

# **Centrifuge Diagnostics version 8**

## **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

G5405-90005

## Edition

Revision 00, August 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](mailto:service.automation@agilent.com)

Customer Service: 1.866.428.9811  
or +1.408.345.8356  
[orders.automation@agilent.com](mailto:orders.automation@agilent.com)

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

Documentation feedback:  
[documentation.automation@agilent.com](mailto:documentation.automation@agilent.com)

Web:  
[www.agilent.com/lifesciences/automation](http://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.**

## Centrifuge Diagnostics Version 8 Quick Reference

The Centrifuge Diagnostics allows you to create a profile and establish communication with the Centrifuge, set bucket teachpoints, and control the motions of the Centrifuge.

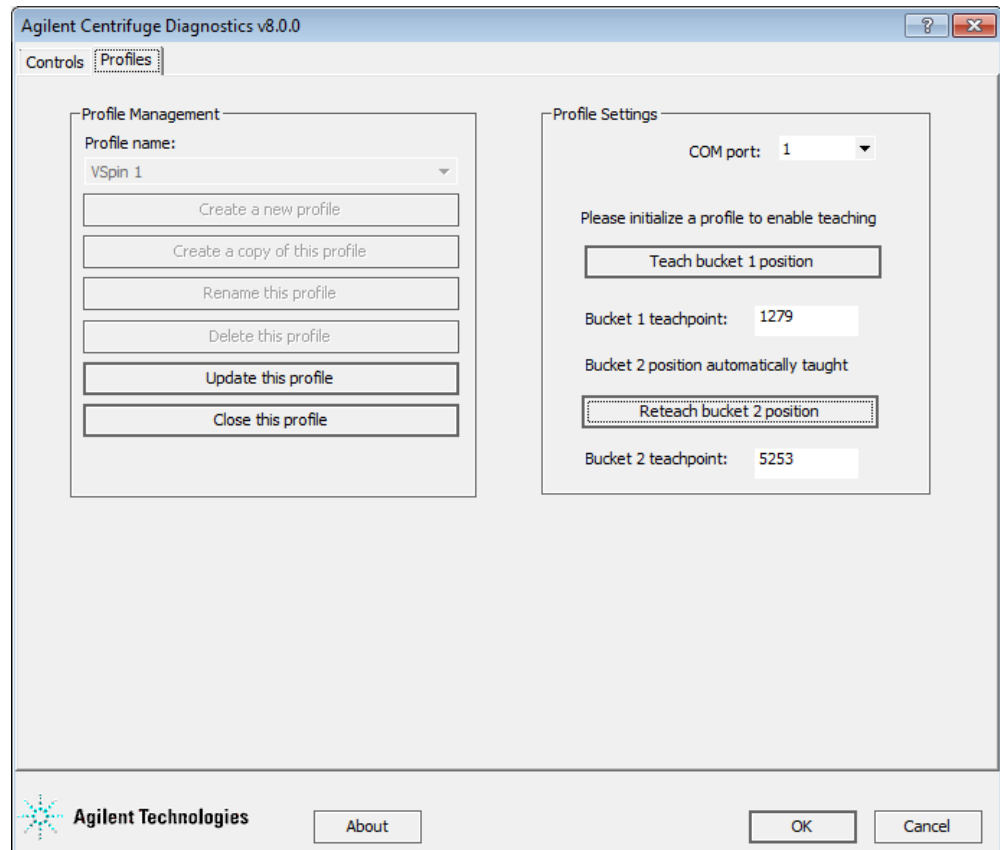
This document provides a quick reference of the commands, selections, and parameters in the Centrifuge Diagnostics dialog box. The topics are:

- “Profiles tab” on page 2
- “Controls tab” on page 4
- “Teach Bucket 1 and Teach Bucket 2 dialog boxes” on page 7

For user information about related 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](http://www.agilent.com/chem/askb).

## Profiles tab



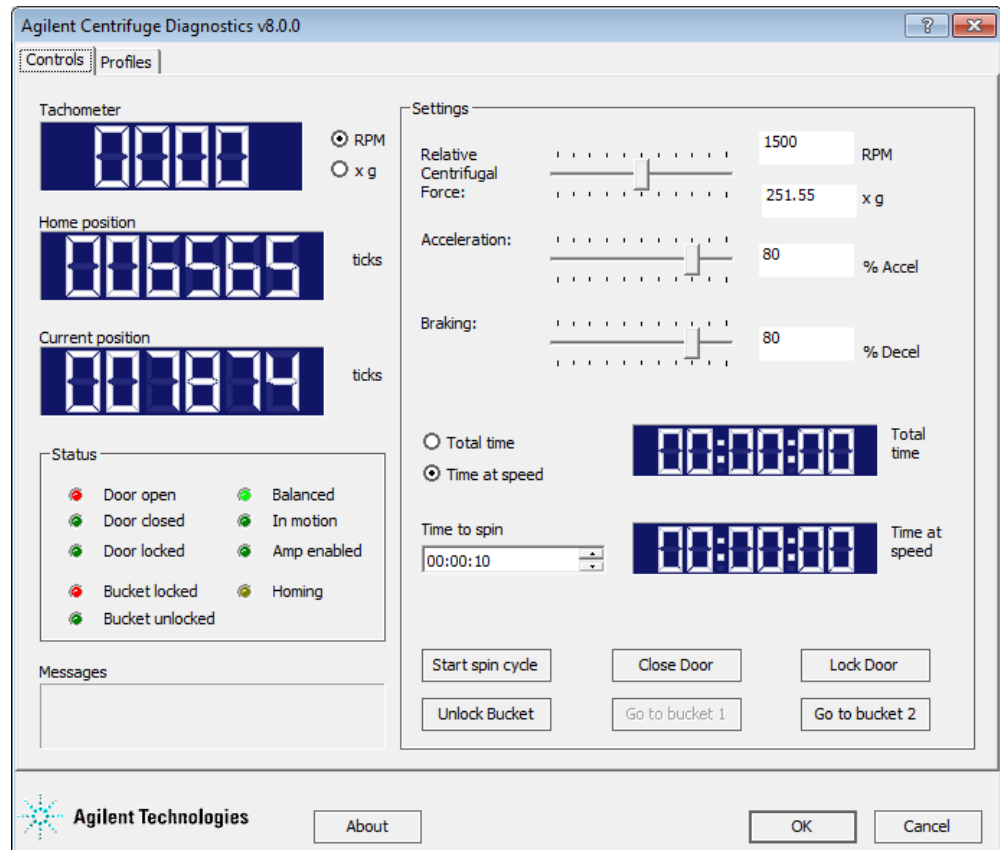
### Profile Management area

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.
Update this profile	Saves changes to the selected profile.
Initialize this profile	Initiates communication with the Centrifuge using the selected profile.

## Profile Settings area

Selection or command	Description
COM port	The controlling computer COM port that is connected to the Centrifuge.
Teach bucket 1 position	<p>The command that opens the Teach Bucket 1 dialog box to allow you to set the bucket-1 teachpoint. See “<a href="#">Teach Bucket 1 and Teach Bucket 2 dialog boxes</a>” on page 7.</p> <p><b>IMPORTANT</b> Make sure you initialize the desired profile before setting the teachpoint.</p> <p><b>IMPORTANT</b> Make sure you unlock the bucket before you set the teachpoint. When the bucket is unlocked, you can move the bucket to the desired position before setting the teachpoint.</p> <p><i>Note:</i> When you set the bucket-1 teachpoint, the software automatically sets the bucket-2 teachpoint. If a bucket-2 teachpoint already exists, you have the option of keeping the existing teachpoint or use the new auto-set bucket-2 teachpoint.</p>
Bucket 1 teachpoint	The bucket-1 teachpoint position. When you set the bucket 1 teachpoint, a value displays to indicate the teachpoint position.
Reteach bucket 2 position	<p>The command that opens the Teach Bucket 2 dialog box to allow you to set the bucket-2 teachpoint. See “<a href="#">Teach Bucket 1 and Teach Bucket 2 dialog boxes</a>” on page 7.</p> <p><b>IMPORTANT</b> Make sure you initialize the desired profile before setting the teachpoint.</p> <p><b>IMPORTANT</b> Make sure you unlock the bucket before you set the teachpoint. When the bucket is unlocked, you can move the bucket to the desired position before setting the teachpoint.</p> <p>You typically do not have to set the bucket-2 teachpoint, because it is set automatically when you set the bucket-1 teachpoint. Use this command only if you want to fine-tune the bucket-2 teachpoint.</p>
Bucket 2 teachpoint	The bucket-2 teachpoint position. When you set the bucket-2 teachpoint (or when it is set automatically), a value displays to indicate the teachpoint position.

## Controls tab



## Display area

Display title	Description
Tachometer	The current spin speed. You can select the desired spin speed unit to display: <b>RPM</b> (revolutions per minute), or <b>x g</b> (a multiple of gravity).
Home position	The home position value. The home position is also the bucket-1 teachpoint.
Current position	The current position of bucket 1.

## Status area

Display title	Description
Door open	Indicates that the Centrifuge door is open.
Door closed	Indicates that the Centrifuge door is closed.
Door locked	Indicates that the door is locked.
Bucket locked	Indicates that the bucket at the door is locked in position to prevent the rotor from moving.
Bucket unlocked	Indicates that the front bucket is unlocked and the rotor can move.
Balanced	Indicates that the buckets are out of balance during spinning, because one of the buckets has more mass than the other bucket.
In motion	Indicates that the buckets are spinning.
Homing	Indicates that the Centrifuge is in the homing process.

## Messages area

Displays error or status messages.

## Settings area

Setting or command	Description
Relative Centrifugal Force	The rotor speed, in RPM (revolutions per minute) and as a multiple of gravity.
Acceleration	The rate of centrifugation, as a percent of maximum acceleration.
Braking	The deceleration of the centrifuge, as a percent of maximum deceleration.
Total time	The specified spin time, including acceleration and braking.
Time at speed	The specified spin time, not including acceleration and braking.
Time to spin	The length of time to spin the microplates in the desired mode.
Start Spin Cycle	The command to start the spin.
Open Door	The command to open the Centrifuge door.
Lock Door	The command to lock the Centrifuge door.

Setting or command	Description
Lock/Unlock Bucket	<p>The command to lock or unlock the front bucket. Locking the front bucket prevents the rotor from moving. Unlocking the front bucket allows the rotor to move.</p> <p>Unlock the bucket when you want to set a teachpoint. Lock the bucket after you finish setting the teachpoints.</p>
Go to Bucket 1	The command to move bucket 1 to the door.
Go to Bucket 2	The command to move bucket 2 to the door.



## Teach Bucket 1 and Teach Bucket 2 dialog boxes

Display or command	Description
Current position	The current position value of the bucket.
Home position	The home position value.
Capture 1	The command that records the current position of the bucket.
Capture 2	<p><b>IMPORTANT</b> Move bucket 1 to the door before capturing its position. Move bucket 2 to the door before capturing its position.</p> <p>The command that records the current position of the bucket a second time.</p> <p>If you do not move the buckets, the second capture value should be the same as or close to the first capture value.</p> <p><i>Note:</i> The final bucket teachpoint is the average of the two captured position values. The software will not allow you to exit this dialog box until you click <b>Capture 2</b>.</p>
Capture average	The average of the two captured position values.
Teachpoint	The position value of the bucket teachpoint, based on the calculated Capture average.







**Quick Reference**

**G5405-90005**

Revision 00, August 2012