



Innovadyne Nanobuilder

Device Driver Quick Reference

Original Instructions

Notices

© Agilent Technologies, Inc. 2012

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.

User Guide Part Number

G5415-90038

Edition

Revision 01, December 2012

Contact Information

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

Technical Support: 1.800.979.4811
or +1.408.345.8011
service.automation@agilent.com

Customer Service: 1.866.428.9811
or +1.408.345.8356
orders.automation@agilent.com

European Service: +44 (0) 845 712 5292
euroservice.automation@agilent.com

Documentation feedback:
documentation.automation@agilent.com

Web:
www.agilent.com/lifesciences/automation

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 sub-contract, 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

 **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.**

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.

Innovadyne Nanobuilder Device Driver Quick Reference

This document contains the following topics:

- [About this document](#)
- [Innovadyne Nanodrop Nanobuilder Diagnostics](#)
- [Innovadyne Nanodrop protocol tasks](#)

About this document

What this guide covers

The Innovadyne Nanobuilder is a software that automates operations in Innovadyne liquid-handling devices, such as the Nanodrop. The Nanodrop liquid handler and the Nanobuilder software can be integrated into a lab automation system that is controlled by the VWorks software. The sequences created in Nanodrop Nanobuilder are used in VWorks protocols.

This document provides a quick reference of the commands, selections, and parameters in the following:

- [Innovadyne Nanodrop Nanobuilder Diagnostics](#)
- [Innovadyne Nanodrop protocol tasks](#)

Assumptions

This document assumes that you are familiar with the VWorks software and that you know how to:

- Add a device to a device file.
- Open the device diagnostics software.
- Create and manage profiles.
- Create a protocol and add protocol tasks.

Software version

The functions described in this document are available in the following software:

- VWorks Automation Control 11.4 or later
- Innovadyne Nanobuilder 4.8 or later

Related documents

Use this document in conjunction with the following:

- [VWorks Automation Control Setup Guide](#)
- [VWorks Automation Control User Guide](#)
- Lab automation system user guide, such as the [BioCel System User Guide](#)
- User documentation for the Innovadyne Nanodrop liquid handler and Nanobuilder software

For user information about other Automation Solutions products, you can search the product knowledge base or download the latest version of a PDF file from the Agilent Technologies website at:

www.agilent.com/chem/askb.

Innovadyne Nanodrop Nanobuilder Diagnostics

About Innovadyne Nanodrop Nanobuilder Diagnostics

You use Innovadyne Nanodrop Nanobuilder Diagnostics to:

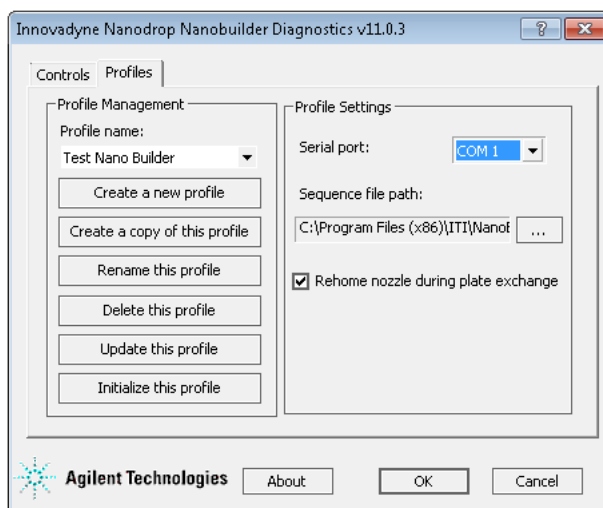
- *Create and manage profiles.* A profile allows you to set up communication between the Nanobuilder and the controlling computer.
- *Run or edit sequence files.* Using the available commands, you can run or edit the selected sequence file.

This topic provides a quick reference of the commands, selections, and parameters in the following Nanodrop Nanobuilder Diagnostics tabs:

- [Profiles tab](#)
- [Controls tab](#)

For instructions on adding the Nanodrop device to your device file and opening Nanodrop Nanobuilder Diagnostics, see the lab automation system user guide, such as the *BioCel System User Guide*.

Profiles tab



Profile Management area

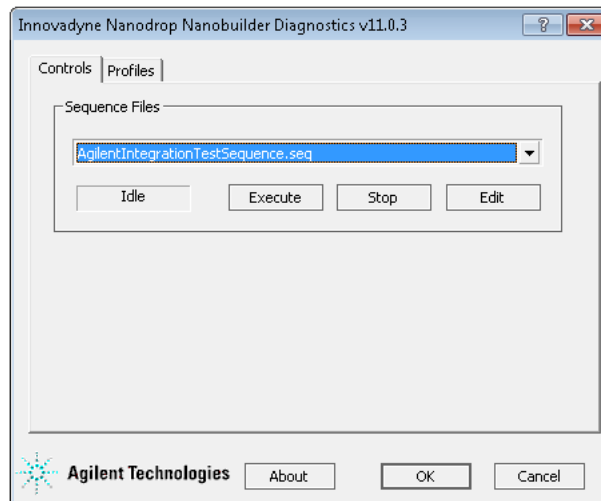
Command	Description
Profile name	Displays the selected profile. Also allows you to select from the list of available profiles.
Create a new profile	Creates a new profile.
Create a copy of this profile	Creates a duplicate copy of the selected profile.
Rename this profile	Renames the selected profile.
Delete this profile	Deletes the selected profile.

Command	Description
Update this profile	Saves changes to the selected profile.
Initialize this profile	Initiates communication with the device using the selected profile.

Profile Settings area

Setting or option	Description
Serial port	The first of the two controlling computer COM ports that is connected to the Nanodrop liquid handler. IMPORTANT The Nanodrop liquid handler requires two consecutive serial ports to operate. The serial port specification in the Profile Settings area should be the first of the two ports. For example, if the Nanodrop liquid handler is connected to COM ports 1 and 2, then you should select COM1 in the Profile Settings area.
Sequence file path	The folder in which the Nanobuilder saves all sequence files.
Rehome nozzle during plate exchange	The option to move the dispense head to the home position while the microplates are being exchanged.

Controls tab



Sequence Files area

Selection or command	Description
Sequence files list	The list of sequence files that are stored in the folder specified in the Profiles tab. Select the sequence file that you want to run or edit. For instructions on how to create and manage sequence files, see the Nanobuilder user documentation.
Status box	The status of the Nanodrop liquid handler. For example, if the device is not performing any task, the status Idle displays in the box.
Execute	Runs the selected sequence file.
Stop	Stops the sequence that is currently running on the Nanodrop liquid handler.
Edit	Opens the sequence file in the Nanobuilder so that you can edit it.

Innovadyne Nanodrop protocol tasks

About the tasks

You can add the following protocol tasks to run sequences on the Nanodrop liquid handler:

- [SubProcess](#)
- [Execute Sequence](#)


The Nanodrop Nanobuilder SubProcess and Execute Sequence tasks are associated with the Nanodrop liquid handlers only. This topic describes the two device-specific tasks.

Requirements

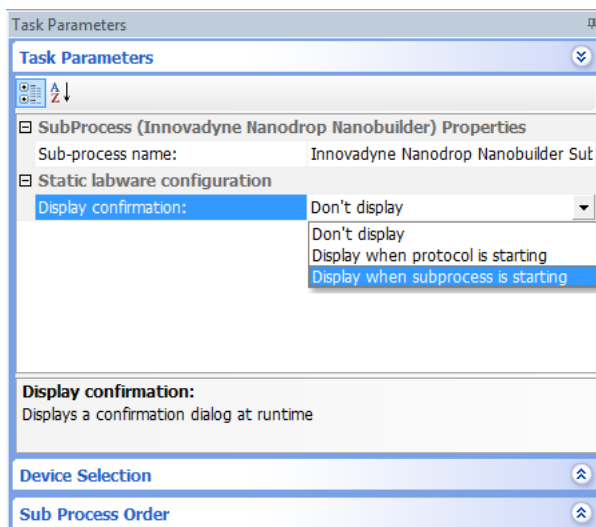
Make sure the desired sequence files are stored in the folder specified in the Nanobuilder Nanodrop Diagnostics Profiles tab.

SubProcess

Description

The SubProcess ( SubProcess (Innovadyne Nanodrop Nanobuilder)) task indicates the start of a protocol subroutine that employs the Nanodrop liquid handler. Within the subprocess, you can add the Execute Sequence task that is unique to the device. You can expand or collapse the subprocess to show or hide the subprocess tasks.

Task parameters



Parameter	Description
Sub-process name	The name of the Nanodrop subprocess. Select from the list of available subprocesses currently in the protocol.
Display confirmation	The option to display a message at the beginning of the protocol run or the subprocess to remind you to verify that the physical locations of the labware match what you specified in the software.

Device selection

You must select the device for the subprocess task.

To select the device for the subprocess task:

- 1 Click **Device Selection**.
- 2 In the **Device Selection** pane, double-click the desired device in the **Devices available to perform** task area to move it to the **Devices involved in task** area.

SubProcess Order


If more than one subprocess uses the same configured labware, and the subprocesses are in different protocol processes, you can specify the sequence in which the subprocesses will be performed.

To specify the sequence in which the subprocesses will be performed on the same configured labware:

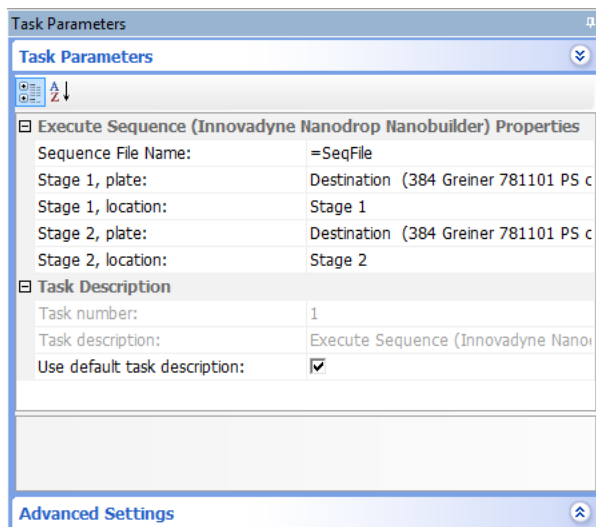
- 1 In the protocol process, select the subprocess that contains the task that uses the configured labware.
- 2 In the **Task Parameters** area, click **Sub Process Order**.
- 3 In the **Sub Process Order** pane, double-click the subprocess names to rearrange the order.

Execute Sequence

Description

The Execute Sequence ( Execute Sequence (Innovadyne Nanodrop Nanobuilder)) task runs the selected sequence on the Nanodrop liquid handler.

Task parameters



Parameter	Description
Sequence file name	The sequence file that the task will run. Select the desired sequence file from the list.
Stage 1, plate	The type of labware that will be placed at the Nanodrop stage.
Stage 1, location	The name of the Nanodrop stage. Select from the list of location names.
Stage 2, plate	The type of labware that will be placed at the second Nanodrop stage, if applicable.
Stage 2, location	The name of the second Nanodrop stage, if applicable. Select from the list of location names.



Quick Reference

G5415-90038

Revision 01, December 2012