



Labcyte Latitude

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

G5415-90005

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.

Labcyte Latitude Device Driver Quick Reference

This document contains the following topics:

- [About this document](#)
- [Labcyte Latitude LX Diagnostics](#)
- [Labcyte Latitude protocol tasks](#)



About this document

What this guide covers

The Labcyte Latitude is a low-volume bulk reagent dispenser device that 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:

- [Labcyte Latitude LX Diagnostics](#)
- [Labcyte Latitude 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 VWorks Automation Control 11.4 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 Labcyte Latitude

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.

Labcyte Latitude LX Diagnostics

About Labcyte Latitude LX Diagnostics

You use Labcyte Latitude LX Diagnostics to:

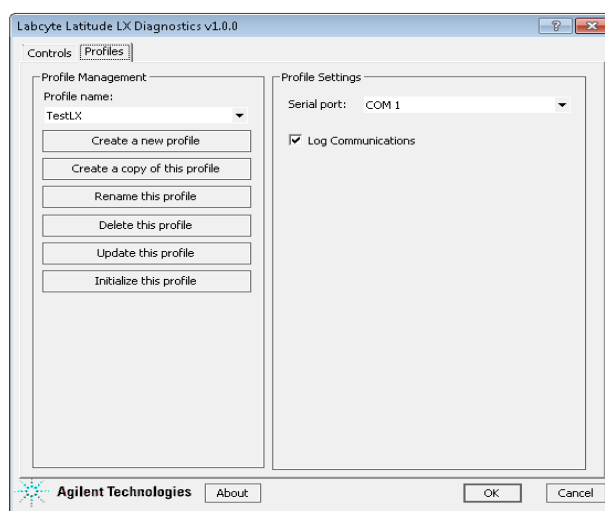
- *Create and manage profiles.* A profile allows you to set up communication between the Latitude and the controlling computer.
- *Home or run methods.* Using the available commands, you can home the device or run liquid transfer methods.

This topic provides a quick reference of the commands, selections, and parameters in the following Labcyte Latitude LX Diagnostics tabs:

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

For instructions on adding the Latitude device to your device file and opening Labcyte Latitude LX 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.

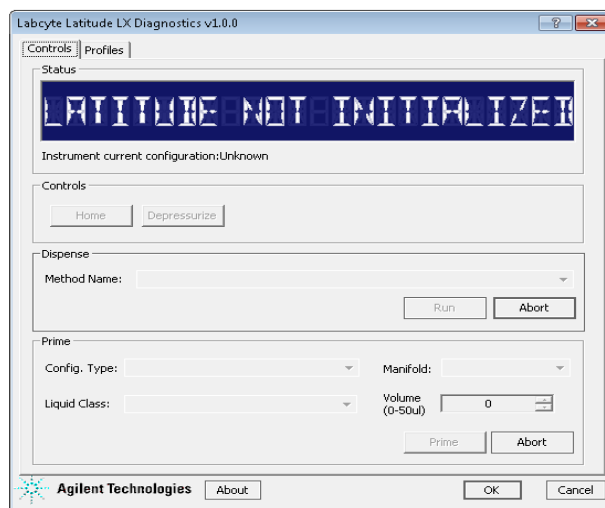
Command	Description
Update this profile	Saves changes to the selected profile.
Initialize this profile	Initiates communication with the device using the selected profile.

Profile Settings area

Setting or option	Description
Serial port	The controlling computer COM port that is connected to the Latitude.
Log communications	The option to record all communication between the Latitude and the controlling computer.

Controls tab

IMPORTANT Before using the commands to operate the Latitude, make sure the Latitude is in the Remote mode. To set the device in the Remote mode, initialize the device from the VWorks software or use the touchscreen on the Latitude.



Status area

Displays the Latitude's current state.

Controls area

Command	Description
Home	Moves the stage to the factory-defined home position.
Depressurize	Releases the pressure in the device. The operation might take up to 30-60 seconds, depending on the amount of pressure currently in the device.

Dispense area

Selection or command	Description
Method name	The method that instructs the Latitude on how to dispense the liquid. Select the desired method from the list.
Run	Pressurizes the Latitude and runs the selected method.
Abort	Aborts the currently running method.

Prime area

Selection or command	Description
Config type	The hardware configuration used to prime the Latitude. Select one of the following: 8x1 (8 tips, 1 liquid) 16x2 (16 tips, 2 liquids) 16x4 (16 tips, 4 liquids)
Manifold	The manifold to use. The manifold selection depends on the configuration type you selected.
Liquid class	The liquid class to use for priming the Latitude. Select from the list.
Volume (0–50 µL)	The volume of liquid to use during priming.
Prime	Pressurizes and primes the Latitude using the specified configuration, manifold, liquid class, and volume.
Abort	Aborts the current priming action.

Labcyte Latitude protocol tasks

About the tasks

You can add the following protocol tasks to run the Labcyte Latitude device:

- [Prime Latitude](#)
- [Run Latitude Method](#)

The Prime Latitude and Run Latitude tasks are associated with the Labcyte Latitude only. This topic describes the two device-specific tasks.

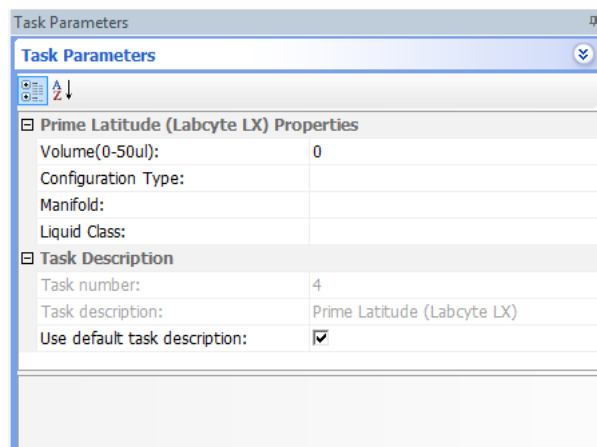
IMPORTANT Before running the protocol that contain these tasks, make sure the Latitude is in the Remote mode. To set the Latitude in the Remote mode, initialize the device from the VWorks software or use the touchscreen on the Latitude.

Prime Latitude

Description

The Prime Latitude ( Prime Latitude (Labcyte LX)) task pressurizes and primes the Latitude device using the specified information.

Task parameters



Task Parameters

Task Parameters

Prime Latitude (Labcyte LX) Properties

Volume(0-50ul): 0

Configuration Type:

Manifold:

Liquid Class:

Task Description

Task number: 4

Task description: Prime Latitude (Labcyte LX)

Use default task description: ☒

Display or command	Description
Volume (0–50 µL)	The volume of liquid used to use during priming.
Configuration type	The hardware configuration used to prime the Latitude. Select one of the following: • 8x1 (8 tips, 1 liquid) • 16x2 (16 tips, 2 liquids) • 16x4 (16 tips, 4 liquids)

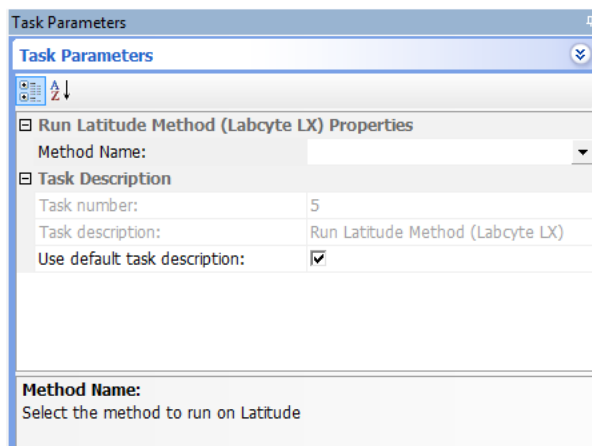
Display or command	Description
Manifold	The manifold to use. The manifold selection depends on the configuration type you selected.
Liquid class	The liquid class to use for priming the Latitude. Select from the list.

Run Latitude Method

Description

The Run Latitude Method ( Run Latitude Method (Labcyte LX)) task pressurizes the Latitude and runs the selected method.

Task parameters



Task Parameters

Task Parameters

Run Latitude Method (Labcyte LX) Properties

Method Name: ▼

Task Description

Task number: 5

Task description: Run Latitude Method (Labcyte LX)

Use default task description: ☒

Method Name:
Select the method to run on Latitude

Display or command	Description
Method name	The method that instructs the Latitude on how to dispense the liquid. Select the desired method from the list.
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.



Agilent Technologies

Quick Reference

G5415-90005

Revision 00, December 2012