AM Cartridge & Tip Seating Station	Agilent G5409-20025	2
96 V11 LT 250 Tip Box 19477-002	Agilent 19477-002	6
96 Eppendorf 30129300, PCR, Full Skirt, PolyPro	Eppendorf 30129300	7, 8
96 Bio-Rad PCR, Hard-Shell, Low-Profile, Full Skirt	Bio-Rad HSP-9611	7, 8
96 Greiner 652270, PCR, Full Skirt, PolyPro	Greiner 652270	7, 8
96 ABgene 1127, 1mL Deep Well, Square Well, Round Bottom	ABgene AB-1127	7, 8
96 Greiner 650201_U-Bottom, Clear PolyPro	Greiner 650201	7, 8
96 Greiner 650207_U-Bottom, White PolyPro	Greiner 650207	7, 8
96 Greiner 651201_V-Bottom, Clear PolyPro	Greiner 651201	7, 8
96 Costar 3363, PP Conical Bottom	Corning Costar 3363	7, 8
96 Greiner 675801, Half Area, Flat-Bottom, UV Star	Greiner 675801	7, 8
12 Column, Low Profile Reservoir, Natural PP	Agilent 201280-100	7, 8
8 Row, Low Profile Reservoir, Natural PP	Agilent 201282-100	7, 8
Reservoir, Seahorse 201254-100, PP, no walls, pyramid bottom	Agilent 201254-100	7, 8
Reservoir, Axygen Scientific RES-SW96-LP, 86mL pyramid bottom	Axygen Scientific RES-SW96-LP	7, 8
96 V11 Manual Fill Reservoir	Agilent G5498B#049	7, 8
96AM Receiver Plate	Included with Agilent cartridge rack	7, 8
96 Eppendorf 96-500 V-bottom, Clear, PolyPro*	Eppendorf 96/500	7, 8
96 Eppendorf 96-1000 U-bottom, Clear, PolyPro*	Eppendorf 96/1000	7, 8
96 Waters 186005837, Clear PolyPro*	Waters 186005837	7, 8

*Pipette tips are required for reagent transfers using these deep-well plate options.

1 Labware options

Labware options for each utility

Reformatting utility v3.0 labware options

Labware	Manufacturer part number	Deck location options*
96 V11 LT 250 Tip Box (250-µL pipette tips)	Agilent 19477-002	3
96AM Cartridge & Tip Seating Station	Agilent G5409-20025	5
96 ABgene 1127, 1mL Deep Well, Square Well, Round Bottom	ABgene AB-1127	2, 6
96 Bio-Rad PCR, Hard-Shell, Low-Profile, Full Skirt	Bio-Rad HSP-9611	2, 6
96 Costar 3363, PP Conical Bottom	Corning Costar 3363	2, 6
96 Eppendorf 30129300, PCR, Full Skirt, PolyPro	Eppendorf 30129300	2, 6
96 Eppendorf 96-500_V-bottom, Clear PolyPro	Eppendorf 96/500	2, 6
96 Greiner 650201_U-Bottom, Clear PolyPro	Greiner 650201	2, 6
96 Greiner 650207_U-Bottom, White PolyPro	Greiner 650207	2, 6
96 Greiner 651201_V-Bottom, Clear PolyPro	Greiner 651201	2, 6
96 Greiner 652270, PCR, Full Skirt, PolyPro	Greiner 652270	2, 6
96 Greiner 675801, Half Area, Flat-Bottom, UV Star	Greiner 675801	2, 6
96 Thermo Matrix 3732, V-bottom, 0.75ml Storage Tubes	Thermo Fisher Scientific 3732	2, 6
96 Thermo Matrix 3735, V-bottom, 500µl Storage Tubes	Thermo Fisher Scientific 3735	2, 6
96 Thermo Matrix 3744, V-bottom, 500µl ScrewTop Storage Tubes	Thermo Fisher Scientific 3744	2, 6
96 Waters 186005837, Clear PolyPro	Waters 186005837	2, 6
12 Column, Low Profile Reservoir, Natural PP	Agilent 201280-100	2, 6
8 Row, Low Profile Reservoir, Natural PP	Agilent 201282-100	2, 6
*At deck locations 2 and 6 the labware must be seated on Plate Risers.		

Serial Dilution utility v3.0 labware options

Labware	Manufacturer part number	Deck location options*
96 V11 LT 250 Tip Box (250-µL pipette tips)	Agilent 19477-002	3
96AM Cartridge & Tip Seating Station	Agilent G5409-20025	5
8 Row, Low Profile Reservoir, Natural PP	Agilent 201282-100	2
12 Column, Low Profile Reservoir, Natural PP	Agilent 201280-100	2
96 V11 Manual Fill Reservoir	Agilent G5498B#049	2
Reservoir, Seahorse 201254-100, PP, no walls, pyramid bottom	Agilent 201254-100	2
96 AbGene 1127, 1mL Deep Well, Square Well, Round Bottom	ABgene AB-1127	2, 6
96 Costar 3363, PP Conical Bottom	Corning Costar 3363	2, 6
96 Eppendorf 30129300, PCR, Full Skirt, PolyPro	Eppendorf 30129300	2, 6
96 Eppendorf 96-500_V-bottom, Clear PolyPro	Eppendorf 96/500	2, 6
96 Bio-Rad PCR, Hard-Shell, Low-Profile, Full Skirt	Bio-Rad HSP-9611	2, 6
96 Greiner 650201_U-Bottom, Clear PolyPro	Greiner 650201	2, 6
96 Greiner 650207_U-Bottom, White PolyPro	Greiner 650207	2, 6
96 Greiner 651201_V-Bottom, Clear PolyPro	Greiner 651201	2, 6
96 Greiner 652270, PCR, Full Skirt, PolyPro	Greiner 652270	2, 6
96 Greiner 675801, Half Area, Flat-Bottom, UV Star	Greiner 675801	6
96 Thermo Matrix 3732, V-bottom, 0.75ml Storage Tubes	Thermo Fisher Scientific 3732	2, 6
96 Thermo Matrix 3735, V-bottom, 500µl Storage Tubes	Thermo Fisher Scientific 3735	2, 6
96 Thermo Matrix 3744, V-bottom, 500µl ScrewTop Storage Tubes	Thermo Fisher Scientific 3744	2, 6
96 Waters 186005837, Clear PolyPro	Waters 186005837	2, 6

*At deck locations 2 and 6 the labware must be seated on Plate Risers.

1 Labware options

Labware options for each utility

Single Liquid Addition utility v2.0 labware options

Labware	Manufacturer part number	Deck location options
96 Greiner 650207_U-Bottom, White PolyPro	Greiner 650207	2, 5, 6, 8, 9
96 Eppendorf 30129300, PCR, Full Skirt, PolyPro	Eppendorf 30129300	4
96 Greiner 650201_U-Bottom, Clear PolyPro	Greiner 650201	4
96 Greiner 652270, PCR, Full Skirt, PolyPro	Greiner 652270	4
96 Costar 3363, PP Conical Bottom	Corning Costar 3363	4
96 ABgene 1127, 1mL Deep Well, Square Well, Round Bottom	ABgene AB-1127	4
12 Column, Low Profile Reservoir, Natural PP	Agilent 201280-100	4
8 Row, Low Profile Reservoir, Natural PP	Agilent 201282-100	4
96 V11 Manual Fill Reservoir	Agilent G5498B#049	4

Syringe Test utility v2.0 labware options

Labware	Manufacturer part number	Deck location options
96AM Cartridge Seating Station (required)	Agilent G5409-20025	2
Reservoir, Seahorse 201254-100, PP, no walls, pyramid bottom	Agilent 201254-100	4, 7
Reservoir, Axygen Scientific RES-SW96-LP, 86mL pyramid bottom	Axygen Scientific RES-SW96-LP	4, 7
12 Column, Low Profile Reservoir, Natural PP	Agilent 201280-100	4, 7
96 V11 Manual Fill Reservoir	Agilent G5498B#049	4, 7
Resin Free cartridges in 96AM Cartridge Rack and Receiver Plate (required)	Agilent G5496-60008	6
96 Greiner 655101 PS Clr Rnd Well Flat Btm	Greiner 655101	8, 9
96 Nunc 269620, Flat Bottom, Polystyrene	Nunc 269620	8, 9
96 Corning 9017, Flat bottom, polystyrene, clear	Corning 9017	8, 9

Syringe Wash utility v3.0 labware options

Labware	Manufacturer part number	Deck location options
96AM Cartridge & Tip Seating Station	Agilent G5409-20025	2
12 Column, Low Profile Reservoir, Natural PP	Agilent 201280-100	3–9
8 Row, Low Profile Reservoir, Natural PP	Agilent 201282-100	3–9
96 Greiner 650201 U-Bottom, Standard, PolyPro	Greiner 650201	3–9
96 ABgene 1127, 1mL Deep Well, Square Well, Round Bottom	ABgene AB-1127	3–9
Reservoir, Seahorse 201254-100, PP, no walls, pyramid bottom	Agilent 201254-100	3–9
Reservoir, Axygen Scientific RES-SW96-LP, 86mL pyramid bottom	Axygen Scientific RES-SW96-LP	3–9
96 V11 Manual Fill Reservoir	Agilent G5498B#049	3–9

System Startup/Shutdown utility v3.0 labware options

Labware	Manufacturer part number	Deck location options
96AM Cartridge & Tip Seating Station	Agilent G5409-20025	2
12 Column, Low Profile Reservoir, Natural PP	Agilent 201280-100	7
8 Row, Low Profile Reservoir, Natural PP	Agilent 201282-100	7
Reservoir, Seahorse 201254-100, PP, no walls, pyramid bottoms	Agilent 201254-100	7
Reservoir, Axygen Scientific RES-SW96-LP, 86mL pyramid bottom	Axygen Scientific RES-SW96-LP	7
96 Greiner 650201, U-Bottom, Standard, PolyPro	Greiner 650201	7
96 ABgene 1127, 1mL Deep Well, Square Well, Round Bottom	ABgene AB-1127	7
96 V11 Manual Fill Reservoir	Agilent G5498B#049	7

1 Labware options

Labware options for each utility

This page is intentionally blank.

2 Labware photo reference

This chapter provides a photograph and description of each labware used in the Protein Sample Prep Workbench.

- “12 Column, Low Profile Reservoir, Natural PP” on page 23
- “8 Row, Low Profile Reservoir, Natural PP” on page 24
- “Reservoir, Seahorse 201254-100, PP, no walls, pyramid bottoms” on page 25
- “Reservoir, Axygen Scientific RES-SW96-LP, 86mL pyramid bottom” on page 26
- “96 AbGene PCR plate” on page 27
- “96 AbGene 1127, 1mL Deep Well, Square Well, Round Bottom” on page 28
- “96 AbGene U-Bottom Thermal Insert + AbGene 1127, Square Well, Round Bottom” on page 29
- “96 Agilent, 2mL Square Deep Well” on page 30
- “96 Bio-Rad PCR, Hard-Shell, Low-Profile, Full Skirt” on page 31
- “96 Red PCR Insert + Bio-Rad PCR, Hard-Shell, Low-Profile, Full Skirt” on page 32
- “96 Corning 9017, Flat Bottom, Polystyrene, Clear” on page 33
- “96 Costar 3363, PP Conical Bottom” on page 34
- “96 Eppendorf 30129300, PCR, Full Skirt, PolyPro” on page 35
- “96 Red PCR Insert + Eppendorf 30129300, PCR, Full Skirt” on page 36
- “96 Eppendorf 96-1000_U-bottom, Clear PolyPro” on page 37
- “96 Eppendorf 96-500_V-bottom, Clear PolyPro” on page 38
- “96 Greiner 650201, U-Bottom Standard, PolyPro” on page 39
- “96 Greiner 650201_U-Bottom, Clear PolyPro” on page 40
- “96 Greiner U-Bottom Thermal Insert + Greiner 650201_U-Bottom, Clear PolyPro” on page 41
- “96 Greiner 650207_U-Bottom, White PolyPro” on page 42
- “96 Greiner 651201_V-Bottom, Clear PolyPro” on page 43
- “96 Greiner V-Bottom Thermal Insert + Greiner 651201, V-Bottom, Clear PolyPro” on page 44
- “96 Greiner 652270, PCR, Full Skirt, PolyPro” on page 45
- “96 Red PCR Insert + Greiner 652270, PCR, Full Skirt, PolyPro” on page 46
- “96 Greiner 655101 PS Clr Rnd Well Flat Btm” on page 47
- “96 Greiner 675801, Half Area, Flat Bottom, UV Star” on page 48
- “96 Nunc 269620, Flat Bottom, Polystyrene” on page 49
- “96 Thermo Matrix 3732, V-bottom, 0.75ml Storage Tubes” on page 50
- “96 Thermo Matrix 3735, V-bottom, 500µl Storage Tubes” on page 51
- “96 Thermo Matrix 3744, V-bottom, 500µl ScrewTop Storage Tubes” on page 52

2 Labware photo reference

- “96 Waters 186005837, Clear PolyPro” on page 53
- “96AM 250 uL Tip Loading Station” on page 54
- “96AM Cartridge Rack and Receiver Plate” on page 55
- “96AM Receiver Plate” on page 56
- “96AM Cartridge Seating Station” on page 57
- “96AM Tip Wash Station” on page 58
- “96 V11 Manual Fill Reservoir” on page 59
- “96 V11 LT 250 Tip Box 19477.002” on page 60
- “Equilibration Holder + Reservoir, Seahorse 201254-100” on page 61
- “Evotip Holder with Evotips” on page 62
- “Bravo Plate Riser” on page 63
- “Universal Lid PS Black” on page 64

12 Column, Low Profile Reservoir, Natural PP



Manufacturer: Agilent Technologies

Part number: 201280-100

Name of labware type in the software:

12 Column, Low Profile Reservoir, Natural PP or
12 Seahorse, Low Profile, Reservoir, Natural PP

General information

One of the most common reagent reservoir options used in the Protein Sample Prep Workbench. The column-wise partitions enable partial-plate runs with the same full-column convention that is used to define the number of full columns of cartridges or pipette tips in most applications and utilities.

Volume recommendations

Number of wells = 12

Maximum volume per well = 7 mL

Recommended overage per well = 3 mL

Additional notes

A dimensionally equivalent alternative is the [Axygen \(Corning\) RES-MW12-LP](#).

Resources

For information about...	See...
Ordering	https://www.agilent.com/store/#quickOrder

8 Row, Low Profile Reservoir, Natural PP



Manufacturer: Agilent Technologies

Part number: 201282-100

Name of labware type in the software:

8 Row, Low Profile Reservoir, Natural PP or

8 Seahorse, Low Profile Reservoir, Natural PP

General information

This is a common reagent reservoir option in the Protein Sample Prep Workbench. The row-wise partitions enable partial-plate runs based on full rows of 12 samples. This reservoir also makes it easy to isolate reagents if experimentation involves testing reagents with different compositions.

Volume recommendations

Number of wells = 8

Maximum volume per well = 10.6 mL

Recommended overage per well = 4.5 mL

Additional notes

A dimensionally equivalent alternative is the [Axygen \(Corning\) RES-MW8-LP](#).

Resources

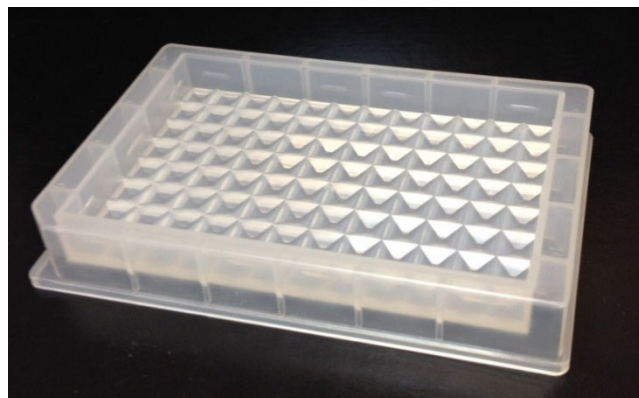
For information about...

See...

Ordering

<https://www.agilent.com/store/#quickOrder>

Reservoir, Seahorse 201254-100, PP, no walls, pyramid bottoms



Manufacturer: Agilent Technologies

Part number: 201254-100

Name of labware type in the software:
Reservoir, Seahorse 201254-100, PP, no walls,
pyramid bottoms

General information

This common reservoir is very useful for AssayMAP operations. It enables easy filling when running a full plate of 96 samples.

Volume recommendations

Number of wells = 1

Maximum volume per well = 86 mL

Recommended overage per well = 20 mL

Additional notes

A dimensionally equivalent alternative is the [Axygen \(Corning\) RES-SW96-LP](#).

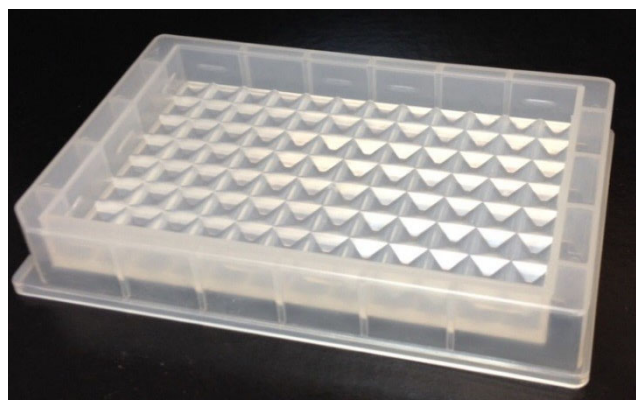
Resources

For information about...	See...
Ordering	https://www.agilent.com/store/#quickOrder

2 Labware photo reference

Reservoir, Axygen Scientific RES-SW96-LP, 86mL pyramid bottom

Reservoir, Axygen Scientific RES-SW96-LP, 86mL pyramid bottom



Manufacturer: Axygen Scientific

Part number: RES-SW96-LP

Name of labware type in the software:

Reservoir, Axygen Scientific RES-SW96-LP, 86mL pyramid bottom

General information

This common reservoir is very useful for AssayMAP operations. It enables easy filling when running a full plate of 96 samples.

Volume recommendations

Number of wells = 1

Maximum volume per well = 86 mL

Recommended overage per well = 20 mL

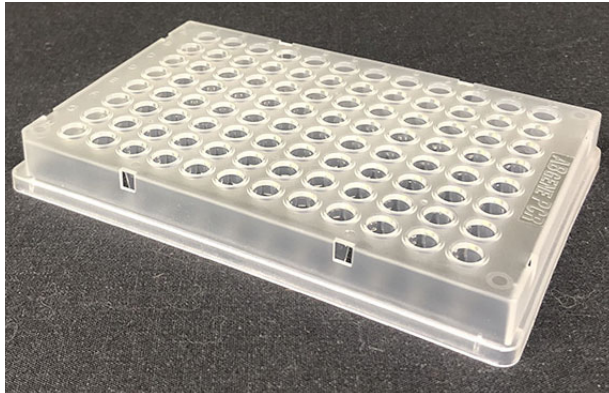
Additional notes

A dimensionally equivalent alternative is the [Reservoir, Seahorse 201254-100, PP, no walls, pyramid bottoms](#).

Resources

For information about...	See...
Ordering	https://ecatalog.corning.com/life-sciences/b2c/US/en/Genomics-%26-Molecular-Biology/Automation-Consumables/Automation-Reservoirs/Axygen%C2%AE-Reagent-Reservoirs/p/RES-SW96-LP

96 AbGene PCR plate



Manufacturer: Fisher Scientific

Part number: AB-0800

Name of labware type in the software:
96 AbGene PCR

General information

The AbGene PCR plate is an option for the sample plate in the AssayMAP Evosep - 96AM, Semi-Automated, Full Plate Workflow.

Volume recommendations

Number of wells = 96

Maximum volume per well = 200 μ L

Recommended overage per well = 10 μ L

Resources

For information about...	See...
Product page	https://www.thermofisher.com/order/catalog/product/AB0800
Evosep loading protocol	https://www.evosep.com

2 Labware photo reference

96 AbGene 1127, 1mL Deep Well, Square Well, Round Bottom

96 AbGene 1127, 1mL Deep Well, Square Well, Round Bottom



Manufacturer: Fisher Scientific

Part number: AB-1127

Name of labware type in the software:

96 AbGene 1127, 1mL Deep Well, Square Well, Round Bottom

General information

The AbGene 1127 is a very important labware type for the AssayMAP Bravo Platform because it is the only known type of deep-well plate that can be used with the AssayMAP head for bare probe operations. The plate has a low enough profile to enable the probes of the head to reach the bottom of the plate without making contact with the cartridges or the head stripper plate.

Volume recommendations

Number of wells = 96

Maximum volume per well = 1050 μ L

Recommended overage per well = 30 μ L

Resources

For information about...	See...
Product page	https://www.thermofisher.com/order/catalog/product/AB1127
Technical drawing	https://tools.thermofisher.com/content/sfs/manuals/MAN0014419_ABgene_Storage_Plate_96well_1_2mL_QR.pdf

96 AbGene U-Bottom Thermal Insert + AbGene 1127, Square Well, Round Bottom

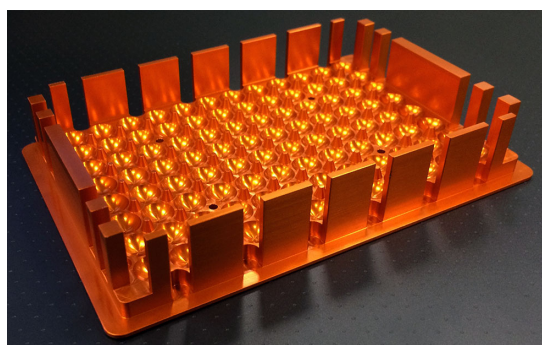


Manufacturer and part number:

- *AbGene plate*: Fisher Scientific AB-1127
- *96 Abgene Deep Well Insert*: Agilent G5498B#127

Name of labware type in the software:

96 AbGene U-Bottom Thermal Insert + AbGene 1127, Square Well, Round Bottom



General information

In applications that require heating of the AbGene 1127 plate, place the AbGene 1127 plate atop the 96 AbGene U-Bottom Thermal Insert to allow efficient heat transfer from the Peltier Thermal Station to the AbGene 1127 plate. The combination of the AbGene 1127 plate stacked on the thermal insert results in a different plate height than the AbGene 1127 plate alone. The combination is designated in the labware selection as **96 AbGene U-Bottom Thermal Insert + AbGene 1127, Square Well, Round Bottom**

Volume recommendations

Number of wells = 96

Maximum volume per well = 1050 μ L

Recommended overage per well = 30 μ L

Resources

For information
about...

See...

AbGene 1127 plate

<https://www.thermofisher.com/order/catalog/product/AB1127>

Thermal Insert

<https://www.agilent.com/cs/library/catalogs/public/5991-0369EN.pdf>

96 Agilent, 2mL Square Deep Well



Manufacturer: Agilent Technologies

Part number: 204353-100

Name of labware type in the software:

96 Agilent, 2mL Square Deep Well

General information

This plate is an option for flow-through and waste collection in the following apps:

- Peptide Cleanup
- Phosphopeptide Enrichment
- Protein Cleanup

Volume recommendations

Number of wells = 96

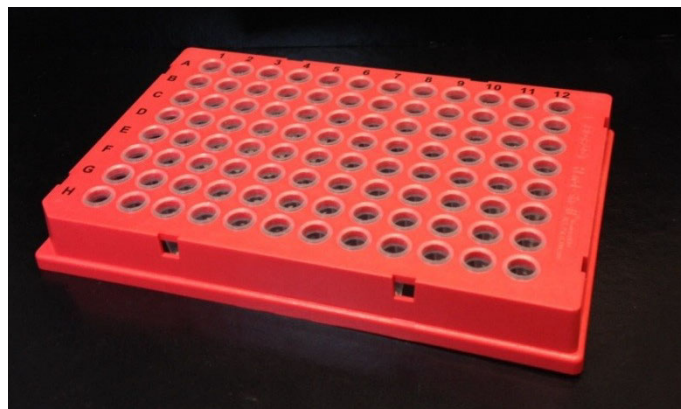
Maximum volume per well = 2200 μ L

Recommended overage per well = Not applicable, not used as a source plate

Resources

For information about...	See...
Ordering	https://www.agilent.com/store/#quickOrder
Data sheet	https://www.agilent.com/cs/library/datasheets/public/ds-cell-analysis-5994-4433en-agilent.pdf

96 Bio-Rad PCR, Hard-Shell, Low-Profile, Full Skirt



Manufacturer: Bio-Rad

Part number: HSP-9611

(See Additional notes.)

Name of labware type in the software:

96 Bio-Rad PCR, Hard-Shell, Low-Profile, Full Skirt

General information

This is one of three types of PCR plates that can be used with the Protein Sample Prep Workbench. Any of the many color varieties of this plate type can be used. The elongated conical geometry of PCR style plates, results in the smallest overage volume requirements.

Volume recommendations

Number of wells = 96

Maximum volume per well = 210 μ L

Recommended overage per well = 10 μ L

Additional notes

Dimensionally equivalent alternatives are the Bio-Rad Hard-Shell, Low-Profile plates, including HSP-9601, 9901, 9621, 9631, 9641, 9661, 9655, 9955, 9665, 9635, 9645, and 9666.

This plate is not compatible with the 25 μ L cartridges at some deck locations. See the relevant application user guide for details.

Note: You can find the application user guides in the Literature Library page of the Protein Sample Prep Workbench.

Resources

For information about...

See...

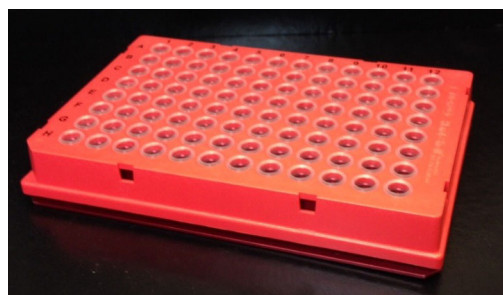
Product page and technical drawing

<https://www.bio-rad.com/en-us/sku/HSP9611-hard-shell-96-well-pcr-plates-low-profile-thin-wall-skirted-red-clear?ID=HSP9611>

2 Labware photo reference

96 Red PCR Insert + Bio-Rad PCR, Hard-Shell, Low-Profile, Full Skirt

96 Red PCR Insert + Bio-Rad PCR, Hard-Shell, Low-Profile, Full Skirt

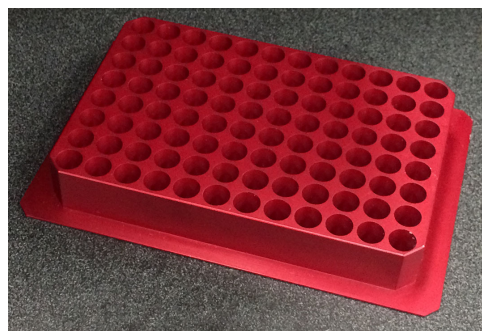


Manufacturer and part number:

- *PCR plate:* Bio-Rad HSP-9611
- *Red PCR Plate Insert:* Agilent G5498B#013

Name of labware type in the software:

96 Red PCR Insert + Bio-Rad PCR, Hard-Shell, Low-Profile, Full Skirt



General information

In applications that require heating the BioRad PCR plate, place the BioRad PCR plate atop the Red PCR Insert to allow efficient heat transfer from the Peltier Thermal Station to the BioRad PCR plate. The combination of the BioRad PCR plate stacked on the thermal insert results in a different plate height than the BioRad PCR plate alone. Therefore, the combination is designated in the labware selection as

96 Red PCR Insert + Bio-Rad PCR, Hard-Shell, Low-Profile, Full Skirt

Volume recommendations

Number of wells = 96

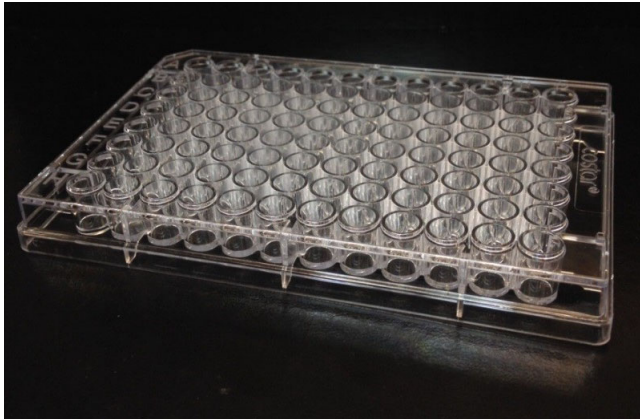
Maximum volume per well = 210 μ L

Recommended overage per well = 10 μ L

Resources

For information about...	See...
Bio-Rad PCR plate	https://www.bio-rad.com/en-us/sku/HSP9611-hard-shell-96-well-pcr-plates-low-profile-thin-wall-skirted-red-clear?ID=HSP9611
Red PCR Plate Insert	https://www.agilent.com/cs/library/catalogs/public/5991-0369EN.pdf

96 Corning 9017, Flat Bottom, Polystyrene, Clear



Manufacturer: Corning

Part number: 9017

Name of labware type in the software: 96 Corning 9017, Flat bottom, polystyrene, clear

General information

This plate is one of three varieties that can be used as a Results Plate and Centrifuge Balance Plate when running the AssayMAP Syringe Test.

Volume recommendations

Number of wells = 96

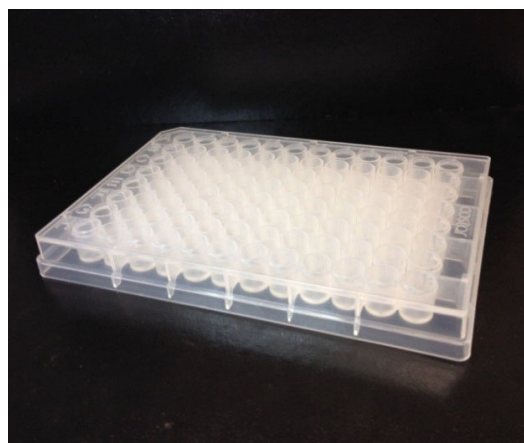
Maximum volume per well = 360 μ L

Recommended overage per well = 50 μ L

Resources

For information about...	See...
Product page	https://ecatalog.corning.com/life-sciences/b2c/US/en/Microplates/Assay-Microplates/96-Well-Microplates/Corning%C2%AE-96-well-Clear-Polystyrene-Microplates/p/9017

96 Costar 3363, PP Conical Bottom



Manufacturer: Corning/Costar

Part number: 3363

Name of labware type in the software:

96 Costar 3363, PP Conical Bottom

General information

The Costar polypropylene conical bottom plate is a good option when attempting to minimize overage volume. The conical bottom forces liquid to drain down to a point, which provides better accessibility for the probes of the AssayMAP Head.

Volume recommendations

Number of wells = 96

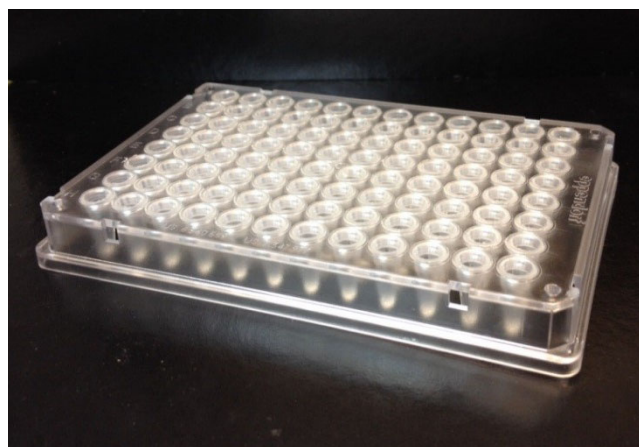
Maximum volume per well = 315 μ L

Recommended overage per well = 20 μ L

Resources

For information about...	See...
Ordering	https://ecatalog.corning.com/life-sciences/b2c/US/en/Microplates/Assay-Microplates/96-Well-Microplates/Corning%C2%AE-96-well-Polypropylene-Microplates/p/3363

96 Eppendorf 30129300, PCR, Full Skirt, PolyPro



Manufacturer: Eppendorf

Part number: 30129300

(See Additional notes.)

Name of labware type in the software:

96 Eppendorf 30129300, PCR, Full Skirt, PolyPro

General information

This is one of three approved types of PCR plates that can be used with the Protein Sample Prep Workbench. While plate movement is not common on AssayMAP applications and utilities, this variety of PCR plate has a rigid skirt, which makes it easier for the Bravo gripper to grip the plate. The elongated conical geometry of PCR style plates, results in the smallest overage volume requirements.

Volume recommendations

Number of wells = 96

Maximum volume per well = 210 μ L

Recommended overage per well = 10 μ L

Additional notes

Any of the Eppendorf 96 well twin.tec skirted plates can be used, such as part numbers 0030129300, 0030129318, 0030129512, 951020401, 951020427, 951020443, 951020460, and 951020486.

The Eppendorf 0030129512 is a low-binding plate that is recommended when low concentrations of proteins or peptides are used to minimize loss due to non-specific binding.

Resources

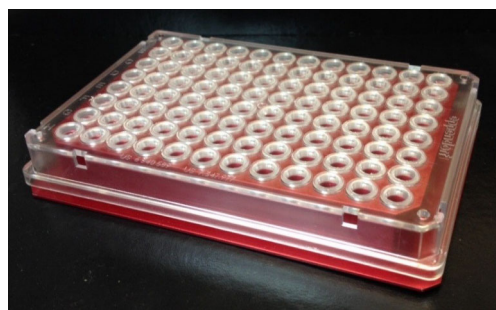
For information about...

See...

Product page and technical drawing

<https://online-shop.eppendorf.us/US-en/Laboratory-Consumables-44512/Plates-44516/Eppendorf-twintec-PCR-Plates-PF-8180.html>

96 Red PCR Insert + Eppendorf 30129300, PCR, Full Skirt



Manufacturer and part number:

- *PCR plate*: Eppendorf 30129300
- *Red PCR Plate Insert*: Agilent G5498B#013

Name of labware type in the software:

96 Red PCR Insert + Eppendorf 30129300, PCR, Full Skirt



General information

In applications that require heating the Eppendorf 30129300 plate, place the Eppendorf 30129300 plate atop the Red PCR Insert to allow efficient heat transfer from the Peltier Thermal Station to the Eppendorf 30129300 plate. The combination of the Eppendorf 30129300 plate stacked on the thermal insert results in a different plate height than the Eppendorf 30129300 plate alone. Therefore, the combination is designated in the labware selection as

96 Red PCR Insert + Eppendorf 30129300, PCR, Full Skirt.

Volume recommendations

Number of wells = 96

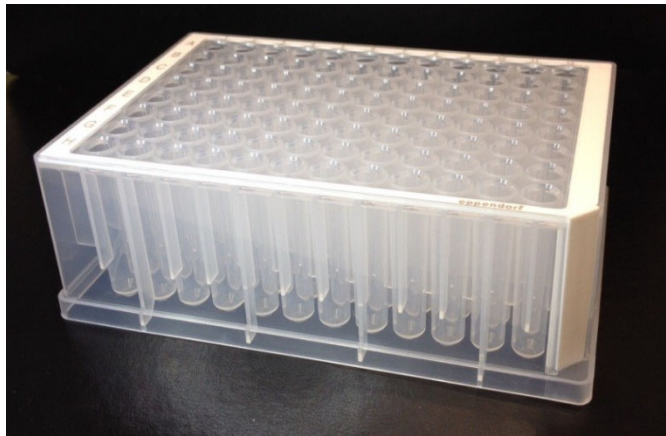
Maximum volume per well = 210 μ L

Recommended overage per well = 10 μ L

Resources

For information about...	See...
Eppendorf 30129300 PCR plate	https://online-shop.eppendorf.us/US-en/Laboratory-Consumables-44512/Plates-44516/Eppendorf-twintec-PCR-Plates-PF-8180.html
Red PCR Plate Insert	https://www.agilent.com/cs/library/catalogs/public/5991-0369EN.pdf

96 Eppendorf 96-1000_U-bottom, Clear PolyPro



Manufacturer: Eppendorf

Part number: Various
(See Additional notes.)

Name of labware type in the software:
96 Eppendorf 96-1000_U-bottom, Clear PolyPro

General information

This type of deep-well plate is an option only in the In-Solution Digestion: Single Plate application and the Reagent Transfer and Reagent Aliquot utilities. Agilent 250- μ L pipette tips are required to use these plates because the AssayMAP Cartridges and the bare probes of the AssayMAP Head cannot reach the well bottoms.

Volume recommendations

Number of wells = 96

Maximum volume per well = 1200 μ L

Recommended overage per well = 25 μ L

Additional notes

This plate comes in many different color options. The following catalog numbers can be used: 951032603, 951032620, 951032662, 951032689, 951032701, 951032727, 951032760, 951032786, 951033006, and 951033103.

Resources

For information about... **See...**

Product page and
technical drawing

<https://online-shop.eppendorf.us/US-en/Laboratory-Consumables-44512/Plates-44516/Eppendorf-Deepwell-Plates-PF-55960.html>

96 Eppendorf 96-500_V-bottom, Clear PolyPro



Manufacturer: Eppendorf

Part number: Various
(See Additional notes.)

Name of labware type in the software: 96
Eppendorf 96-500_V-bottom, Clear PolyPro

General information

This type of deep-well plate is an option only in the In-Solution Digestion: Single Plate application and the Reagent Transfer, Reagent Aliquot, Reformatting, and Serial Dilution utilities. Agilent 250- μ L pipette tips are required to use these plates because the AssayMAP Cartridges and the bare probes of the AssayMAP Head cannot reach the well bottoms.

Volume recommendations

Number of wells = 96

Maximum volume per well = 700 μ L

Recommended overage per well = 25 μ L

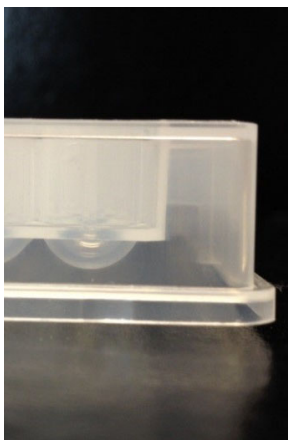
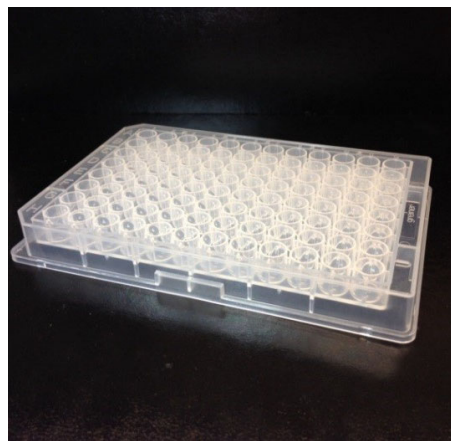
Additional notes

This plate comes in many different color options, which can be used, including the following part numbers: 951031801, 951031828, 951031861, 951031887, 951031909, 951031925, 951031968, 951031984, 951032204, and 951032301.

Resources

For information about...	See...
Product page and technical drawing	https://online-shop.eppendorf.us/US-en/Laboratory-Consumables-44512/Plates-44516/Eppendorf-Deepwell-Plates-PF-55960.html

96 Greiner 650201, U-Bottom Standard, PolyPro



Manufacturer: Greiner Bio-One

Part number: 650201

Name of labware type in the software:

96 Greiner 650201, U-Bottom Standard, PolyPro

General information

The 96 Greiner 650201 is a very common plate type for AssayMAP applications and utilities. It provides a relatively high liquid capacity, and the round bottom helps minimize overage by forcing liquid to drain to the center of the well.

Volume recommendations

Number of Wells = 96

Maximum Volume per Well = 315 μ L

Recommended Overage per Well = 20 μ L

Additional notes

This plate type has two different labware definitions in the software:

1 96 Greiner 650201, U-Bottom Standard, PolyPro

2 [96 Greiner 650201_U-Bottom, Clear PolyPro](#)

Only one of the two naming varieties is available in each Protein Sample Prep Workbench application or utility.

Resources

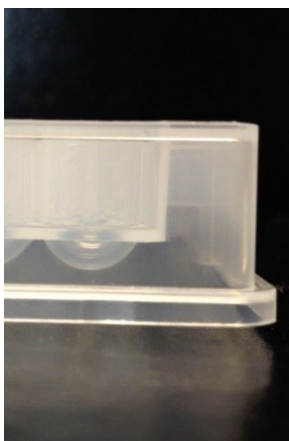
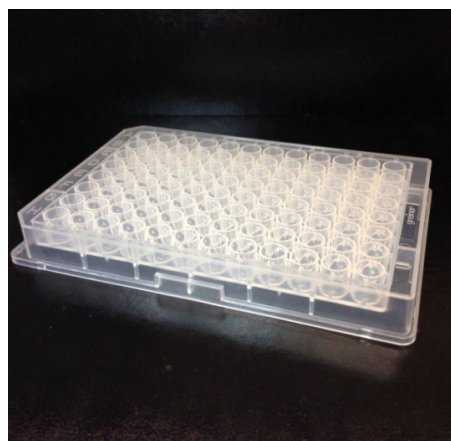
For information about...

See...

Product page and technical drawing

https://shop.gbo.com/en/usa/products/bioscience/microplates/96-well-microplates/96-well-polypropylene-microplates/650201.html?sword_list%5B0%5D=650201&no_cache=1

96 Greiner 650201_U-Bottom, Clear PolyPro



Manufacturer: Greiner Bio-One

Part number: 650201

Name of labware type in the software:

96 Greiner 650201_U-Bottom, Clear PolyPro

General information

This is a very common plate type for AssayMAP applications and utilities. It provides a relatively high liquid capacity, and the round bottom helps minimize overage by forcing liquid to drain to the center of the well.

Volume recommendations

Number of Wells = 96

Maximum Volume per Well = 315 μ L

Recommended Overage per Well = 20 μ L

Additional notes

This plate type has two different labware definitions in the software:

1 96 Greiner 650201_U-Bottom, Clear PolyPro

2 [96 Greiner 650201, U-Bottom Standard, PolyPro](#)

Only one of the two naming varieties is available in each Protein Sample Prep Workbench application or utility.

Resources

For information about...

See...

Product page and technical drawing

https://shop.gbo.com/en/usa/products/bioscience/microplates/96-well-microplates/96-well-polypropylene-microplates/650201.html?sword_list%5B0%5D=650201&no_cache=1

96 Greiner U-Bottom Thermal Insert + Greiner 650201_U-Bottom, Clear PolyPro

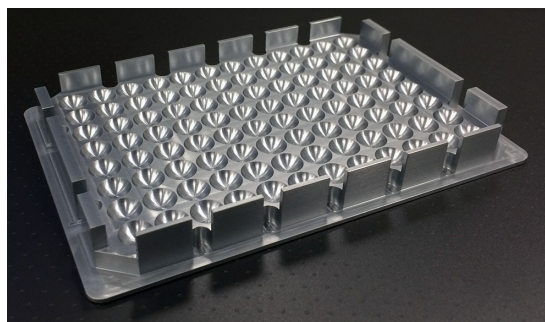


Manufacturer and part number:

- *Greiner plate:* Greiner Bio-One 650201
- *96 Greiner U- or V-Bottom Insert:* Agilent G5498B#126

Name of labware type in the software:

96 Greiner U-Bottom Thermal Insert +
Greiner 650201_U-Bottom, Clear PolyPro



General information

In applications that require heating the Greiner 650201 plate, place the plate atop the 96 Greiner U-Bottom Thermal Insert to allow efficient heat transfer from the Peltier Thermal Station to the Greiner 650201 plate. The combination of the Greiner 650201 plate stacked on the thermal insert results in a different plate height than the Greiner 650201 plate alone. Therefore, the combination is designated in the labware selection as **96 Greiner U-Bottom Thermal Insert + Greiner 650201_U-Bottom, Clear PolyPro**

Volume recommendations

Number of Wells = 96

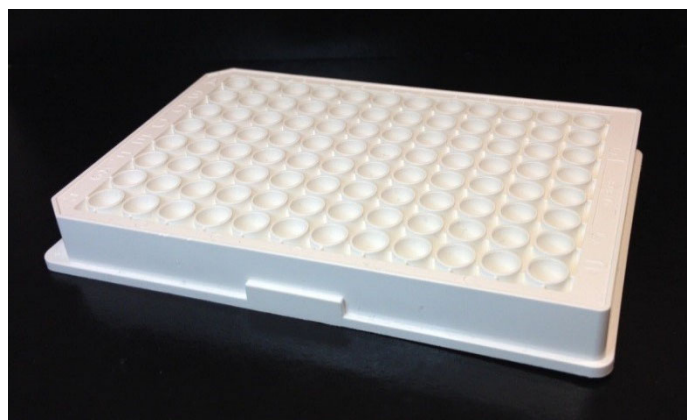
Maximum Volume per Well = 315 µL

Recommended Overage per Well = 20 µL

Resources

For information about...	See...
Greiner plate	https://shop.gbo.com/en/usa/products/bioscience/microplates/96-well-microplates/96-well-polypropylene-microplates/650201.html?sword_list%5B0%5D=650201&no_cache=1
Thermal Insert	https://www.agilent.com/cs/library/catalogs/public/5991-0369EN.pdf

96 Greiner 650207_U-Bottom, White PolyPro



Manufacturer: Greiner Bio-One

Part number: 650207

Name of labware type in the software:
96 Greiner 650207_U-Bottom, White PolyPro

General information

This is a very common plate type for AssayMAP applications and utilities. It provides a relatively high liquid capacity, and the round bottom helps minimize overage by forcing liquid to drain to the center of the well. The opaque white coloration also provides additional light protection for sensitive reagents.

Volume recommendations

Number of wells = 96

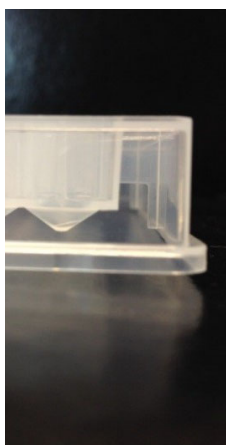
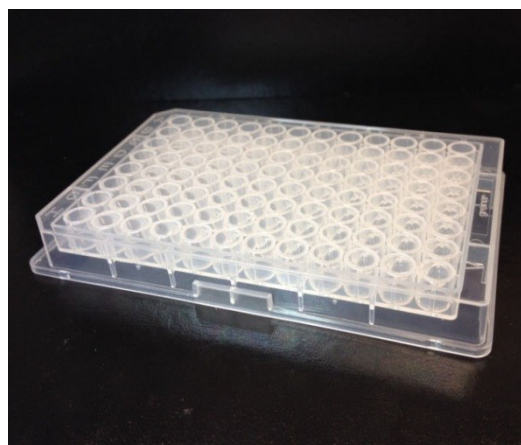
Maximum volume per well = 315 μ L

Recommended overage per well = 20 μ L

Resources

For information about...	See...
Product page and technical drawing	https://shop.gbo.com/en/usa/products/bioscience/microplates/us-made-microplates/0110_0040_0200_0030/650207.html?sword_list%5B0%5D=650207&no_cache=1

96 Greiner 651201_V-Bottom, Clear PolyPro



Manufacturer: Greiner Bio-One

Part number: 651201

Name of labware type in the software:
96 Greiner 651201_V-Bottom, Clear PolyPro

General information

The Greiner polypropylene V-bottom plate is a good option when attempting to minimize overage volume. The conical bottom forces liquid to drain down to a point, providing good accessibility for the bare probes of the AssayMAP Head or for the cartridges.

Volume recommendations

Number of wells = 96

Maximum volume per well = 315 μ L

Recommended overage per well = 20 μ L

Resources

For information about...

See...

Product page and technical drawing

https://shop.gbo.com/en/usa/products/bioscience/microplates/96-well-microplates/96-well-polypropylene-microplates/651201.html?sword_list%5B0%5D=651201&no_cache=1

96 Greiner V-Bottom Thermal Insert + Greiner 651201, V-Bottom, Clear PolyPro

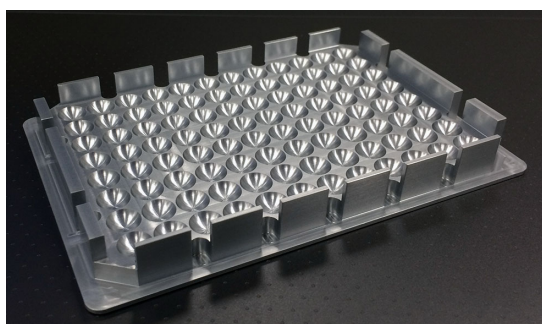


Manufacturer and part number:

- Greiner plate: Greiner Bio-One 651201
- 96 Greiner U- or V-Bottom Insert: Agilent G5498B#126

Name of labware type in the software:

96 Greiner V-Bottom Thermal Insert +
Greiner 651201, V-Bottom, Clear PolyPro



General information

In applications that require heating the Greiner 651201 plate, place the plate atop the 96 Greiner V-Bottom Thermal Insert to allow efficient heat transfer from the Peltier Thermal Station to the Greiner 651201 plate. The combination of the Greiner 651201 plate stacked on the thermal insert results in a different plate height than the Greiner 651201 plate alone. Therefore, the combination is designated in the labware selection as **96 Greiner V-Bottom Thermal Insert + Greiner 651201, V-Bottom, Clear PolyPro**.

Volume recommendations

Number of wells = 96

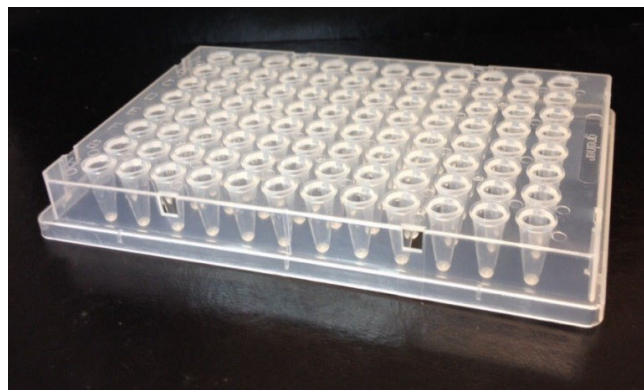
Maximum volume per well = 315 μ L

Recommended overage per well = 20 μ L

Resources

For information about...	See...
Greiner plate	https://shop.gbo.com/en/usa/products/bioscience/microplates/96-well-microplates/96-well-polypropylene-microplates/651201.html?sword_list%5B0%5D=651201&no_cache=1
Thermal Insert	https://www.agilent.com/cs/library/catalogs/public/5991-0369EN.pdf

96 Greiner 652270, PCR, Full Skirt, PolyPro



Manufacturer: Greiner Bio-One

Part number: 652270

Name of labware type in the software:
96 Greiner 652270, PCR, Full Skirt, PolyPro

General information

This is one of three approved types of PCR plates that can be used with the Protein Sample Prep Workbench. The elongated conical geometry of PCR style plates, results in the smallest overage volume requirements.

Volume recommendations

Number of wells = 96

Maximum volume per well = 210 μ L

Recommended overage per well = 10 μ L

Additional notes

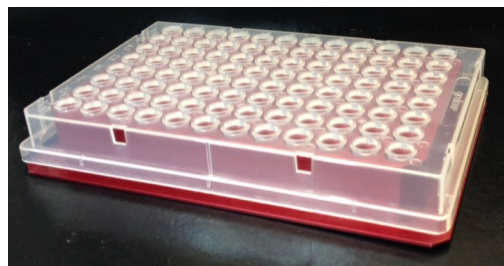
This plate is not compatible with the 25 μ L cartridges at some deck locations. See the relevant application user guide for details.

Note: You can find the application user guides in the Literature Library page of the Protein Sample Prep Workbench.

Resources

For information about...	See...
Product page and technical drawing	https://shop.gbo.com/en/usa/products/bioscience/molecular-biology/pcr-microplates/652270.html?sword_list%5B0%5D=652270&no_cache=1

96 Red PCR Insert + Greiner 652270, PCR, Full Skirt, PolyPro

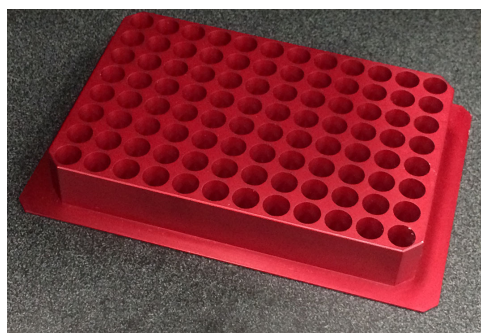


Manufacturer and part number:

- *PCR plate:* Greiner Bio-One 652270
- *Red PCR Plate Insert:* Agilent G5498B#013

Name of labware type in the software:

96 Red PCR Insert + Greiner 652270, PCR, Full Skirt, PolyPro



General information

In applications that require heating the Greiner 652270 plate, place the Greiner 652270 plate atop the Red PCR Insert to allow efficient heat transfer from the Peltier Thermal Station to the Greiner 652270 plate. The combination of the Greiner 652270 plate stacked on the thermal insert results in a different plate height than the Greiner 652270 plate alone. Therefore, the combination is designated in the labware selection as **96 Red PCR Insert + Greiner 652270, PCR, Full Skirt, PolyPro**.

Volume recommendations

Number of wells = 96

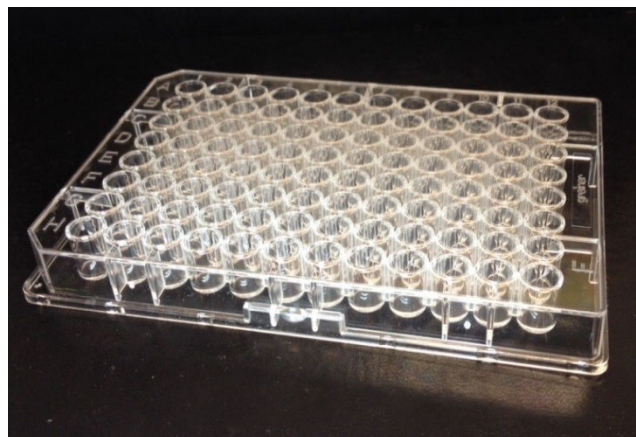
Maximum volume per well = 210 µL

Recommended overage per well = 10 µL

Resources

For information about...	See...
Greiner plate	https://shop.gbo.com/en/usa/products/bioscience/molecular-biology/pcr-microplates/652270.html?sword_list%5B0%5D=652270&no_cache=1
Red PCR Plate Insert	https://www.agilent.com/cs/library/catalogs/public/5991-0369EN.pdf

96 Greiner 655101 PS Clr Rnd Well Flat Btm



Manufacturer: Greiner Bio-One

Part number: 655101

Name of labware type in the software:
96 Greiner 655101 PS Clr Rnd Well Flat Btm

General information

This plate is one of three varieties that can be used as a Results Plate and Centrifuge Balance Plate when running an AssayMAP Syringe Test.

Volume recommendations

Number of wells = 96

Maximum volume per well = 360 μ L

Recommended overage per well = 50 μ L

Additional notes

Two of these plates are included with the AssayMAP Syringe Test kit.

Resources

For information about...	See...
Product page and technical drawing	https://shop.gbo.com/en/usa/products/bioscience/microplates/96-well-microplates/96-well-microplates-clear/655101.html?sword_list%5B0%5D=655101&no_cache=1

96 Greiner 675801, Half Area, Flat Bottom, UV Star



Manufacturer: Greiner Bio-One

Part Number: 675801

Name of labware type in the software:

96 Greiner 675801, Half Area, Flat Bottom, UV Star

General information

This plate is generally used for the purification and normalization applications, when protein concentration will need to be determined using protein A280. The flat bottom on this Greiner plate results in a very high overage volume, which is an important consideration when transitioning from this labware into another type.

Volume recommendations

Number of wells = 96

Maximum volume per well = 190 μ L

Recommended overage per well = 25 μ L

Resources

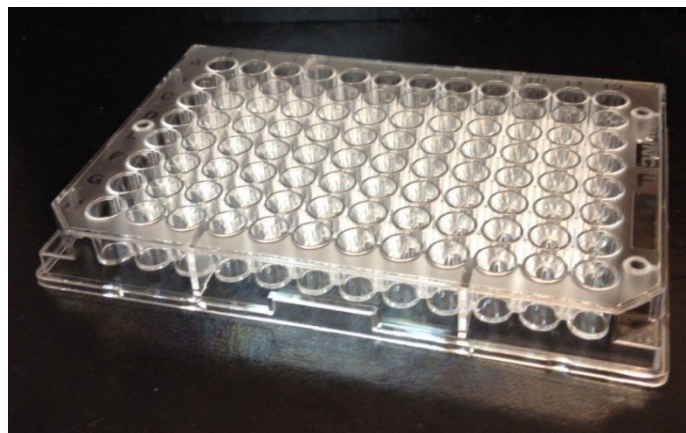
For information about...

See...

Product page and technical drawing

https://shop.gbo.com/en/usa/products/bioscience/microplates/uv-star-microplates/96-well-uv-star-microplate/675801.html?sword_list%5B0%5D=675801&no_cache=1

96 Nunc 269620, Flat Bottom, Polystyrene



Manufacturer: Nunc (Thermo-Fisher)

Part number: 269620

Name of labware type in the software:
96 Nunc 269620, Flat Bottom, Polystyrene

General information

This plate is one of three varieties that can be used as a Results Plate and Centrifuge Balance Plate when running the AssayMAP Syringe Test utility.

Volume recommendations

Number of wells = 96

Maximum volume per well = 360 μ L

Recommended overage per well = 50 μ L

Resources

For information about...

See...

Product page and
technical drawing

<https://www.thermofisher.com/order/catalog/product/269620?SID=srch-hj-269620>

96 Thermo Matrix 3732, V-bottom, 0.75ml Storage Tubes



Manufacturer: ThermoFisher Scientific

Part number: 3732

Name of labware type in the software:

96 Thermo Matrix 3732, V-bottom, 0.75ml
Storage Tubes

General information

This storage tube rack is one of three varieties that can be used for the following:

- Diluent reservoir and Serial Dilution plate for the Serial Dilution utility
- Source and Destination plates for the Reformatting utility

Volume recommendations

Number of wells = 96

Maximum volume per tube = 0.75 mL

Recommended overage per well = 20 μ L

Resources

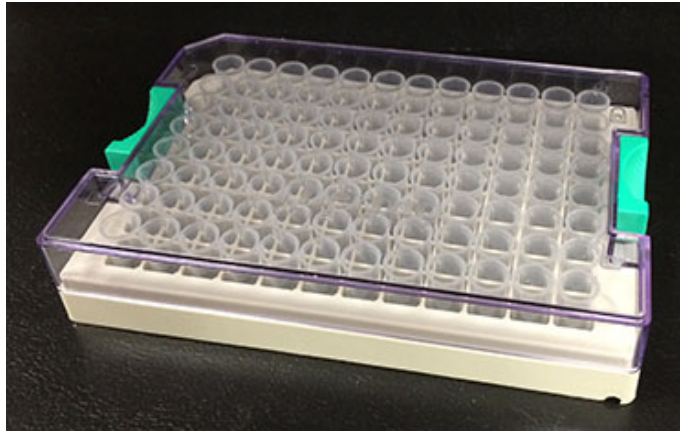
For information about...

See...

Product page

<https://www.thermofisher.com/order/catalog/product/3732?SID=srch-srp-3732>

96 Thermo Matrix 3735, V-bottom, 500µl Storage Tubes



Manufacturer: ThermoFisher Scientific

Part number: 3735

Name of labware type in the software:
96 Thermo Matrix 3735, V-bottom, 500µl
Storage Tubes

General information

This storage tube rack is one of three varieties that can be used for the following:

- Diluent reservoir and Serial Dilution plate for the Serial Dilution utility
- Source and Destination plates for the Reformatting utility

Volume recommendations

Number of wells = 96

Maximum volume per tube = 500 µL

Recommended Overage per Well = 20 µL

Resources

For information about...	See...
Product page	https://www.thermofisher.com/order/catalog/product/3735?SID=srch-srp-3735

96 Thermo Matrix 3744, V-bottom, 500µl ScrewTop Storage Tubes



Manufacturer: ThermoFisher Scientific

Part number: 3744

Name of labware type in the software:

96 Thermo Matrix 3744, V-bottom, 500µl
ScrewTop Storage Tubes

General information

This storage tube rack is one of three varieties that can be used for the following:

- Diluent reservoir and Serial Dilution plate for the Serial Dilution utility
- Source and Destination plates for the Reformatting utility

Volume recommendations

Number of wells = 96

Maximum volume per tube = 500 µL

Recommended Overage per Well = 20 µL

Resources

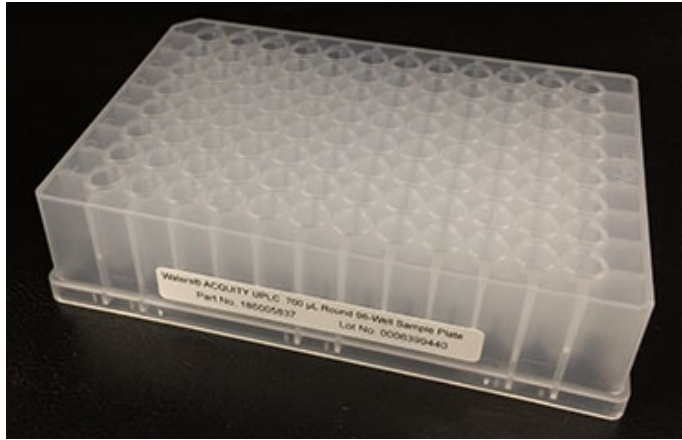
For information about...

See...

Product page

<https://www.thermofisher.com/order/catalog/product/3744?SID=srch-srp-3744>

96 Waters 186005837, Clear PolyPro



Manufacturer: Waters

Part number: 186005837

Name of labware type in the software:
96 Waters 186005837, Clear PolyPro

General information

This plate can be used for the following:

- Reagent Aliquot, Reagent Transfer, Reformatting, and Serial Dilution utilities
- In-Solution Digestion: Single Plate application

Volume recommendations

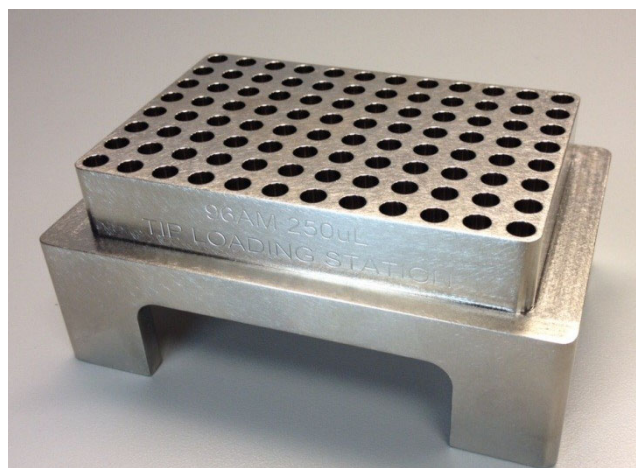
Number of wells = 96

Maximum volume per tube = 500 µL

Resources

For information about...	See...
Product page	https://www.waters.com/nextgen/us/en/shop/vials-containers--collection-plates/186005837-96-well-sample-collection-plate-700--l-round-well-25-pk.html

96AM 250 uL Tip Loading Station



Manufacturer: Agilent Technologies

Part number: G5409-20025

Name of labware type in the software:
96AM 250 uL Tip Loading Station

General information

The Agilent 96AM Cartridge & Tip Seating Station accessory is included with all AssayMAP Bravo purchases. This accessory is required when seating pipette tips on the probes of the AssayMAP Bravo head.

Volume recommendations

Number of wells = 96

Maximum volume per well = Not applicable

Recommended overage per well = Not applicable

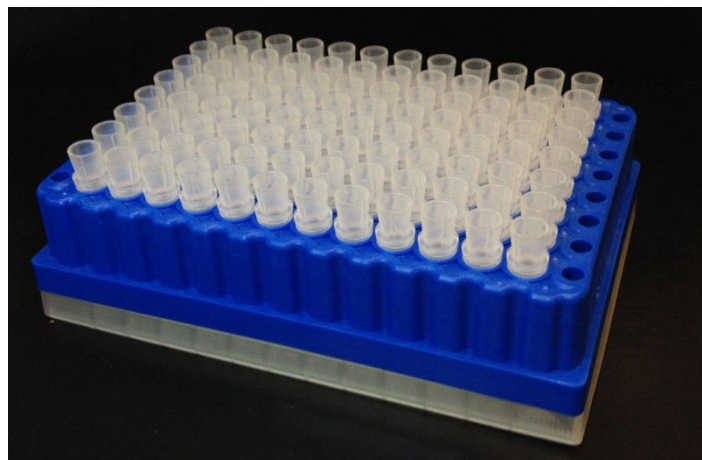
Additional notes

This accessory has two labware names in the software:

- 1 96AM 250 uL Tip Loading Station, which is used for seating tips on the AssayMAP Bravo head
- 2 [96AM Cartridge Seating Station](#), which is used for seating cartridges on the AssayMAP Bravo head

The engraving, 96AM 250uL TIP LOADING STATION, appears on the side of the accessory.

96AM Cartridge Rack and Receiver Plate



Manufacturer: Agilent Technologies

Part number: See the Agilent [AssayMAP Cartridge Selection Guide](#).

Name of labware type in the software:
96AM Cartridge Rack and Receiver Plate

General information

The AssayMAP Cartridges are shipped in the 96AM Cartridge Rack and Receiver Plate. The labware consists of two distinct pieces: the blue cartridge rack and the white receiver plate. The cartridges are transferred from this labware into the [96AM Cartridge Seating Station](#) before use on the AssayMAP Bravo Platform.

Volume recommendations

Number of wells = 96

Maximum volume per well = Not applicable

Recommended overage per well = Not applicable

Additional notes

Partial plate operations with the head positioned at an offset to the deck location are possible with this piece of labware, for example, using the Cartridge Transfer utility.

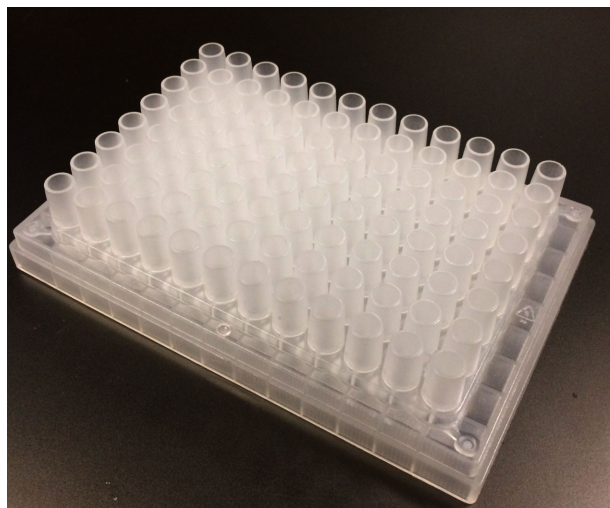
One known issue is that the AssayMAP Bravo head might pick up the cartridge rack unintentionally while attempting to seat a full rack of 96 cartridges. When using a full cartridge rack, pay close attention to the AssayMAP Bravo head during a run of the Cartridge Transfer utility.

Resources

[For information about...](#) [See...](#)

Product details [Agilent AssayMAP Cartridge Selection Guide](#)

96AM Receiver Plate



Manufacturer: Agilent Technologies

Part number: Not applicable

Name of labware type in the software: 96AM Receiver Plate

General information

The 96AM Receiver Plate is the bottom piece of the [96AM Cartridge Rack and Receiver Plate](#). The cartridge rack stacks on the receiver plate during shipment. The cartridges are transferred from the 96AM Cartridge Rack and Receiver plate into the [96AM Cartridge Seating Station](#) for automation tasks on the AssayMAP Bravo Platform. If the cartridge rack is used during automation tasks, the receiver plate is required for support.

Volume recommendations

Number of wells = 96

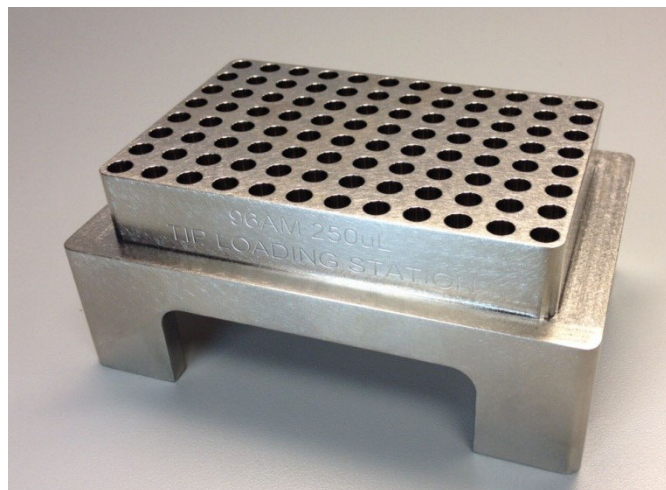
Maximum volume per well = 200 μ L when a cartridge is present

Recommended overage per well = Not applicable

Additional notes

The wells of the receiver plate can be filled with 200 μ L of liquid and used as a temporary storage location for used cartridges, for example, after running the immobilization protocol and before using the cartridges in the affinity purification protocol. When the cartridges are put in the blue rack in a position where the well below it has 200 μ L of liquid, the cartridge tips will be submerged in the liquid and thereby avoid drying out. Cartridges should never be allowed to dry out.

96AM Cartridge Seating Station



Manufacturer: Agilent Technologies

Part number: G5409-20025

Name of labware type in the software:
96AM Cartridge Seating Station

General information

The Agilent 96AM Cartridge & Tip Seating Station accessory is included with all AssayMAP Bravo purchases. This accessory is the “labware” used for seating the cartridges on the probes of the AssayMAP Bravo head.

Volume recommendations

Number of wells = 96

Maximum volume per well = Not applicable

Recommended overage per well = Not applicable

Additional notes

This accessory has two labware type names in the software:

- 1 96AM Cartridge Seating Station, which is used for seating cartridges on the AssayMAP Bravo head
- 2 [96AM 250 uL Tip Loading Station](#), which is used for seating tips on the AssayMAP Bravo head

The engraving, 96AM 250uL TIP LOADING STATION, appears on the side of the accessory.

96AM Tip Wash Station



Manufacturer: Agilent Technologies

Part number:

G5498B#90

96 Channel Wash Station (G5409-60125)

or

G5498B#057

96AM Wash Station (11961-311)

Name of labware type in the software:

96AM Tip Wash Station

General information

The wash station is installed on deck location 1 of the AssayMAP Bravo Platform. The wash station is used for internally and externally washing syringes, cartridges, and pipette tips. The wash station also serves as a waste collection location.

The AssayMAP Bravo Platform requires one of the following Agilent wash stations:

- **96 Channel Wash Station G5498B#90.** Equipped with 96 wide-bore chimneys (part number G5409-68004) and a bracketed platepad to secure it.
- **96AM Wash Station G5498B#057.** Originally equipped with 96 standard-bore chimneys (part number 15586-002) This model also has a bracketed platepad to secure it. This wash station may be retrofitted with the wide-bore chimneys.

Volume recommendations

Number of wells = 96

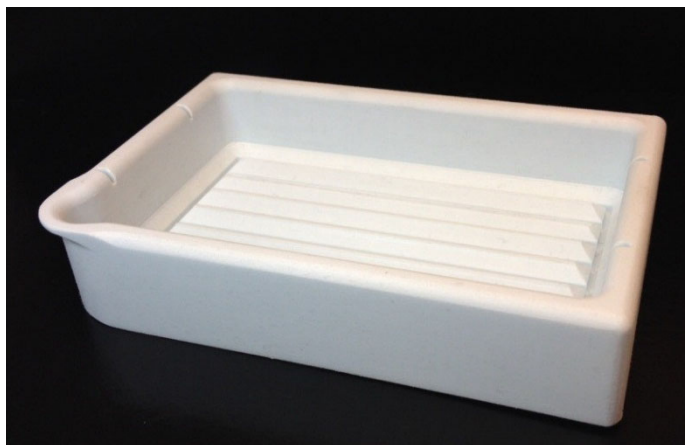
Maximum volume per well = Not applicable

Recommended overage per well = Not applicable

Resources

For information about...	See...
Wash station chimneys	Agilent 96 Channel Wash Station Maintenance Guide

96 V11 Manual Fill Reservoir



Manufacturer: Agilent Technologies

Part Number: G5498B#049
(Internal part number: 08105-001)

Name of labware type in the software:
96 V11 Manual Fill Reservoir,
96 V11 Reservoir (Manual Fill)

General information

The 96 V11 Manual Fill Reservoir is an open-bath reservoir that can be used for full plate runs of 96 samples or cartridges.

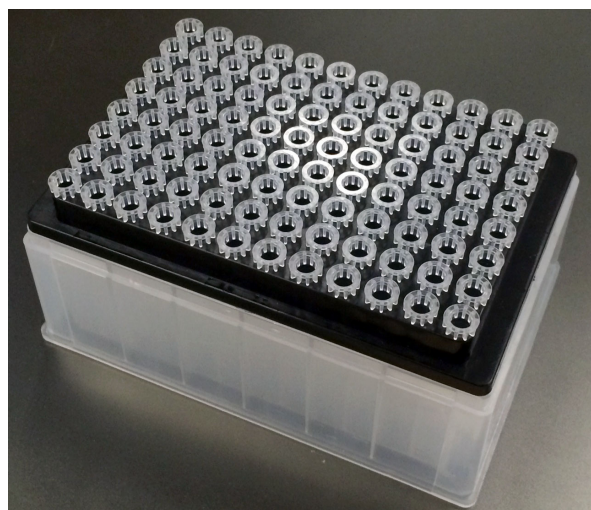
Volume recommendations

Number of wells = 1

Maximum volume per well = 150 mL

Recommended overage per well = 35 mL

96 V11 LT 250 Tip Box 19477.002



Manufacturer: Agilent Technologies

Part number: 19477-002

Name of labware type in the software:
96 V11 LT 250 Tip Box

General information

The pipette tips are transferred from the tip box into the [96AM 250 uL Tip Loading Station](#) before use in an AssayMAP application or utility.

Volume recommendations

Number of wells = 96

Maximum volume per well = 140 μ L can be pulled into a mounted tip before the excess fluid overflows into the syringe.

Recommended overage per well = Not applicable

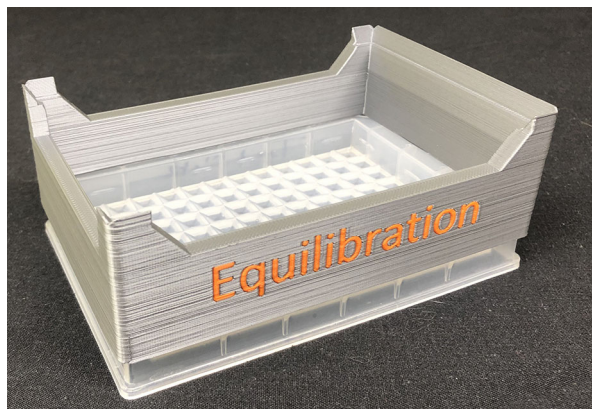
Additional notes

The [96AM 250 uL Tip Loading Station](#) is required to seat the pipette tips on the Bravo 96AM Head.

Resources

For information about...	See...
Ordering	https://www.chem.agilent.com/store/

Equilibration Holder + Reservoir, Seahorse 201254-100

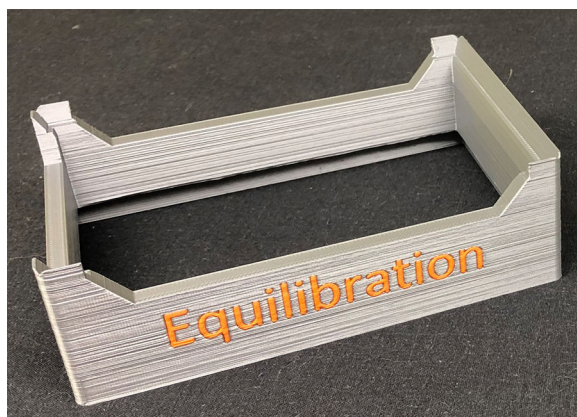


Manufacturer and part number:

- *Equilibration Holder*: Evosep EV1142
- *Reservoir*: One of the following:
 - Agilent 201254-100
 - Agilent 201280-100
 - Agilent 201282-100

Name of labware type in the software:

Equilibration Holder + Reservoir, Seahorse 201254-100



General information

The Equilibration Holder with reservoir is required for the Agilent Evosep - 96AM, Semi-Automated, Full Plate Workflow. The following reservoir types are options for this configuration:

- [Reservoir, Seahorse 201254-100, PP, no walls, pyramid bottoms](#)
- [12 Column, Low Profile Reservoir, Natural PP](#)
- [8 Row, Low Profile Reservoir, Natural PP](#)

You manually place the reservoir in the Equilibration Holder when setting up the protocol run. The reservoir fit is snug, so you must ensure that it is well seated within the holder. During the run, the AssayMAP Bravo moves the Evotip carrier with Evotips to and from the top of the reservoir that is seated in the Equilibration Holder.

Resources

For information about... See...

Ordering

<https://www.evosep.com>

Evotip Holder with Evotips



Manufacturer: Evosep

Part number:

- *Evotip Holder:* EV1142
- *Evotips in tip box:* EV2001

Name of labware type in the software:

Evotip Holder with Evotips

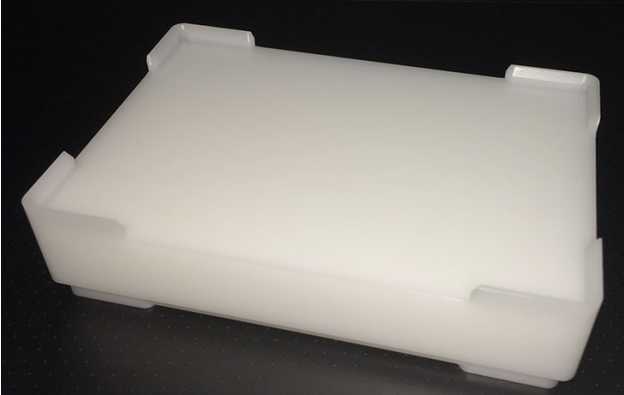
General information

The Evotip Holder with Evotips is required for the Agilent Evosep - 96AM, Semi-Automated, Full Plate Workflow. The Evotip Holder is designed to hold an Evotip carrier (yellow) with 96 Evotips. You manually position the carrier with Evotips in the holder when setting up the protocol run. During the run, the AssayMAP Bravo moves the carrier with Evotips to and from the top of the reservoir that is seated in the Equilibration Holder ("[Equilibration Holder + Reservoir, Seahorse 201254-100](#)" on [page 61](#)).

Resources

For information about...	See...
Ordering	https://www.evosep.com

Bravo Plate Riser



Vendor: Agilent Technologies

Part number: G5498B#061

Name of labware type in the software:
Plate Riser

General information

The Bravo plate riser is required for certain AssayMAP applications and utilities where pipette tips are used for the liquid handling. The 2.84-cm tall riser lifts the labware to a height to the Bravo 96AM Head to be offset when using pipette tips. For example, the Normalization, Serial Dilution, and Reformatting utilities require a pair of plate risers. One riser is placed on the platepad under the source plate and one riser is placed on the platepad under the destination plate.

Universal Lid PS Black



Vendor: Agilent Technologies

Part number: 200858-100

Name of labware type in the software:
Plate Lidded

General information

Lidded plates, not the lid by itself, are allowed for the In-Solution Digestion: Single-Plate application. If you plan to use a lidded plate, ensure that you use this universal microplate lid, which is the only approved lid type.

Additional notes

Several different vendors offer the exact same lid under different part numbers, for example, Phenix Research ML-5009 or ML-5010, E&K Scientific EK-2079, EK-2080, or EK-2081.

Three color varieties are available for this lid: opaque black, opaque white, and transparent. All three colors are acceptable for use with the In-Solution Digestion: Single-Plate application.

Resources

For information about...	See...
Ordering	https://www.chem.agilent.com/store/

In This Book

This guide provides a labware reference for the AssayMAP Protein Sample Prep Workbench, including

- Labware options and default settings
- Labware photo reference

www.agilent.com

© Agilent Technologies, Inc. 2022

October 2022



D0012628 Revision A

