

Filtration Microplates

For protein purification, sample preparation, and DNA/RNA/oligo isolation

Agilent filtration microplates incorporate most media types, such as glass fiber, polyethylene, polypropylene, nylon, polyethersulfone (PES), and polyvinylidene fluoride (PVDF). They are available in 4 well configurations (24, 48, 96 and 384) with well volumes ranging from 140 μ l to 15 ml and are suitable for vacuum and positive pressure manifolds, as well as centrifugation. All Agilent filter plates can be paired with a corresponding collection plate.

To add items to your “My Favorites”* list at the [Agilent online store](#), simply click the part number. Then, enter the quantities for the products you need.



The filtration microplates listed below are organized by pore size and well volume.

Description	Pore Size	Well Volume (ml)	Part Number
96-well polypropylene with polyethersulfone membrane, short drip, pack of 25, recommended receiver plate 204600-100 (204601-100 irradiated)	10 KDa	0.4	201013-100
96-well polypropylene with polyethersulfone membrane, short drip, pack of 25, recommended receiver plate 204600-100 (204601-100 irradiated)	30 KDa	0.4	201015-100
96-well polypropylene with polyethersulfone membrane, short drip, pack of 25, recommended receiver plate 204600-100 (204601-100 irradiated)	100 KDa	0.4	201596-100
96-well polypropylene with polyethersulfone membrane, short drip, pack of 25, recommended receiver plate 204600-100 (204601-100 irradiated)	300 KDa	0.4	201598-100
24-well polypropylene with polyvinylidene fluoride membrane, long drip, pack of 25, recommended receiver plate 201272-100 (204023-100 irradiated)	0.2 μ m	10	204465-100
24-well polypropylene with polyethersulfone membrane, long drip, pack of 25, recommended receiver plate 201272-100 (204023-100 irradiated)	0.2 μ m	10	204469-100
96-well polypropylene with polyethersulfone membrane, long drip, pack of 25, recommended receiver plate 201240-100 (204353-100 irradiated)	0.2 μ m	2	204510-100

Description	Pore Size	Well Volume (ml)	Part Number
96-well polypropylene with polyvinylidene fluoride membrane, long drip, pack of 25, recommended receiver plate 201276-100 (204355-100 irradiated)	0.2 µm	0.8	203982-100
96-well polypropylene with polyethersulfone membrane, long drip, pack of 25, recommended receiver plate 204600-100 (204601-100 irradiated)	0.2 µm	0.4	203940-100
96-well polypropylene with polyvinylidene fluoride membrane, long drip, pack of 25, recommended receiver plate 204600-100 (204601-100 irradiated)	0.2 µm	0.4	203980-100
24-well polypropylene with polyvinylidene fluoride membrane, long drip, pack of 25, recommended receiver plate 201272-100 (204023-100 irradiated)	0.45 µm	10	204463-100
24-well polypropylene with polyethersulfone membrane, long drip, pack of 25, recommended receiver plate 201272-100 (204023-100 irradiated)	0.45 µm	10	204467-100
96-well polypropylene with polypropylene membrane, long drip, pack of 25, recommended receiver plate 201240-100 (204353-100 irradiated)	0.45 µm	2	201009-100
96-well polypropylene with polyethersulfone membrane, long drip, pack of 25, recommended receiver plate 201240-100 (204353-100 irradiated)	0.45 µm	2	201011-100
96-well polypropylene with polyvinylidene fluoride membrane, long drip, pack of 25, recommended receiver plate 201240-100 (204353-100 irradiated)	0.45 µm	2	201731-100
96-well polypropylene with polyvinylidene fluoride membrane, long drip, pack of 25, recommended receiver plate 201276-100 (204355-100 irradiated)	0.45 µm	0.8	200931-100
96-well polypropylene with polypropylene membrane, long drip, pack of 25, recommended receiver plate 201276-100 (204355-100 irradiated)	0.45 µm	0.8	200933-100
96-well polypropylene with polypropylene membrane, short drip, pack of 25, recommended receiver plate 201276-100 (204355-100 irradiated)	0.45 µm	0.8	200961-100
96-well polypropylene with polyvinylidene fluoride membrane, short drip, pack of 25, recommended receiver plate 201276-100 (204355-100 irradiated)	0.45 µm	0.8	200959-100
96-well polypropylene with polyvinylidene fluoride membrane, long drip, pack of 25, recommended receiver plate 204600-100 (204601-100 irradiated)	0.45 µm	0.4	200943-100
96-well polypropylene with polyvinylidene fluoride membrane, short drip, pack of 25, recommended receiver plate 204600-100 (204601-100 irradiated)	0.45 µm	0.4	200971-100
96-well polypropylene with polypropylene membrane, long drip, pack of 25, recommended receiver plate 204600-100 (204601-100 irradiated)	0.45 µm	0.4	200945-100
96-well polypropylene with polypropylene membrane, short drip, pack of 25, recommended receiver plate 204600-100 (204601-100 irradiated)	0.45 µm	0.4	200973-100
96-well polypropylene with polyvinylidene fluoride membrane, long drip, pack of 50, recommended receiver plate 203942-100 (204602-100 irradiated)	0.45 µm	0.3	200981-100
96-well polypropylene with polyvinylidene fluoride membrane, short drip, pack of 50, recommended receiver plate 203942-100 (204602-100 irradiated)	0.45 µm	0.3	200995-100
96-well polypropylene with polypropylene membrane, long drip, pack of 50, recommended receiver plate 203942-100 (204602-100 irradiated)	0.45 µm	0.3	200983-100
96-well polypropylene with polypropylene membrane, short drip, pack of 50, recommended receiver plate 203942-100 (204602-100 irradiated)	0.45 µm	0.3	200997-100
96-well polypropylene with nylon membrane, short drip, pack of 50, recommended receiver plate 203942-100 (204602-100 irradiated)	0.45 µm	0.3	201738-100
96-well polypropylene with glass fiber membrane, long drip, pack of 25, recommended receiver plate 201240-100	0.7 µm	2	201007-100
96-well polypropylene for protein precipitation, long drip, pack of 5, recommended receiver plate 201240-100	0.7 µm	1.8	203845-100
96-well polypropylene for protein precipitation, long drip, includes a receiver plate 201240-100 and lid 200856-100, pack of 1 kit	0.7 µm	1.8	203846-100
96-well polypropylene for protein precipitation, long drip, pack of 1, recommended receiver plate 201240-100	0.7 µm	1.8	203847-100
96-well polypropylene half skirt with 4 layers of fiber membrane, long drip with turbine, pack of 25, recommended receiver plate 201276-100	0.7 µm	1	201719-100
96-well polypropylene half skirt with polyethylene frit plus 2 layers of glass fiber membrane, long drip with turbine, pack of 25, recommended receiver plate 201276-100	0.7 µm	1	201720-100
96-well polypropylene with glass fiber membrane, long drip, pack of 25, recommended receiver plate 201276-100	0.7 µm	0.8	200937-100

Description	Pore Size	Well Volume (ml)	Part Number
96-well polypropylene with glass fiber membrane, short drip, pack of 25, recommended receiver plate 201276-100	0.7 µm	0.8	200965-100
96-well polypropylene with 4X glass fiber membranes, long drip, pack of 25, recommended receiver plate 201276-100	0.7 µm	0.8	201537-100
96-well polypropylene with 65 µm polyethylene frit plus glass fiber membrane, long drip, pack of 25, recommended receiver plate 201276-100	0.7µm	0.8	201019-100
96-well polypropylene with glass fiber membrane, long drip, pack of 25, recommended receiver plate 201276-100	0.7 µm	0.4	200951-100
96-well polypropylene with glass fiber membrane, short drip, pack of 25, recommended receiver plate 204600-100	0.7 µm	0.4	200977-100
96-well polypropylene with glass fiber membrane, long drip, pack of 50, recommended receiver plate 203942-100	0.7 µm	0.3	200987-100
96-well polypropylene with glass fiber membrane, short drip, pack of 50, recommended receiver plate 203942-100	0.7 µm	0.3	201001-100
384-well polypropylene with glass fiber membrane, long drip, pack of 10, recommended receiver plate 204586-100	0.7 µm	0.14	201027-100
96-well polypropylene with glass fiber membrane, long drip, pack of 25, recommended receiver plate 201276-100	1 µm	0.8	200935-100
96-well polypropylene with glass fiber membrane, , short drip, pack of 25, recommended receiver plate 201276-100	1 µm	0.8	200963-100
96-well polypropylene with glass fiber membrane, long drip, pack of 25, recommended receiver plate 201276-100	1 µm	0.4	200949-100
96-well polypropylene with glass fiber membrane, short drip, pack of 25, recommended receiver plate 204600-100	1 µm	0.4	200975-100
96-well polypropylene with glass fiber membrane, long drip, pack of 50, recommended receiver plate 203942-100	1 µm	0.3	200985-100
96-well polypropylene with glass fiber membrane, short drip, pack of 50, recommended receiver plate 203942-100	1 µm	0.3	200999-100
96-well polypropylene with hydrophobic polyvinylidene fluoride membrane on a polypropylene support, short drip, pack of 50, recommended receiver plate 203942-100	1 µm	0.3	201023-100
24-well polypropylene with glass fiber membrane, long drip, pack of 25, recommended receiver plate 201272-100	1.2 µm	10	204474-100
96-well polypropylene with glass fiber membrane, long drip, pack of 25, recommended receiver plate 201276-100	1.2 µm	0.8	200929-100
96-well polypropylene with glass fiber membrane, short drip, pack of 25, recommended receiver plate 201276-100	1.2 µm	0.8	200957-100
96-well polypropylene with glass fiber membrane, long drip, pack of 25, recommended receiver plate 204600-100	1.2 µm	0.4	200941-100
96-well polypropylene with glass fiber membrane, short drip, pack of 25, recommended receiver plate 204600-100	1.2 µm	0.4	200969-100
96-well polypropylene with glass fiber membrane, short drip, pack of 50, recommended receiver plate 203942-100	1.2 µm	0.3	200993-100
384-well polypropylene with glass fiber membrane, long drip, pack of 10, recommended receiver plate 204586-100	1.2 µm	0.14	201025-100
96-well polypropylene with glass fiber + 25 µm polyethylene support membranes, long drip, pack of 25, recommended receiver plate 201240-100	3 µm	2	204495-100
96-well polypropylene with glass fiber membrane, long drip, pack of 50, recommended receiver plate 203942-100	3 µm	0.3	201021-100
96-well polypropylene with glass fiber membrane, long drip, pack of 25, recommended receiver plate 201240-100	5 µm	2	201702-100
384-well polypropylene with glass fiber membrane, long drip, pack of 10, recommended receiver plate 204586-100	5 µm	0.14	201033-100
96-well polypropylene with polyethylene frit, short drip, pack of 25, recommended receiver plate 201276-100	7 µm	0.8	202501-100
96-well polypropylene with glass fiber membrane, long drip, pack of 25, recommended receiver plate 204600-100	7 µm	0.4	202400-100

Description	Pore Size	Well Volume (ml)	Part Number
24-well polypropylene with polyethylene frit, long drip, pack of 15, recommended receiver plate 202061-100	25 µm	15	201415-100
24-well polypropylene with polyethylene membrane, long drip, pack of 25, recommended receiver plate 201272-100	25 µm	10	204471-100
48-well polypropylene with polyethylene frit, long drip, pack of 15, recommended receiver plate 201236-100	25 µm	7.5	201003-100
48-well polypropylene with polyethylene frit, EXTRA long drip, pack of 25, recommended receiver plate 201238-100	25 µm	5	200925-100
48-well polypropylene with polyethylene frit, long drip, pack of 25, recommended receiver plate 201238-100	25 µm	5	200923-100
96-well polypropylene with polyethylene frit, extra long drip, pack of 25, recommended receiver plate 201240-100	25 µm	2	200921-100
96-well polypropylene with polyethylene frit, long drip, pack of 25, recommended receiver plate 201240-100	25 µm	2	200919-100
96-well polypropylene with ultrahigh molecular weight polyethylene membrane, long drip, pack of 25, recommended receiver plate 201240-100	25 µm	2	201005-100
96-well polypropylene half skirt with polyethylene frit, long drip with turbine, pack of 25, recommended receiver plate 201276-100	25 µm	1	201718-100
96-well polypropylene with high molecular weight polyethylene membrane, long drip, pack of 25, recommended receiver plate 201236-100	25 µm	0.8	200927-100
96-well polypropylene with high molecular weight polyethylene membrane, short drip, pack of 25, recommended receiver plate 201276-100	25 µm	0.8	200955-100
96-well polypropylene with polyethylene frit, long drip, pack of 25, recommended receiver plate 201276-100	25 µm	0.8	200953-100
96-well polypropylene with high molecular weight polyethylene membrane, long drip, pack of 25, recommended receiver plate 204600-100	25 µm	0.4	200939-100
96-well polypropylene with high molecular weight polyethylene membrane, short drip, pack of 25, recommended receiver plate 204600-100	25 µm	0.4	200967-100
96-well polypropylene with high molecular weight polyethylene membrane, long drip, pack of 50, recommended receiver plate 204600-100	25 µm	0.3	200989-100
96-well polypropylene with high molecular weight polyethylene membrane, short drip, pack of 50, recommended receiver plate 204600-100	25 µm	0.3	200991-100
384-well polypropylene with polyethylene frit, long drip, pack of 10, recommended receiver plate 204586-100	25 µm	0.14	201035-100
384-well polypropylene with no filter, long drip, pack of 10, recommended receiver plate 204586-100	NA	0.14	201031-100
96-well polypropylene with no filter, extra long drip, pack of 25, recommended receiver plate 201240-100	NA	2	202951-100

*First time using "My Favorites?" You will be asked to enter your email address for account verification. If you have an existing Agilent account, you will be able to log in. However, if you don't have a registered Agilent account, you will need to [register](#) for one. This feature is valid only in regions that are e-commerce enabled. All items also can be ordered through your regular sales and distributor channels.

See the difference for yourself. Request your free sample:

explore.agilent.com/microplate-sample

Find microplates for any workflow:

www.agilent.com/en/promotions/microplate-for-any-workflow

DE44280.4359722222

This information is subject to change without notice.

© Agilent Technologies, Inc. 2021
 Printed in the USA, April 6, 2021
 5994-3223EN

