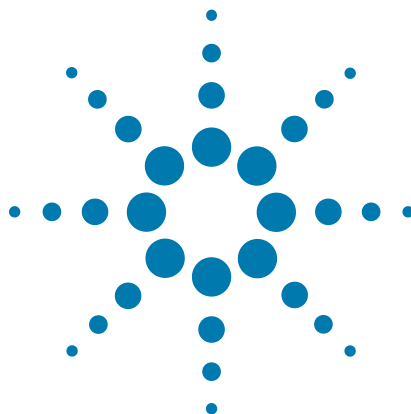




Roche Applied Science Cellavista Analyzer

Device Driver Quick Reference

Original Instructions



Agilent Technologies

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

G5415-90007

Edition

Revision 00, 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.

Roche Applied Science Cellavista Analyzer Device Driver Quick Reference

This document contains the following topics:

- [About this document](#)
- [Roche Applied Science Cellavista Diagnostics](#)
- [Cellavista Analyzer protocol tasks](#)
- [Cellavista XML RPC Server](#)

About this document

What this guide covers

The Roche Applied Science Cellavista Analyzer is a microplate reader for cell-based applications. The device can be integrated into a lab automation system that is controlled by the VWorks software.

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

- [Roche Applied Science Cellavista Diagnostics](#)
- [Cellavista Analyzer protocol tasks](#)
- [Cellavista XML RPC Server](#)

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

This document describes the functions that are available in the following software:

- VWorks Automation Control 11.4 or later
- Cellavista software 2.0 or later
- Cellavista XML RPC Server 1.0 or later

The VWorks software runs on a Windows 7 computer. Because the Cellavista software cannot run on a Windows 7 computer, it is installed on a different computer that is connected to the Cellavista Analyzer. This dual-computer setup requires a program, the Cellavista XML RPC Server, to manage communication between the Cellavista software and the VWorks software. The Cellavista XML RPC Server is installed on the Cellavista computer.

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 Roche Applied Science Cellavista Analyzer

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.

Roche Applied Science Cellavista Diagnostics

About Roche Applied Science Cellavista Diagnostics

You use Roche Applied Science Cellavista Diagnostics to:

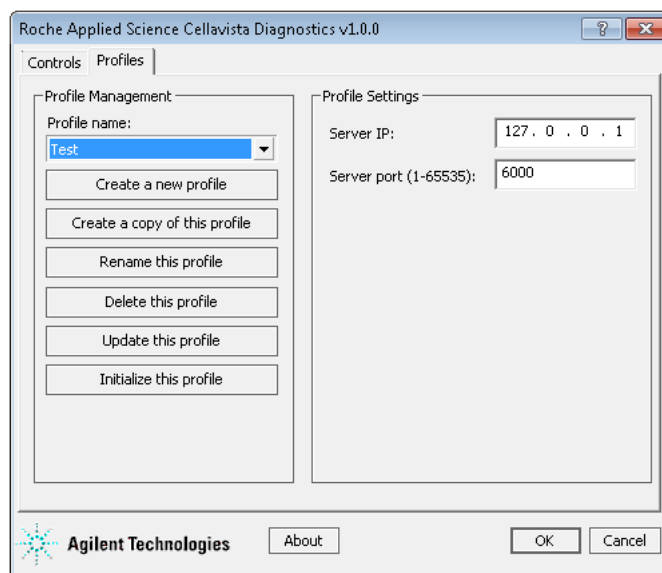
- *Create and manage profiles.* A profile allows you to set up communication between the Cellavista Analyzer and the controlling computer.
- *Read a microplate.* Using the available commands, you can select a measurement file, start reading a microplate, and stop the reading process.

This topic provides a quick reference of the commands, selections, and parameters in the following Roche Applied Science Cellavista Diagnostics tabs:

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

For instructions on adding the Cellavista Analyzer device to your device file and opening Roche Applied Science Cellavista 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.

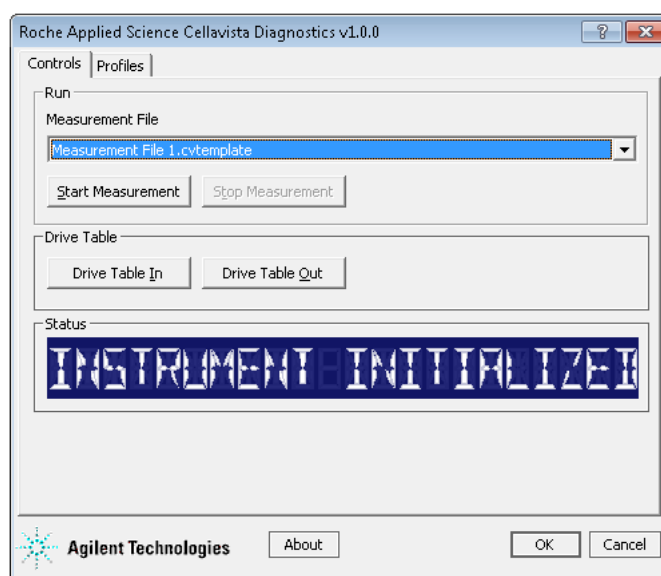
Command	Description
Create a new profile	Creates a new profile.
Create a copy of this profile	Creates a duplicate copy of the selected profile.
Rename this profile	Renames the selected profile.
Delete this profile	Deletes the selected profile.
Update this profile	Saves changes to the selected profile.
Initialize this profile	Initiates communication with the device using the selected profile.

Profile Settings area

Setting or option	Description
Server IP	The IP address of the computer that is running the Cellavista software and the Cellavista XML RPC Server.
Server port	The IP address of the computer that is connected to the computer running the Cellavista software and the Cellavista XML RPC Server.

Controls tab

IMPORTANT Before using the commands in diagnostics, make sure the Cellavista computer is turned on, and the Cellavista software and the Cellavista XML RPC Server are both running. Make sure the Cellavista XML RPC Server has been configured correctly. See [Cellavista XML RPC Server](#).



Run area

Selection or command	Description
Measurement File	The measurement file that you want to use when reading a microplate. Select from the list of available measurement files. Notes: - The measurement files are created in the Cellavista software. For instructions on how to create the files, see the Cellavista software user documentation. - The location of the files is specified in the Cellavista XML RPC Server. See Cellavista XML RPC Server.
Start Measurement	Starts the microplate reading process on the Cellavista Analyzer.
Stop Measurement	Stops the current reading process on the Cellavista Analyzer.

Drive Table area

Command	Description
Drive Table In	Moves the drive table into the Cellavista Analyzer. The drive table hosts the microplate. Moving it into the analyzer prepares the microplate for the reading process.
Drive Table Out	Moves the drive table out so that the automation system robot can access the microplate.

Status area

Displays the Cellavista Analyzer's current state.

Cellavista Analyzer protocol tasks

About the tasks

You can add the following protocol tasks to read microplates in the Cellavista Analyzer:

- [Start Measurement](#)
- [Stop Measurement](#)

The Start Measurement and Stop Measurement tasks are associated with the Cellavista Analyzer only. This topic describes the two device-specific tasks.

IMPORTANT Before running the protocol that contain these tasks, make sure the Cellavista computer is turned on, and the Cellavista software and the Cellavista XML RPC Server are both running. Make sure the Cellavista XML RPC Server has been configured correctly. See [Cellavista XML RPC Server](#).

Start Measurement

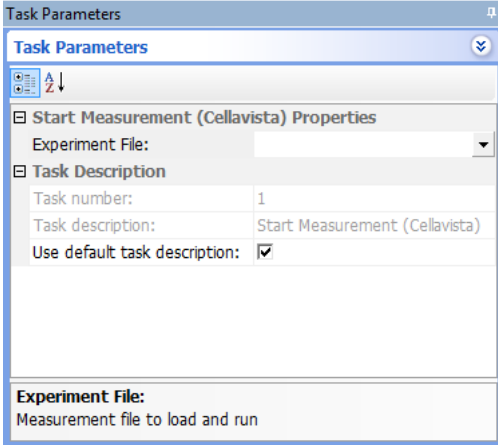
Description

The Start Measurement ( Start Measurement (Cellavista)) task starts the microplate reading process in the Cellavista Analyzer.

When the protocol run reaches the Start Measurement task, the Cellavista Analyzer drive table will automatically move out to receive the microplate, and then move in to prepare the microplate for the reading process.

Note: The Start Measurement task is only available in the Main Protocol.

Task parameters



The screenshot shows the 'Task Parameters' dialog box for the 'Start Measurement (Cellavista)' task. The dialog has a title bar 'Task Parameters' and a close button. Below the title bar is a section 'Task Parameters' with a dropdown arrow. Underneath is a section 'Start Measurement (Cellavista) Properties' with a dropdown arrow. This section contains three fields: 'Experiment File:' with a dropdown menu, 'Task number:' with a text box containing '1', and 'Task description:' with a text box containing 'Start Measurement (Cellavista)'. Below these is a checkbox 'Use default task description:' which is checked. At the bottom of the dialog is a section 'Experiment File:' with a text box containing 'Measurement file to load and run'.

Parameter	Description
Experiment File	The measurement file to use when reading a microplate. Select from the list of available measurement files. The measurement files are created in the Cellavista software. For instructions on how to create the files, see the Cellavista software user documentation.
Task number	The number that indicates the position of the task in the protocol.
Task description	The description of the task.
Use default task description	The option to use the default task description or provide your own description for the task. Select the check box to use the default description. Clear the check box to provide your own description.

Stop Measurement

Description

The Stop Measurement ( Stop Measurement (Cellavista)) task stops the microplate reading process in the Cellavista Analyzer.

When the protocol run reaches the Stop Measurement task, the Cellavista Analyzer stops the reading process, and the drive table will automatically move out so that the automation system robot can access the microplate.

Note: The Stop Measurement task is only available in the Main Protocol.

Task parameters

The Stop Measurement task does not have any parameters.

Cellavista XML RPC Server

About the Cellavista XML RPC Server

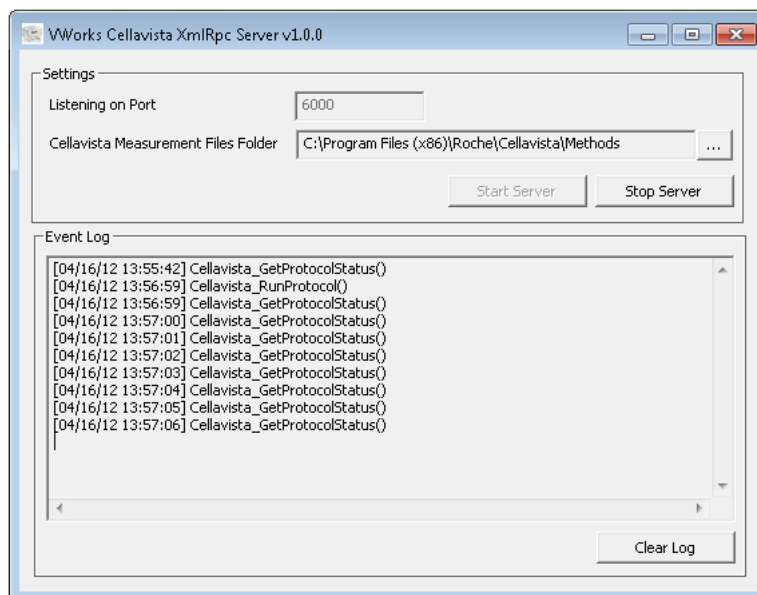
The Cellavista XML RPC Server is a software that is installed on the Cellavista computer and manages the communication between the Cellavista software and the VWorks software. To ensure successful communication, the following information is required when configuring the server:

- The IP port number that the VWorks software uses to communicate with the Cellavista software.

Note: The IP port number is also specified in the Roche Applied Science Cellavista Diagnostics Profiles tab.

- The folder that contains the measurement files.

To start the Cellavista XML RPC Server, double-click the Cellavista XML RPC Server icon on the Cellavista computer. The Cellavista XML RPC Server window opens.



Settings area

Selection or command	Description
Listening on Port	The IP port number that the VWorks software uses to communicate with the Cellavista software. <i>Note:</i> The IP port number is also specified in the Roche Applied Science Cellavista Diagnostics Profiles tab. Make sure they match.
Cellavista Measurement Files Folder	The location of the measurement files on the Cellavista computer. Click the browse button, then locate and select the folder. During startup, the server will retrieve the list of Cellavista measurement files and send it to the VWorks software.
Start Server	Starts serving the requests from the VWorks software and the Cellavista software.
Stop Server	Stops serving the requests from the VWorks software and the Cellavista software.

Event Log area

Display area or command	Description
Log display area	Displays the actions that occur in the Cellavista software.
Clear Log	Removes the existing events listed in the log display area.



Quick Reference

G5415-90007

Revision 00, December 2012