



400-DS Apparatus 7

Operator's Manual

Original Instructions

Notices

Manual Part Number

70-9064Rev D
March 2018

Copyright

© Agilent Technologies, Inc. 2018

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.

Agilent Technologies, Inc.
3501 Stevens Creek Blvd.
Santa Clara, CA 95052 USA

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

U.S. Government Restricted Rights. Software and technical data rights granted to the federal government include only those rights customarily provided to end user customers. Agilent provides this customary commercial license in Software and technical data pursuant to FAR 12.211 (Technical Data) and 12.212 (Computer Software) and, for the Department of Defense, DFARS 252.227-7015 (Technical Data - Commercial Items) and DFARS 227.7202-3 (Rights in Commercial Computer Software or Computer Software Documentation).

Safety Notices

CAUTION

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.

WARNING

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.

Content

1 Safety 7

General Safety Information 8

Safety Standards 8

General 8

Warning 9

Caution 9

Note 9

A 'Note' appears in the manual to give advice or information. 9

Electrical Hazards 10

Before Applying Power 10

Ground the Instrument 11

Do Not Operate in an Explosive Atmosphere 11

Do Not Remove Instrument Cover 11

Do Not Modify the Instrument 12

In Case of Damage 12

Disconnecting Device 12

Fuses and Fuse Replacement 13

Leak and Waste Handling 14

Information Symbols 15

2 Introduction 17

Conventions Used in this Manual 19

Serial Number Format 19

3 Setting Up the 400-DS (G7975A) 21

Initial Setup 22

Unpacking Procedure 22

Environmental Requirements for Installation 23

Clearance 23

Parts and Accessories 24

Leveling Your Apparatus 25

Electrical Connection 25

Content

Cable Connections	27
Heater Jacket / Sample Holder Installation	28
Tubing Connections	29
PC Setup	30
Initial Power Up	30
Software Installation and Setup	31
400-DS Workstation Software Installation	31
Local Security Policy	32
Starting 400-DS Workstation	33
400-DS Workstation Logon	34
Adding Users to the Application	34
4 Operating the 400-DS	37
400-DS Workstation Logon	38
System Configuration	41
Initial Setup of 400-DS Configuration	41
Copying a System Configuration	45
Deleting a System Configuration	45
System Configuration Report	46
Editing an Existing System Configuration	46
Recovering a Deleted Configuration	47
Show Audit Trail	47
Verify Integrity	48
Import/Export XML File	48
400-DS System Calibration	50
Temperature Calibration	50
Volume Calibration	54
Manual Control/Diagnostics	56
Verifying Sample Tray Control	56
Verifying Media Control	57
Verifying the Cell Temperature	58

Verifying Reciprocation	58
Evaporation Cover Control	59

Method Editor 60

Creating a Method	60
Copying Methods	67
Deleting Methods	67
Editing an Existing Method	67
Method Report	68
Recovering a Method	69
Method Audit Trail	69
Import/Export XML File	70

Active Channel Selection Valves 72

Running a Method 73

Test Reports 77

Electronic Signatures	78
-----------------------	----

Clean System 79

5 Troubleshooting and Maintenance 81

Preventive Maintenance 82

Calibration	82
Routine Maintenance (to be Performed Between Each Dissolution Method)	82
Sample Holders	83

Troubleshooting 84

Obtaining Warranty and Other Services	84
---------------------------------------	----

This page was intentionally left blank, except for this message.

1

Safety

General Safety Information 8

Electrical Hazards 10

Leak and Waste Handling 14

The 400-DS Apparatus 7 (G7975A) has been designed and tested so that when used properly you have an accurate, fast, flexible, and safe instrument.

If the equipment is used in a manner not specified by the manufacturer, the protection provided by the equipment may be impaired.

Operation of an Agilent 400-DS (G7975A) involves the use of aqueous liquids (or methanol or ethanol) and various pharmaceutical dosage forms. Unskilled, improper, or careless use of this instrument can create shock hazards, fire hazards, or other hazards which can cause death, serious injury to personnel, or severe damage to equipment and property.

Information on safety practices is provided with your instrument and operation manuals. Before using your instrument or accessories, you must thoroughly read these safety practices.

Observe all relevant safety practices at all times.

General Safety Information

The following general safety precautions must be observed during all phases of operation, service, and repair of this instrument. Failure to comply with these precautions or with specific warnings elsewhere in this manual violates safety standards of design, manufacture, and intended use of the instrument. Agilent Technologies assumes no liability for the customer's failure to comply with these requirements.

Safety Standards

This is a Safety Class I instrument (provided with terminal for protective earthing) and has been manufactured and tested according to international safety standards.

General

Do not use this product in any manner not specified by the manufacturer.

The protective features of this product may be impaired if it is used in a manner not specified in the operation instructions.

Warning

WARNING

A 'Warning' message appears in the manual when failure to observe instructions or precautions could result in death or injury.

Read all warnings and cautions carefully and observe them at all times.

Caution

CAUTION

A 'Caution' message appears in the manual when failure to observe instructions could result in damage to equipment (Agilent supplied and / or other associated equipment).

Note

NOTE

A 'Note' appears in the manual to give advice or information.

Electrical Hazards

The 400-DS (G7975A) contains electrical circuits, devices, and components operating at dangerous voltages. Contact with these circuits, devices, and components can cause death, serious injury, or painful electric shock.

Panels or covers that are retained by fasteners which require the use of a tool for removal may be opened only by Agilent-trained, Agilent-qualified, or Agilent-authorized service engineers. Consult the manuals or product labels supplied with the 400-DS to determine which parts are operator-accessible.

Application of the wrong supply voltage, connection of the instrument to an incorrectly wired supply outlet, or lack of proper electrical grounding can create a fire hazard or a potentially serious shock hazard and could seriously damage the instrument and any attached ancillary equipment.

Always use a three-wire outlet with ground connection which is adequately rated for the load. The installation must comply with local, state, and federal safety regulations.

Do not connect the instrument to the main power supply until you have made sure that the operating voltage is correctly set for the main power supply in the specific outlet in your laboratory to which the equipment will be connected.

Before Applying Power

WARNING

Wrong voltage range, frequency or cabling - personal injury or damage to the instrument.

Verify that the voltage range and frequency of your power distribution matches to the power specification of the individual instrument. Never use cables other than the ones supplied by Agilent Technologies to ensure proper functionality and compliance with safety or EMC regulations. Make all connections to the unit before applying power.

Ground the Instrument

WARNING

Missing electrical ground - electrical shock.

If your product is provided with a grounding type power plug, the instrument chassis and cover must be connected to an electrical ground to minimize shock hazard. The ground pin must be firmly connected to an electrical ground (safety ground) terminal at the power outlet. Any interruption of the protective (grounding) conductor or disconnection of the protective earth terminal will cause a potential shock hazard that could result in personal injury.

Do Not Operate in an Explosive Atmosphere

WARNING

Presence of flammable gases or fumes - explosion hazard.

Do not operate the instrument in the presence of flammable gases or fumes.

Do Not Remove Instrument Cover

WARNING

Instrument covers removed - electrical shock.

Do not remove the instrument covers or panels. Only Agilent authorized personnel are allowed to remove instrument covers or panels. The power cables and any external circuits should be disconnected before removing the instrument cover.

Do Not Modify the Instrument

Do not install substitute parts or perform any unauthorized modification to the product. Contact an Agilent Repair Center for service and repair to ensure that safety features are maintained.

In Case of Damage

WARNING

Damage to the module - personal injury (e.g., electrical shock, intoxication).

Instruments that appear damaged or defective should be made inoperative and secured against unintended operation until they can be repaired by qualified service personnel.

Disconnecting Device

WARNING

Power removal.

The module is partially energized when switched off, as long as the power cord is connected. In case of emergency, it must be possible to disconnect the instrument from the power line at any time. Make sure the power connector of the instrument can be easily reached and unplugged. Provide sufficient space behind the power socket of the instrument to unplug the cable.

Fuses and Fuse Replacement

WARNING

Double-pole neutral protection.

The instrument is provided having 2 mains fuses, so that one is in the active connector and one in the neutral connector.

WARNING

Externally accessible fuses, located in the fuse holder of the power inlet.

For exchanging the fuses, disconnect the power cord. Exchange the fuses only with fuses specified in this manual.

Leak and Waste Handling

WARNING

The handling of toxic, flammable and hazardous solvents, samples and reagents can hold health and safety risks.

- When working with these substances observe appropriate safety procedures (for example by wearing goggles, safety gloves and protective clothing) as described in the material handling and safety data sheet supplied by the vendor, and follow good laboratory practice.
- The volume of substances should be reduced to the minimum required for the analysis.
- Do not operate the instrument in an explosive atmosphere.
- Operate the instrument in a fume cupboard or in a well ventilated room.
- The residual free volume in the appropriate waste container must be large enough to collect the waste liquid.
- Check the filling level of the waste container regularly.
- To achieve maximal safety, check the correct installation regularly.
- The 400-DS (G7975A) may be used with up to 100% concentrations of Methanol and Ethanol. Other organic solvents have not been tested/validated for use.

Information Symbols

I

Switches main power on

0

Switches main power off



Indicates single-phase alternating current



Indicates the product complies with the requirements of one or more European Union (EU) directives.



Indicates an electrical hazard.



Indicates a warning.



Indicates specific equipment meets standards of safety. These products are safe for use in the workplace for North America.



Indicates that this product must not be disposed of as unsorted municipal waste.

All Agilent products that are subject to the WEEE directive shipped after August 13, 2005 are compliant with the WEEE marking requirements. Such products are marked with the “crossed out wheeled bin” WEEE symbol in accordance with European Standard EN 50419.

For more information on collection, reuse, and recycling systems, please contact your local/regional waste administration, your local distributor, or Agilent.



Indicates the product complies with regulatory compliance requirements of New Zealand and Australia.



2

Introduction

Conventions Used in this Manual 19

Serial Number Format 19

The Agilent 400-DS Dissolution Apparatus (G7975A) is the first compendial small-volume dissolution apparatus for testing medical devices such as drug-eluting stents (DES). This integrated, small-footprint dissolution apparatus features several first-in-class innovations that greatly increase the convenience of release rate testing of combinatorial drug products. The 400-DS is ideal for dissolution testing of extended release products that either contain and / or release small amounts of API over time.

All parameters are controlled using the 400-DS Dissolution Workstation software via the PC. Up to 4 units can be fully programmed to execute all dissolution method and sampling parameters from a single PC. Samples can be collected in pre-capped HPLC vials for direct transfer to an HPLC system. Septa are pierced using an innovative needle manifold, which lowers and raises at each sample point. Sample trays are available in 1.5 mL and 4 mL sizes.

Conventions Used in this Manual

- Items you are asked to click are in bold. For example, “Click **OK**.”

Serial Number Format

The serial number contains 10 characters and follows this syntax:

CC1234xxxx

Syntax Code	Meaning	Description
CC	Country of origin	2 alpha characters matching the required trade designation for the country of origin
12	Year of manufacture	'09' for 2009, '10' for 2010, etc.
34	Week of manufacture	'01' for week 1 to '52' for week 52

This page was intentionally left blank, except for this message.



3

Setting Up the 400-DS (G7975A)

Initial Setup	22
PC Setup	30
Initial Power Up	30
Software Installation and Setup	31

Initial Setup

Complete the following sections to initially set up the Agilent 400-DS and all other system components.

Unpacking Procedure

The Agilent 400-DS is shipped in one crate. The crate contains the apparatus and accessories.

- 1 Open the crate and check the contents for damage, which may have occurred during shipping. Shipping damage rarely occurs, but if it does contact both the carrier who delivered the instruments and the Dissolution Systems Service Department. Though claims for damage should be filed with the carrier, we can help you file a claim.
- 2 Remove the top of the crate.
- 3 Remove one side of the crate and lift the accessories box out of the crate.
- 4 Disassemble the other three sides of the crate.
- 5 Remove any cushioning material and tape. If any item is missing, contact the Dissolution Systems Service Department or your local representative for replacements.
- 6 Carefully remove the apparatus from the base of the shipping crate.

WARNING

Because of its heavy weight (approximately 59 kg), two people should lift the apparatus. Lift by holding underneath the base of the unit on each side. Support feet are located at each corner of the unit. Hands should be positioned between the support feet to ensure they are not trapped when lowering.

- 7 Place the apparatus on a clean, dry, and level section of the bench.

NOTE

The electrical connection at the back of the apparatus is the primary disconnect for the instrument. The apparatus should be positioned to allow accessibility to the power cords for easy disconnection.

Environmental Requirements for Installation

- Humidity: max relative humidity 80% for temperatures up to 31 °C decreasing linearly to 50% relative humidity at 40°C
- Indoor use only
- Pollution Degree: 2
- Installation Category: II
- Altitude: 2000 m
- Temperature: 5 °C to 40°C
- Power: 115/230 V ac, 50/60 Hz, 8 A

Main supply voltage fluctuations are not to exceed $\pm 10\%$ of the nominal supply voltage

Clearance

- 1 Prepare the area where the equipment is to be located.
- 2 Ensure a minimum clearance of 40 cm above the unit and 10 cm at the rear and on both sides of the Agilent 400-DS. Approximately 65 cm x 65 cm total bench space is required per apparatus.

NOTE

Ensure space is also available for the PC, media containers, and waste.

- 3 Orient the Agilent 400-DS and all other system components appropriately.

Parts and Accessories

Locate the following items, as applicable for your system configuration:

Item	Quantity
Agilent 400-DS Automated Apparatus 7	1
AC power cord	1
DB-9 cable	1
6-pin Mini-DIN cable	1
Sample cells	13
Cell heater jackets	13
Sample holders	13
Sample holder extraction tool	1
Roll of tubing (25 feet)	1
Sample tray with 12 rows	1
Additional sample tray for calibration	1
Magnetic plate	1
Evaporation cover with 2 guide rod stabilizer plates	1
2.5 mm Allen wrench located on rear of 400-DS	1
9/64" Allen key screwdriver	1
1.5 mm Allen key screwdriver	1
Modified temperature probe (calibration certificate included)	1
Sample vials, 12 x 32, Qty. 100	2
Sample vials, 15 x 35, Qty. 100	1
Level	1
400-DS Workstation software	1
License dongle	1
Technical Documentation CD	1

Leveling Your Apparatus

- 1 Place the level provided in the accessory kit on the vessel table.
- 2 Check the level of the instrument in the front and rear center of the unit as well as the left and right sides.
- 3 The apparatus has adjustable feet under each corner. These may be screwed in or out to raise or lower each leg. Adjust the feet so the bubble inside the level is within the circle at all four points.

Electrical Connection

NOTE

The electrical connection at the back of the apparatus is the primary disconnect for the instrument. The apparatus should be positioned to allow accessibility to the power cords.

- 1 Verify the Agilent 400-DS power switch, located in the lower left portion of the front of the instrument, is in the off position.
- 2 Connect the power cord between the receptacle on the drive unit rear panel and an outlet of the appropriate voltage.

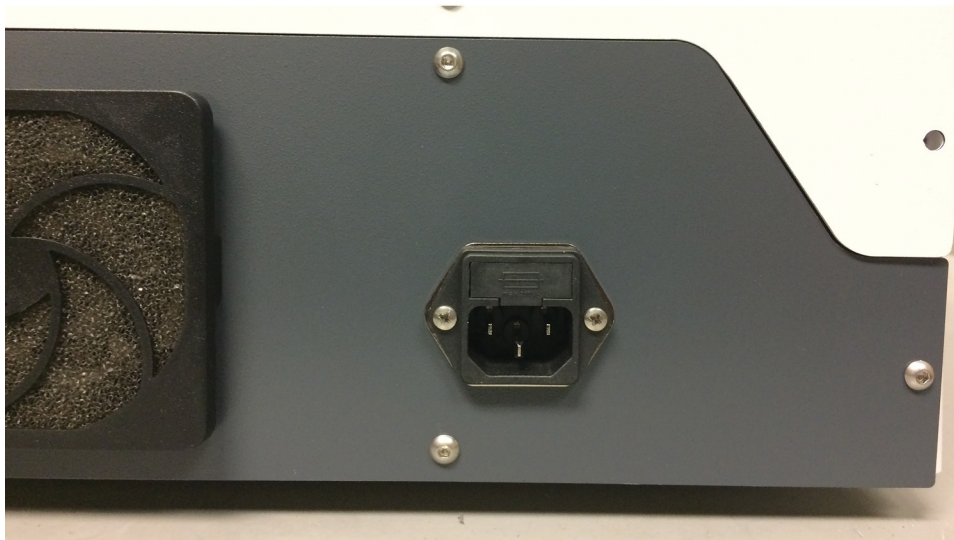
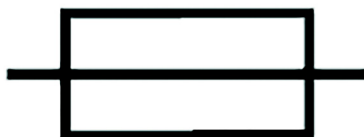


Figure 1. Power outlet located on the rear panel of the 400-DS.

WARNING

For replacing the line fuses, disconnect the power cable and replace the fuses only by the following:



T 10A H 250V (PN 2110-1004)

Consult your monitor and printer manuals for details of their individual cabling requirements.

Cable Connections

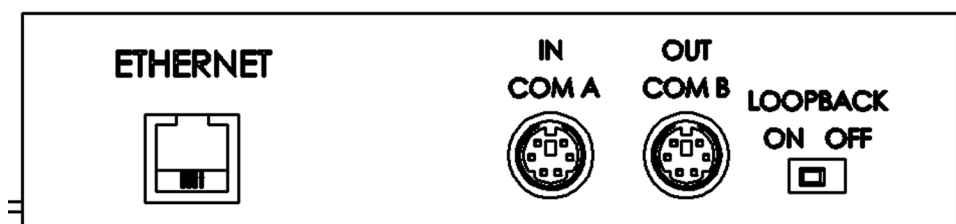


Figure 2. Diagram of cable connections

Single System Setup

- 1 Connect a cable between COM1 on the PC and the left serial port (IN COM A) on the rear of the Agilent 400-DS.
- 2 Ensure the loop back switch (located to the right of the right serial port) is switched to the left (ON).

Multiple System Setup

- 1 Connect a cable between COM1 on the PC and the left serial port (IN COM A) on the rear of the Agilent 400-DS.
- 2 Ensure the loop back switch (located to the right of the right serial port) is switched to the right (OFF).
- 3 Connect a 6-pin mini-DIN cable between the right serial port (OUT COM B) of the first apparatus and the left serial port of the second apparatus.
- 4 If this is the final Agilent 400-DS in the series, ensure the loop back switch (located to the right of the right serial port) is switched to the left (ON).

If this is not the final Agilent 400-DS in the series, ensure the loop back switch (located to the right of the right serial port) is switched to the right (OFF).

- 5 Repeat steps 3 and 4 for each additional apparatus.

Heater Jacket / Sample Holder Installation

- 1 Locate the 1.5 mm Allen wrench to install the sample heater jackets.
- 2 Install the first cell heater by placing it over the thermistor at the desired cell position of the Agilent 400-DS.
- 3 While holding the jacket in place, use the Allen wrench to tighten the screw. Do not overtighten the screw.
- 4 Insert the 2-pin plug for the heater in the appropriate location on the sample heater jacket.
- 5 Repeat steps 2 - 4 to install the sample heater jackets for the remaining cell positions.
- 6 Install the magnetic plate on the reciprocation guide rod, which is the shorter rod in the middle of the 3 rods on the Agilent 400-DS.
- 7 Tighten the bolt on each end of the magnetic plate using the 9/64" Allen wrench.
- 8 Remove the guide rod stabilizers from the bottom of the evaporation cover.
- 9 Install the guide rod stabilizers on top of the magnetic plate using the two longer guide rods for alignment.
- 10 Tighten the two bolts on each stabilizer plate using the 9/64" Allen wrench.
- 11 Locate the sample cells and sample holders provided with your instrument.
- 12 Remove the sample cells and holders from the packing material to ensure no damage occurred during shipment.
- 13 To install a sample cell in the heater jacket, slide one end through the magnetic lift plate. Use the finger tabs on the heater jacket to spread it open.
- 14 Continue to lower the sample cell until it is secure and completely covers the bottom cell seal.

CAUTION

Be careful when sliding the sample cells over the O-ring seals as damage can occur. Ensure the cell and seal are properly aligned before applying pressure.

- 15 With the magnet at the top, insert the sample holder into the sample cell.
- 16 Repeat steps 13 - 15 to install the remaining sample cells and holders.
- 17 Replace the evaporation cover and tighten the two thumb screws.

Tubing Connections

- 1 If necessary, install the tubing connectors at each location of the sample valve, rinse, and waste lines.
- 2 Connect an adequate length of tubing to each location of the sample valve, rinse, and waste lines.
- 3 Place each media / rinse / waste line into an appropriate media / rinse / waste container.

CAUTION

If one of the provided media bottle caps is used as a waste container, remove the dip tubes before attaching the waste line. If an alternative waste container is used, ensure that the waste connection is not capable of siphoning back into the instrument. The inlet tubing must never fall below the level of waste in the container.

The media bottles should be located at the same bench level as the 400-DS.

PC Setup

400-DS Workstation software requires Microsoft Windows 7 or Windows 10 to be installed on the computer. For instructions on installing the operating system, refer to the documentation supplied with Windows.

Minimum PC Configuration

Windows 7 or Windows 10	4 GB minimum RAM (8 GB recommended) 1024 x 768 pixel display resolution
10 GB free hard disk drive space	2x USB port

If the PC is not equipped with a standard serial port, a USB-to-serial converter must be purchased. Refer to the manufacturer's documentation for installation instructions.

If Agilent is installing the 400-DS Workstation for you, note that installation of the operating system is not included as part of the standard instrument installation.

Initial Power Up

- 1 Turn on the power switch on the front panel and ensure the integrated red LED illuminates.
- 2 Repeat Step 1 for each Agilent 400-DS.
- 3 Turn on the PC, monitor, and printer. Verify the PC power light illuminates and the operating system initializes.

Software Installation and Setup

NOTE

To complete the installation of 400-DS Workstation, you must log on to the PC as an administrator. Attaching the Database and Local Security Policy are included as part of the software installation procedures. The CD contains all of the requirements for an installation.

400-DS Workstation Software Installation

- 1 Insert the 400-DS Workstation CD and access the files.
- 2 Right-click **setup.exe** and select **Run as Administrator**. Follow the on-screen prompts.
- 3 When asked if you want to install the Microsoft .Net Framework, you *must* click **Yes**.
- 4 After the installation has successfully completed, install the license dongle by plugging it into an open USB port.



Figure 3. License dongle

Local Security Policy

For 21 CFR Part 11 compliance purposes, you *must* ensure that the following minimum requirements are met by your system's security policy.

NOTE

It may be necessary to coordinate the following configurations with the domain administrator.

- 1 Click **Start > Run** to run the Local Security Settings Manager. Type `secpol.msc` and press **Enter**. The Local Security Settings manager displays.
- 2 Click **Security Settings > Account Policies > Password Policy** and set the following security policy configuration:

Policy	Security Setting
Enforce Password History	3 passwords remembered
Maximum password age	30 days
Minimum password length	6 characters
Password Must Meet Complexity Requirements	Enabled

- 3 Click and set the following configuration:

Policy	Security Setting
Account lockout duration	0 minutes (infinite)
Account lockout threshold	3 invalid login attempts
Reset account lockout counter	99999 minutes

- Click **Security Settings > Local Policies > Audit Policy** and set the following configuration:

Policy	Security Setting
Audit account logon events	Success, Failure
Audit account management	Success, Failure
Audit logon events	Success, Failure
Audit policy change	Success, Failure

- Close the Local Security Settings Management screen.

Starting 400-DS Workstation

NOTE

You must log on to the computer as an administrator to set up the software and run it for the first time.

- Double-click the **400-DS Workstation** icon on the Windows desktop to start the software.
- If your system has Windows Firewall enabled, the Windows Security Alert screen displays. Click **Unblock** to enable the program.

400-DS Workstation Logon

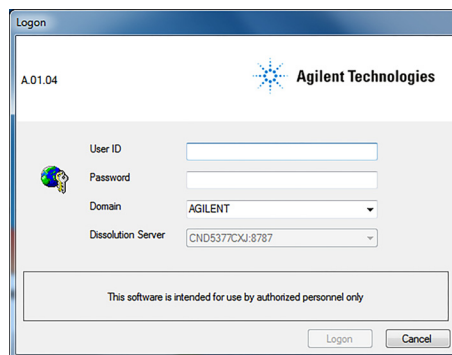


Figure 4. Logon

- 1 From the Logon screen, enter your credentials in the User ID and Password boxes.
- 2 Ensure the domain selected is appropriate and click **Logon** to initiate the software.

Adding Users to the Application

- 1 After successfully logging on to the software, click **Tools > Options**. The Configuration Dialog screen displays.
- 2 To add a user to a group, select the **Security tab** on the Configuration Dialog screen.

NOTE

To complete this section, you must be logged on as an administrator

- 3 Click **User Administration** at the bottom of the screen. The Local Users and Groups screen displays.

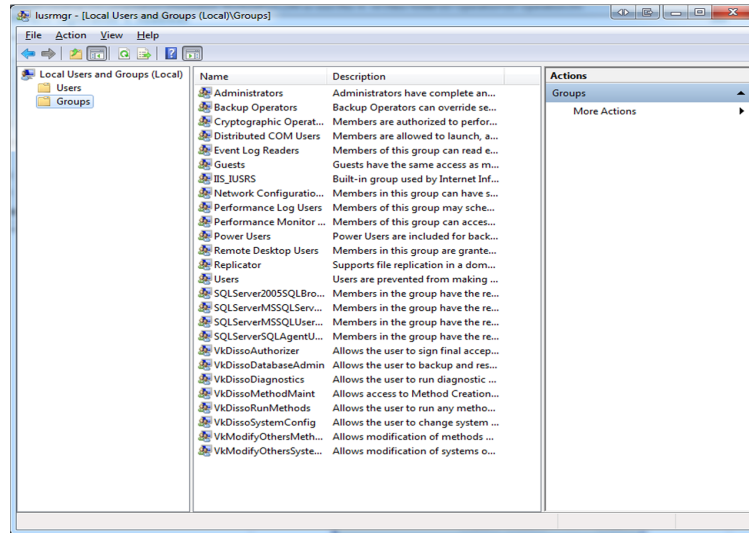


Figure 5. Local Users and Groups

- 4 Double-click the **Groups** folder to expand the list of groups.
- 5 Double-click all eight of the groups that begin with **Vk** and ensure that your username is logged in and is identified as a member of these groups.
- 6 To add a user to a group, click **Add...** from the respective group screen. The Select Users, Computers, or Group screen displays.
- 7 Enter your user identification in the empty box and click **Check Names**. Ensure your user identification and domain populate the empty field. Click **OK**.
- 8 Close the Local Users and Groups screen.
- 9 Click **OK** to close the Configuration Dialog screen.

NOTE

It may be necessary to log in again to the application to reveal the newly granted permissions for the user.

- 10 Click **Unblock** to enable the program.

This page was intentionally left blank, except for this message.



4

Operating the 400-DS

400-DS Workstation Logon	38
System Configuration	41
400-DS System Calibration	50
Manual Control/Diagnostics	56
Method Editor	60
Active Channel Selection Valves	72
Running a Method	73
Test Reports	77
Clean System	79

400-DS Workstation Logon

- 1 Double-click the **Agilent 400-DS Workstation** icon on your desktop. The Logon screen displays.

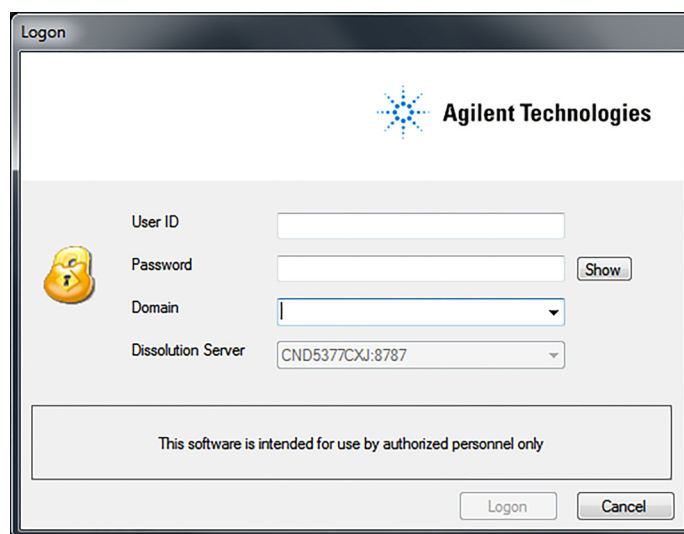


Figure 6. Logon Screen

- 2 Enter your user identification and password. Verify the domain is correct and click **Logon**. The 400-DS Workstation screen displays.

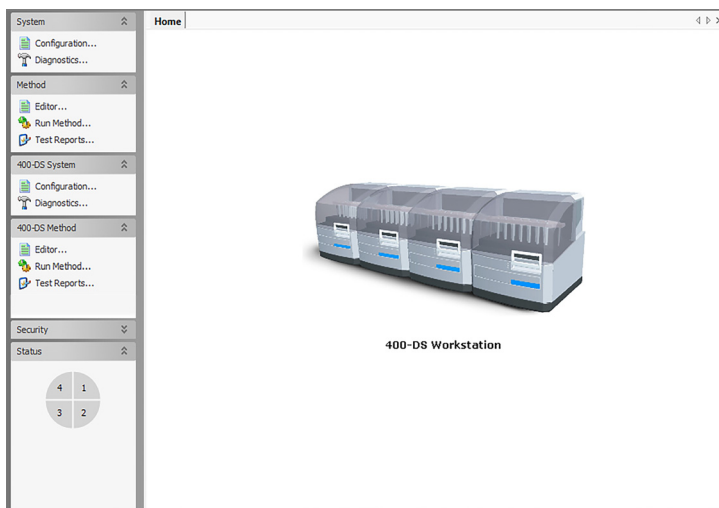


Figure 7. 400-DS Workstation Screen

NOTE

Options in the left pane may be grayed out depending on user permissions.

Following is a description of the screen options listed on the navigation bar of the 400-DS Workstation screen:

Table 1 400-DS Workstation Screen Options

Option		Description
System	Configuration	Use this option to configure the system. See "System Configuration" on page 41.
	Diagnostics	Use this option to check the diagnostics of a dissolution apparatus or accessory within the system. See "Manual Control/Diagnostics" on page 56.
Method	Editor	Use this option to create a method, change or delete method parameters, run a report of the method setup, view the method audit trail, and verify the integrity of the method. See "Method Editor" on page 60.
	Run Method	Use this option to run a method. See "Running a Method" on page 73.
	Test Reports	Use this option to run a report of the completed method. See "Test Reports" on page 77.
Security	Change User	To change the user, click Change User . The Logon screen displays. Enter the appropriate user identification and password and click Logon .
	Lock Application	To lock the 400-DS Workstation screen, click Lock Application . The 400-DS Workstation Locked screen displays. Click the lock or the link to unlock the screen. The Logon screen displays. Enter the appropriate user identification and password as applicable and click Logon.
	Audit Trail	Click Audit Trail . The Security Audit Trail screen displays. Click Show Report to display the report. Change the start date, end date, and/or user identification as desired and click Retrieve Records to change the parameters for the information displayed.
	Permissions	Use this option to view permissions assigned to the current user and the descriptions of the corresponding privileges. Click Permissions . The User Group Membership screen displays.
Status		Click one of the pie-shaped graphics under status to view the status of the corresponding dissolution system.

System Configuration

It is necessary to set up the 400-DS systems that will be used. Systems can be added, modified, and removed from the database. All system configuration activity is recorded in the system audit log. The dissolution software allows the configuration of multiple systems. A maximum of four systems can be running methods at one time on a single PC.

System configuration entails selecting the appropriate equipment and setting the communication and other physical properties of the system. Serial numbers are stored for each system to allow tracking of physical system changes.

Initial Setup of 400-DS Configuration

- 1 From the navigation bar, click **Configuration**. The System Configuration screen displays.

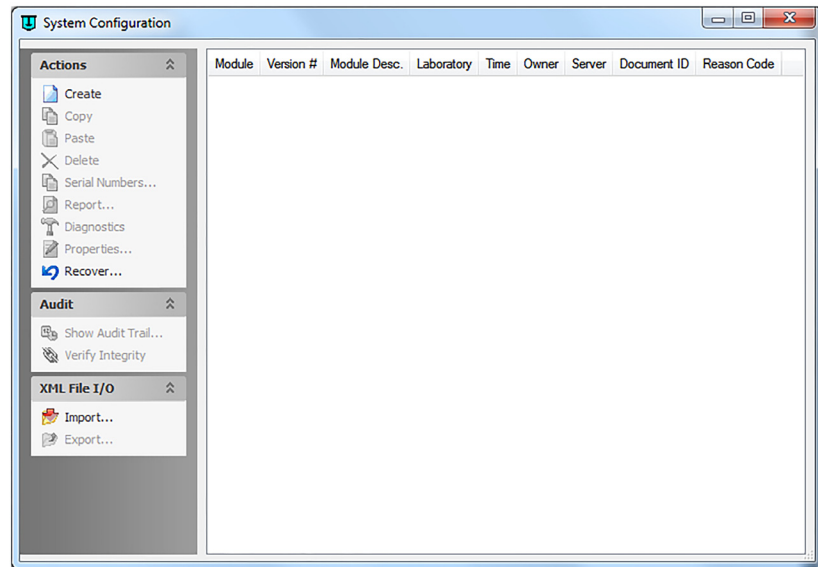



Figure 8. System Configuration Screen

- 2 Click **Create**. The System Editor screen displays (see **Figure 9**, “System Editor Screen,” on page 42).

Figure 9. System Editor Screen

Following is a description of the System Editor screen options:

Table 2 System Editor Screen 1 Options

Screen Section	Parameter	Description
System	System Name	Enter a name for your system.
	Laboratory	Enter a laboratory name.
Restrict Execution To Specified Server		Click  to display the name of the workstation or enter the name of the workstation connected to the system. Click No Restriction to allow the system to be run from any workstation. Note: The system must be physically connected.
Change Management		If applicable, select the box under Change Management in order to restrict the system editing rights to the current user. Note: This option prevents system configuration modification by anyone other than the system owner or a user with VkModifyOthersSystems privilege.

- 3 Click **Next**. The next System Editor screen displays.

The screenshot shows the 'System Editor' window with the following sections:

- Automated Apparatus 7:**
 - Serial Number: US*****
 - Cell Size: 10ml
 - Vial Size: 1.5ml
 - Syringe Size: 10ml
- Volume Calibration Coefficients:**
 - Offset:
 - Slope:
- Cell Calibration Temperatures:**
 - Manufacturer: Agilent
 - Model Name: 280-DS
 - Serial Number: MY*****
 - Setpoint Temperature 1 (C): 32.0
 - Setpoint Temperature 2 (C): 37.0
 - Setpoint Temperature 3 (C): 45.0
- Table:**

Channel	Temperature 1 (C)	Temperature 2 (C)	Temperature 3 (C)
1	32.0	37.0	45.0
2	32.0	37.0	45.0
3	32.0	37.0	45.0
4	32.0	37.0	45.0
5	32.0	37.0	45.0
6	32.0	37.0	45.0
7	32.0	37.0	45.0
8	32.0	37.0	45.0
9	32.0	37.0	45.0
10	32.0	37.0	45.0
11	32.0	37.0	45.0
12	32.0	37.0	45.0
13	32.0	37.0	45.0

At the bottom of the window are four buttons: Cancel, < Back, Next >, and Finish.

Figure 10. System Editor Screen

Following is a description of the System Editor screen options:

Table 3 System Editor Screen 2 Options

Screen Section	Parameter	Response
Agilent 400-DS	Serial Number	Click ... to display the serial numbers of all connected instruments. Highlight the appropriate serial number and click Select .
	Cell Size	Click the drop-down arrow to indicate the appropriate cell size.
	Vial Size	Click the drop-down arrow to indicate the appropriate vial size.
	Syringe Size	Click the drop-down arrow to indicate the appropriate syringe size.
Cell Calibration Temperatures	Manufacturer	Enter the manufacturer of the calibrated temperature probe.
	Model Name	Enter the model name of the calibrated temperature probe.
	Serial Number	Enter the serial number of the calibrated temperature probe.
	Set Point Temperatures (°C)	Enter the 3 set points for the cell temperatures. Enter the temperature for each cell at each set point as determined from the diagnostics screen (see "Close the System Configuration screen." on page 55).
Volume Calibration Coefficient	Offset	The current offset displays in this area.
	Slope	The current slope displays in this area.

- 4 Click **Next**. The Serial Numbers screen displays.
- 5 Enter the serial number for the sample cells, sample holders, etc. and click **Add**.
- 6 Click **Finish**.
- 7 Repeat all the sections under **"System Configuration"** on page 41 for each additional system.
- 8 Close the System Configuration screen.

Copying a System Configuration

To copy a system configuration, complete the following steps:

- 1 From the navigation bar, click **Configuration**. The System Configuration screen displays (see **Figure 8**, “System Configuration Screen,” on page 41).
- 2 Select the desired **system configuration**.
- 3 Click **Copy**.
- 4 Click **Paste**. A new system configuration displays. The description of the new system configuration is *Copy of....*
- 5 Close the System Configuration screen.
- 6 To edit the system configuration, see **“Editing an Existing System Configuration”** on page 46.

Deleting a System Configuration

To delete a system configuration, complete the following steps:

- 1 From the navigation bar, click **Configuration**. The System Configuration screen displays (see **Figure 8**, “System Configuration Screen,” on page 41).
- 2 Select the desired **system configuration**.
- 3 Click **Delete**.
- 4 Click **Yes**.
- 5 Close the System Configuration screen.

NOTE

The system configuration is never permanently deleted. It is marked as deleted in the database.

System Configuration Report

To display a report of the system configuration, complete the following steps:

- 1 From the navigation bar, click **Configuration**. The System Configuration screen displays (see **Figure 8**, “System Configuration Screen,” on page 41).
- 2 Select the desired **system configuration**.
- 3 Click **Report**. By default, the most recent report version displays.

NOTE

Alternately, you can right-click the desired **system configuration** and select **Report**. The Version Selection screen displays as a result of either of these actions.

- 4 Use the up and down arrows to indicate the desired report version and click **OK**.
- 5 The system report displays. The report can be printed, exported, searched, or verified.

NOTE

PCL6 or PPD printer drivers are required for printing a report.

Editing an Existing System Configuration

To display and edit the properties of an existing system configuration, complete the following steps:

- 1 From the navigation bar, click **Configuration**. The System Configuration screen displays (see **Figure 8**, “System Configuration Screen,” on page 41).
- 2 Select the desired **system configuration**.
- 3 Click **Properties**. The System Editor screen displays (see **Figure 9**, “System Editor Screen,” on page 42).

NOTE

Alternately, you can double-click the desired **system configuration** or right-click the desired **system configuration** and select **Properties**. The System Editor screen displays as a result of any of these actions.

- 4 Select the appropriate **tabs** and change the relevant information in the same manner that the system was created.
- 5 Close the System Configuration screen.

Recovering a Deleted Configuration

To recover a deleted configuration, complete the following steps:

- 1 From the navigation bar, click **Configuration**. The System Configuration screen displays.
- 2 Click **Recover**. A list of the deleted systems is displayed.
- 3 Select the **system** to recover and click **OK**.
- 4 Verify the system is now in the current system list.

Show Audit Trail

To display the audit trail for a system configuration, complete the following steps:

- 1 From the navigation bar, click **Configuration**. The System Configuration screen displays (see [Figure 8](#), "System Configuration Screen," on page 41).
- 2 Select the **system configuration** and click **Show Audit Trail**. The System Audit Trail screen displays.

NOTE

Alternately, you can right-click the desired **system configuration** and select **Show Audit Trail**. The System Audit Trail screen displays as a result of either of these actions.

- 3 While holding the Ctrl key, select two or more versions and click **Differences** in the navigation bar. A change report displays. The report can be printed or exported.

NOTE

PCL6 or PPD printer drivers are required for printing a report.

Verify Integrity

To verify that the system configuration has not been changed outside of the 400-DS Workstation program, complete the following steps:

- 1 From the navigation bar, click **Configuration**. The System Configuration screen displays (see **Figure 8**, “System Configuration Screen,” on page 41).
- 2 Select the **system configuration** and click **Verify Integrity**. Either the data is verified successfully or the user is directed to contact their system administrator.

NOTE

Alternately, you can right-click the desired **system configuration** and select **Verify Integrity**. The integrity of the system configuration is checked as a result of either of these actions.

- 3 Click **OK** to close the Data Verification screen.
- 4 Close the System Configuration screen.

Import/Export XML File

To use an existing system configuration from one 400-DS Workstation computer on a different 400-DS Workstation computer, you can export and import the system configuration as an XML file.

To export the system configuration, complete the following steps:

- 1 From the navigation bar, click **Configuration**. The System Configuration screen displays (see **Figure 8**, “System Configuration Screen,” on page 41).
- 2 Select the desired **system configuration**.
- 3 Click **Export**. The Version Selection screen displays.

NOTE

Alternately, you can right-click the desired **system configuration** and select **Export**. The Version Selection screen displays as a result of either of these actions.

- 4 If applicable, indicate which version to export and click **OK**. The Export System to XML File screen displays.

NOTE

Each time a system configuration is saved, a new version is created. To export a version other than the most recently saved, indicate the appropriate version number on the Version Selection screen.

- 5 Indicate the directory and file name and click **Save**. The code is saved as an XML file.

To import the XML file at another 400-DS Workstation, complete the following steps:

- 1 From the navigation bar, click **Configuration**. The System Configuration screen displays (see **Figure 8**, "System Configuration Screen," on page 41).
- 2 Click **Import**. The Import System from XML File screen displays.

NOTE

Alternately, you can right-click the desired **system configuration** and select **Import**. The Import System from XML File screen displays as a result of either of these actions.

- 3 Select the appropriate **directory** and **file name** and click **Open**. The system configuration displays on the System Configuration screen.
- 4 Close the System Configuration screen.

400-DS System Calibration

Temperature Calibration

NOTE

If a previous temperature calibration has been performed, it is necessary to reset the cell temperature values before executing this procedure:

Step 1. Click **Configuration**.

Step 2. Select the appropriate **system** from System Configuration.

Step 3. Click **Next** (from System Editor).

All channels should be set to the desired temperature (°C).

- 1 Place the rinse media line into a suitable container holding deionized water.
- 2 Ensure sample holders are installed in each sample cell.
- 3 From the navigation bar, click **Diagnostics**. The Select Item screen displays.

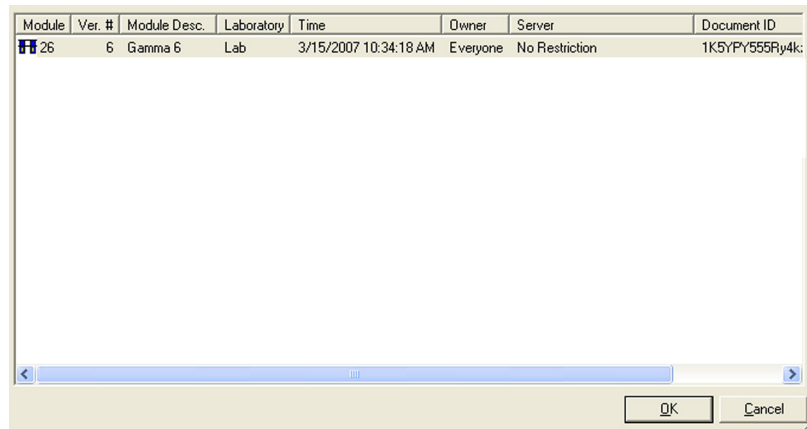


Figure 11. Select Item Screen

- 4 Select the desired **system** and click **OK**. The System Diagnostics screen displays.

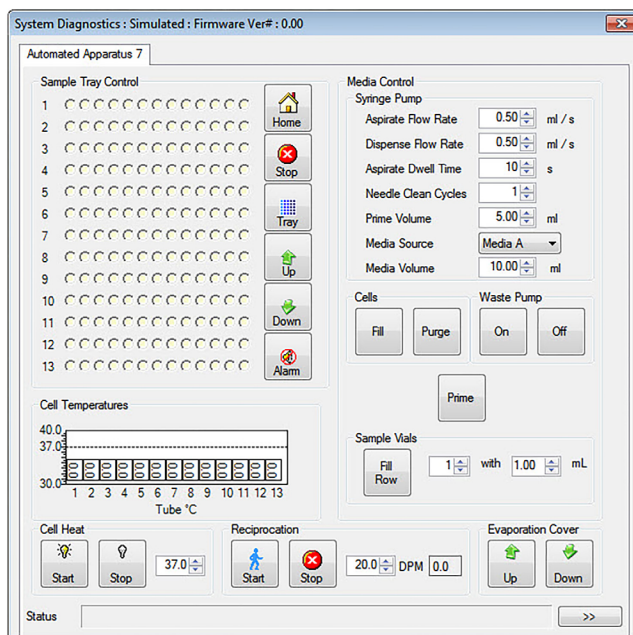




Figure 12. System Diagnostics

- 5 Ensure the media source option selected is Rinse Media. In the Cells box under Media Control, click **Fill**. Media moves into the cells.
- 6 In the Cell Heat box, set the cell temperature to the desired temperature.
- 7 Click  (start heating). DO THE SAMPLE CELLS CONTAIN LIQUID? displays.
- 8 Click **Yes**. The temperatures display in the Cell Temperature box.
- 9 Set the reciprocation rate to at least 20 DPM and click .
- 10 Allow the media to equilibrate to the set temperature.
- 11 Fill a beaker with a sufficient amount of deionized water and verify the temperature using a calibrated thermometer. Record the value.



NOTE

If using the standard QA Station probe, it will be necessary to first verify the temperature with the QA Station probe and then replace it with the Agilent 400-DS modified probe.

- 12 Using the modified temperature probe, confirm the temperature of the beaker of deionized water. Record the value. Temperatures should match ± 0.2 °C.
- 13 In the Evaporation Cover box, click the **up arrow** to raise the evaporation cover.
- 14 Remove the two thumb screws from the evaporation cover and remove the evaporation cover.
- 15 Remove the sample holder for Cell 1 and replace it with the modified temperature probe.
- 16 Close the top cover of the Agilent 400-DS. Allow the temperature to stabilize and record the temperature.
- 17 Repeat steps 14 - 16 for each additional cell.
- 18 Reinstall the sample holders and the evaporation cover.
- 19 Close the Diagnostics screen.
- 20 From the navigation bar, click **Configuration**. The System Configuration screen displays.
- 21 Select the appropriate **system** and double-click to open. The System Editor screen displays.
- 22 Click **Next**. Enter the Setpoint temperature.
- 23 Enter the temperature values for each cell (as recorded in the table beneath Step 40) into the cell calibration field.
- 24 Click **Finish**. The Reason for Change screen displays.
- 25 Enter a reason code and change description and click **OK**.
- 26 Close the System Configuration screen.

NOTE

To update 400-DS Workstation Software with the new temperature values, initiate a new method. See "Creating a Method" on page 60. and "**Running a Method**" on page 73.

- 27 From the navigation bar, click **Diagnostics**. The Select Item screen displays.
- 28 Highlight the system being qualified and click **OK**.
- 29 In the Cell Heat box, set the cell temperature.
- 30 Click  (start heating). DO THE SAMPLE CELLS CONTAIN LIQUID? displays.
- 31 Click **Yes**. The temperatures display in the Cell Temperature box.
- 32 If necessary, set the reciprocation rate and click .
- 33 Allow the media to equilibrate to the set temperature.

- 34 Record the temperatures displayed for each position.
- 35 Remove the two thumb screws from the evaporation cover and remove the cover.
- 36 Remove the sample holder for Cell 1 and replace it with the modified temperature probe.
- 37 Close the top cover of the Agilent 400-DS. Allow the temperature to stabilize.
- 38 After stabilization, record the temperature displayed on the QA Station.
- 39 Repeat steps 36 - 38 for each additional sample cell position.

NOTE

The temperatures displayed in 400-DS Workstation Software should match the modified QA Station temperature probe ± 0.2 °C to ensure accurate temperatures.

- 40 Replace the sample holders and the evaporation cover.
- 41 Click **Purge** to empty the sample cells.
- 42 Close the Diagnostics screen.

Volume Calibration

- 1 Place the tubing from the Media A valve into a suitable container with deionized water.
- 2 From the navigation bar, click **Configuration**.
- 3 Select desired configuration.
- 4 Click **Calibration**. The System Calibration screen displays.
- 5 Enter the type of media used during the calibration procedure.
- 6 Enter the density of the media used at ambient temperature.

Label 26 empty sample vials (13 x 4ml and 13 x 1.5ml) with IDs (Row 1 Vial 1 - Row 2 Vial 13). Weigh each of the vials and record their empty masses in the assigned cells in the table below. Place the sample vials in rows 1 and 2 of the sample tray. Click on the "Fill Calibration Volumes" button to fill the vials in the first row with with 4ml and the second row with 1ml of specified media. After the fill process is complete, weigh each of the vials and record their filled masses in the table below.

Media

Name Density g/ml

Balance Calibration Details

Manufacturer Model Name Calibration Date

Serial Number Calibration Due Date

System Serial Number

ID	Empty Vial Mass (g)	Filled Vial Mass (g)	Vial Volume (ml)
Row 1 Vial 1	4.7839	8.6882	3.916
Row 1 Vial 2	4.7985	8.7060	3.920
Row 1 Vial 3	4.7920	8.7050	3.925
Row 1 Vial 4	4.7912	8.6956	3.917
Row 1 Vial 5	4.8231	8.7380	3.927
Row 1 Vial 6	4.7867	8.6845	3.910
Row 1 Vial 7	4.7483	8.6586	3.922
Row 1 Vial 8	4.7595	8.6665	3.919
Row 1 Vial 9	4.8196	8.7265	3.919
Row 1 Vial 10	4.7611	8.6631	3.914
Row 1 Vial 11	4.7631	8.6578	3.901
Row 1 Vial 12	4.8344	8.7433	3.921
Row 1 Vial 13	4.8024	8.7097	3.919
Row 2 Vial 1	2.4220	3.3913	0.972
Row 2 Vial 2	2.3921	3.3756	0.987
Row 2 Vial 3	2.2003	3.1829	0.986
Row 2 Vial 4	2.4437	3.4257	0.979
Row 2 Vial 5	2.2084	3.1903	0.985
Row 2 Vial 6	2.2285	3.2029	0.977
Row 2 Vial 7	2.3778	3.3546	0.980
Row 2 Vial 8	2.2527	3.2353	0.986

Figure 13. System Calibration Screen

- 7 Label 13 empty 4 mL sample vials (Row 1 Vial 1 - Row 1 Vial 13).
- 8 Label 13 empty 1.5 mL sample vials (Row 2 Vial 1 - Row 2 Vial 13).
- 9 Weigh each of the vials and record the empty masses in the assigned cells in the table.
- 10 Place the sample vials in row 1 (4 mL) and row 2 (1.5 mL) of the sample tray.

- 11 Click **Fill Calibration Volumes** to fill the vials in the first row with 4 mL and the second row with 1 mL of deionized water.
- 12 After the fill process is complete, weigh each of the vials and record their filled masses in the table. The vial volume calculates and displays in the table automatically.
- 13 Click **OK**. The Reason for Change screen displays.

NOTE

The software performs a Least Squares Fit analysis. The volume deviation from the curve is 1.5% of the vial volume. An automated error message is generated if the calibration fails.

- 14 Enter a reason code and change description and click **OK**.
- 15 Highlight the appropriate system.
- 16 Click **Properties**.
- 17 Click **Next** and ensure the Volume Calibration Coefficients are visible.
- 18 Click **Finish**.
- 19 Close the System Configuration screen.

Manual Control/Diagnostics

To check the diagnostics of a dissolution apparatus, complete the procedures on the following pages. These procedures are performed on one system at a time. Repeat the procedures as applicable for each additional system.

- 1 From the navigation bar, click **Diagnostics**. The Select Item screen displays.
- 2 Select the desired **system** and click **OK**. The System Diagnostics screen displays.

Verifying Sample Tray Control

- 1 Click any **vessel position** corresponding to the desired row in the Sample Tray Control box. The tray moves and the indicated row raises.

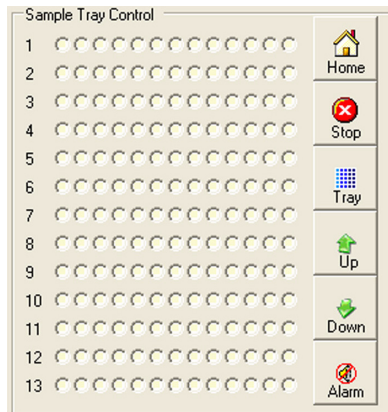








Figure 14. Sample Tray Control Box

- 2 Click  under the Sample Tray Control box. The drive unit returns to the home position.
- 3 Click  to stop the movement.
- 4 Click  to eject or retract the sample tray. Select **Ship** to place the tray in ship mode.
- 5 Click  to raise the sample tray row.

- 6 Click  to lower the sample tray row.
- 7 Click  to turn off any audible alarm.

Verifying Media Control

- 1 Under the Media Control box, set the following parameters:

Table 4 Media Control Options

Option	Description
Aspirate Flow Rate (mL/s)	Use the up and down arrows to indicate the rate at which the media is pulled into the syringe.
Dispense Flow Rate (mL/s)	Use the up and down arrows to indicate the rate at which the media is dispensed from the syringe.
Aspirate Dwell Time (s)	Use the up and down arrows to indicate the duration of time the plunger holds at the top of the stroke before dispensing. Note: The time specified is for full syringe volume. For smaller syringe volumes, it is calculated proportionately.
Needle Clean Cycles	Use the up and down arrows to indicate the number of times the unit performs the needle cleaning cycle.
Prime Volume (mL)	Use the up and down arrows to indicate the volume to be pulled before taking measurement.
Media Source	Use the drop-down arrow to indicate the source of the media.
Media Volume	Use the up and down arrows to indicate the desired volume of the media.

- 2 Under the Media Control box, click **Prime**.
- 3 Under the Cells box, click **Fill**. The media moves into the cells.

CAUTION

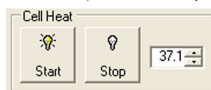
Do not click Fill more than one time without purging in between. The media will overflow the cells.

- 4 Under the Sample Vials box, use the up and down arrows to indicate the desired row number to sample into and the desired volume.
- 5 Ensure the selected row contains vials in all positions. Under the Sample Vials box, click **Fill Row**. The media moves from the cell into the sample vial.

- 6 Under the Waste Pump box, click **On** to turn on the waste pump.
- 7 Under the Waste Pump box, click **Off** to turn off the waste pump.

Verifying the Cell Temperature

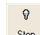
- 1 In the Cell Heat box, use the up and down arrows to set the media cell temperature to an appropriate temperature (see the sample screen below).



- 2 Click . Do THE SAMPLE CELLS CONTAIN LIQUID? displays.


CAUTION

Do not begin heating the sample cells if they are empty as this can damage the heating jacket. Ensure the each cell is filled with media before activating the heater(s).

- 3 Click **Yes**. The temperatures display in the Cell Temperature box.
- 4 As applicable, click .
- 5 Under the Cells box, click **Purge** to purge the remaining media from the tube.

Verifying Reciprocation

To set the dips per minute, complete the following steps:

- 1 In the Reciprocation box, use the up and down arrows in the box that corresponds to DPM, to set the dips per minute to a desired value and click . Dipping begins.

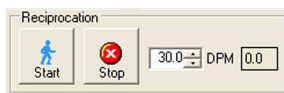


Figure 15. Reciprocation box

- 2 Click . The dipping stops.

Evaporation Cover Control

To raise the evaporation cover, click  in the Evaporation Cover box.

To lower the evaporation cover, click  in the Evaporation Cover box.

Method Editor

- 1 From the navigation bar, click **Editor**. The Methods screen displays. All test parameters are entered via the Method Editor screen.

Function	Procedure
Create a new method	See "Creating a Method" on page 60.
Copy a method	See "Copying Methods" on page 67.
Delete a method	See "Deleting Methods" on page 67.
Edit an existing method	See "Editing an Existing Method" on page 67.
Run a report of the method setup	See "Method Report" on page 68.
View the method audit trail	See "Show Audit Trail" on page 69.
Verify the integrity of a method	See "Verify Integrity" on page 69.
Import or export the method between two 400-DS Workstations	See "Import/Export XML File" on page 70.

Creating a Method

To create a new method, complete the following steps:

- 1 From the navigation bar, click **Editor**. The Methods screen displays.

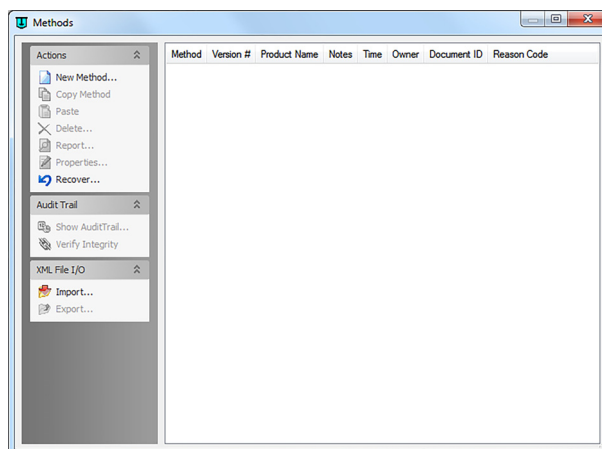


Figure 16. Methods Screen

- 2 Click **New Method**. The Method Editor screen displays.

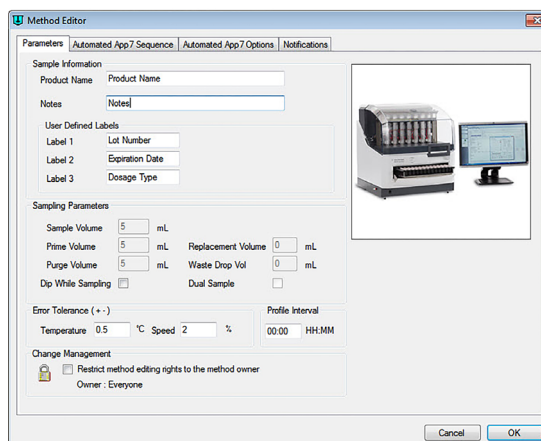


Figure 17. Method Editor Screen

3 Select the **Parameters** tab.

Following is a description of the Parameters tab options:

Table 5 Parameters Tab Options

Option	Description	
Sample Information	Product Name	Enter the name of the product.
	Notes	Enter any relevant notes regarding the sample.
	User Defined Labels	The fields in this area have default values of LOT, BATCH, and GROUP. These fields are customizable; enter the information that best serves the needs of the method parameters.
	Label 1	
	Label 2	
Error Tolerance (\pm)	Label 3	The fields in this area have default values of LOT, BATCH, and GROUP. These fields are customizable; enter the information that best serves the needs of the method parameters.
	Temperature	
	Speed	
Profile Interval		Enter the timepoint in hh:mm format at which the temperature and speed settings are recorded. Note: Profile measurements are optional. Values are always recorded at sample timepoints independent of this setting.
Change Management		If applicable, select the box under Change Management in order to restrict the method editing rights to the current user or any user with VkModifyOthersMethod.

- 4 Select the **Automated App7 Sequence** tab (see [Figure 18](#), “Automated App7 Sequence tab,” on page 63).

The screenshot shows the 'Method Editor' window with the 'Automated App7 Sequence' tab selected. The 'Parameters' sub-tab is active, showing various input fields for sample and media settings. Below these fields is a table with three rows of sequence data. The table columns are: Seq#, Sample, Sample Vol, Media Type, Media Vol, Replace Media, Tube °C, Hold Time, Dip Interval, and Speed(DPM). The first row is highlighted with a green checkmark icon, the second with a green checkmark icon, and the third with a red X icon.

Seq#	Sample	Sample Vol	Media Type	Media Vol	Replace Media	Tube °C	Hold Time	Dip Interval	Speed(DPM)
1	1	1.00	Media A	10.00	Full	37.0	00:00	001:00:00	20.0
2	2	1.00	Media B	10.00	Full	37.0	00:00	001:00:00	20.0
3	3	1.00	Media C	10.00	Partial	37.0	00:00	001:00:00	20.0

Figure 18. Automated App7 Sequence tab

Following is a description of the Automated App7 Sequence tab options:

Table 6 Automated App7 Sequence Options

Option	Description
Sample Row	Use the up and down arrows to indicate the sample tray row number for which the parameters are being set. After clicking Add to List , use the up and down arrows to indicate the next row. Note: The sample row will increment automatically from the previous timepoint.
Media	Use the drop-down arrows to indicate Media A, Media B, Media C, Media D, Media E, or Rinse media.
Sampled Volume (mL)	Use the up and down arrows to indicate appropriate volume based on your vial size.
Dissolution Volume (mL)	Use the up and down arrows to indicate appropriate volume based on the system cell size.
Cell Temperature	Enter the desired tube temperature in degrees Celsius for the first timepoint.
Hold Time (mm:ss)	Enter the desired duration for the sample holder to remain stationary prior to dipping in mm:ss format.
Dip Speed (DPM)	Use the up and down arrows to indicate the desired dips per minute (DPM) for each applicable row of the 400-DS.
Dip Interval (hhh:mm:ss)	Enter the desired duration for dipping in hhh:mm:ss format.

- 5 Select the **Automated App7 Options** tab.

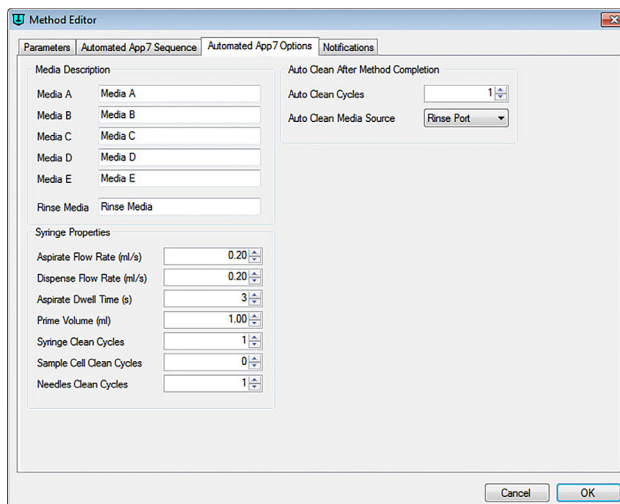


Figure 19. Automated App7 Options tab

Following is a description of the Automated App7 Options tab options:

Table 7 Automated App7 Options

Screen Section	Parameter	Description
Media Description	Media A	Enter descriptions as appropriate for the different medias.
	Media B	
	Media C	
	Media D	
	Media E	
	Rinse Media	
Syringe Properties	Aspirate Flow Rate (mL/s)	Use the up and down arrows to indicate the rate at which the media is pulled into the syringe.
	Dispense Flow Rate (mL/s)	Use the up and down arrows to indicate the rate at which the media is dispensed from the syringe.
	Aspirate Dwell Time (s)	Use the up and down arrows to indicate the duration of time the plunger holds at the top of the stroke before dispensing. Note: The time specified is for full syringe volume. For smaller syringe volumes, it is calculated proportionately.
	Prime Volume (mL)	Use the up and down arrows to indicate the volume to be pulled before taking a sample.
	Syringe Clean Cycles	Use the up and down arrows to indicate the number of times to rinse the syringe between samples.
	Sample Cell Clean Cycles	Use the up and down arrows to indicate the number of times to rinse the sample cell between samples.
	Needles Clean Cycles	Use the up and down arrows to indicate the number of times to rinse the needles between samples
Auto Clean After Method Completion	Auto Clean Cycles	Specify the number of cleaning cycles to be performed automatically at the conclusion of each test. Each cycle will flush the specified solution through the entire system - including syringes, sample cells and needles - and then dispense to waste.
	Auto Clean Media Source	Indicate the source of the solution to be pumped through the system during for the cleaning cycle. Options include Media A, B, C, D, E and Rinse Port.

- 6 Select the **Notifications** tab.
- 7 Enter appropriate address(es) to receive notification of system operation (if applicable).
- 8 Indicate which options would require notification.
- 9 Click **OK** to close the Method Editor screen.

- 10 Close the Methods screen.

Copying Methods

To copy a method, complete the following steps:

- 1 From the navigation bar, click **Editor**. The Methods screen displays (see **Figure 16**, “Methods Screen,” on page 61).
- 2 Select the desired **method**.
- 3 Click **Copy Method**.
- 4 Click **Paste**. A new method displays. The description of the new method is *Copy of....*
- 5 To change any of the parameters of the method, see “**Editing an Existing Method**” on page 67.

Deleting Methods

To delete a method, complete the following steps:

- 1 From the navigation bar, click **Editor**. The Methods screen displays (see **Figure 16**, “Methods Screen,” on page 61).
- 2 Select the desired **method**.
- 3 Click **Delete**.
- 4 Click **Yes**.

NOTE

A deleted method can be recovered at any time. See “**Recovering a Method**” on page 69.

Editing an Existing Method

To edit a method already entered on 400-DS Workstation, complete the following steps:

- 1 From the navigation bar, click **Editor**. The Methods screen displays (see **Figure 16**, “Methods Screen,” on page 61).
- 2 Click the desired **method**.
- 3 Click **Properties** on the navigation bar. The Method Editor screen displays (see **Figure 17**, “Method Editor Screen,” on page 61).

NOTE

Alternately, you can double-click the desired **method** or right-click the desired **method** and select **Properties**. The Method Editor screen displays as a result of any of these actions.

- 4 Select the appropriate **tabs** and change the relevant information in the same manner that the method was created.

Method Report

To display a report of the method parameters, complete the following steps:

- 1 From the navigation bar, click **Editor**. The Methods screen displays (see **Figure 16**, “Methods Screen,” on page 61).
- 2 Select the desired **system configuration**.
- 3 Click **Report**. The Version Selection screen displays.
- 4 If applicable, indicate which version and click **OK**.

NOTE

Each time a method is changed, a new version is created. To create a report of a version other than the most recently saved, indicate the appropriate version number on the Version Selection screen.

- 5 The method report displays. The report can be printed, exported, searched, verified, and/or signed.

Recovering a Method

To recover a method, complete the following steps:

- 1 From the navigation bar, click **Editor**. The Methods screen displays (see **Figure 16**, “Methods Screen,” on page 61).
- 2 Click **Recover**. The Select Deleted Method(s) to Recover, screen displays.
- 3 Select the **method** to recover and click **OK**.

Method Audit Trail

Once a method has completed, the results are available for review, audited modification, and electronic signature. The software maintains complete history for all runs executed on the system. Results can be previewed and printed.

Show Audit Trail

To display the audit trail for a method, complete the following steps:

- 1 From the navigation bar, click **Editor**. The Methods screen displays (see **Figure 16**, “Methods Screen,” on page 61).
- 2 Select the **method** and click **Show Audit Trail**. The Method Audit Trail screen displays.

NOTE

Alternately, you can right-click the desired **method** and select **Show Audit Trail**. The Method Audit Trail screen displays as a result of either of these actions.

- 3 While holding the Ctrl key, select two or more **versions** and click **Differences** in the navigation bar. A change report displays. The report can be printed or exported.

Verify Integrity

To verify that the method has not been changed outside of the application, complete the following steps:

- 1 From the navigation bar, click **Editor**. The Methods screen displays (see **Figure 16**, “Methods Screen,” on page 61).
- 2 Select the **method** and click **Verify Integrity**. Either the data is verified successfully or the user is directed to contact their system administrator.

NOTE

Alternately, you can right-click the desired **method** and select **Verify Integrity**. The integrity of the method is checked as a result of either of these actions.

Import/Export XML File

To use an existing method from one 400-DS Workstation on a different 400-DS Workstation, you can export and import the method as an XML file.

To export the method, complete the following steps:

- 1 From the navigation bar, click **Editor**. The Methods screen displays (see **Figure 16**, “Methods Screen,” on page 61).
- 2 Select the desired **method**.
- 3 Click **Export**. The Version Selection screen displays.

NOTE

Alternately, you can right-click the desired **method** and select **Export**. The Version Selection screen displays as a result of either of these actions.

- 4 If applicable, indicate which version to export and click **OK**. The Export Method to XML File screen displays.

NOTE

Each time a method is saved, a new version is created. To export a version other than the most recently saved, indicate the appropriate version number on the Version Selection screen.

- 5 Indicate the directory and file name and click **Save**. The code is saved as an XML file which displays in Windows Notepad.
- 6 Close the file.

To import the XML file at another 400-DS Workstation, complete the following steps:

- 1 From the navigation bar, click **Editor**. The Methods screen displays (see **Figure 16**, “Methods Screen,” on page 61).
- 2 Click **Import**. The Import Method from XML File screen displays.

NOTE

Alternately, you can right-click the desired method and select **Import**. The Import Method from XML File screen displays as a result of either of these actions.

- 3 Select the appropriate **directory** and **file name** and click **Open**. The method displays on the Method screen.

Active Channel Selection Valves

The Agilent 400-DS is equipped with channel selection valves which allow the you to isolate and shut off any unused sample cells. For example, if an experiment only required the use of eight sample vessels, Vessels 9 - 13 could be shut off by closing the red channel selection valves.

- 1 Identify the sample cells you want to disable and their corresponding valves.
- 2 Turn the red knob(s) clockwise one quarter turn.
- 3 To re-enable a valve, turn the applicable knob one quarter turn counterclockwise.
- 4 The number of active channels is programmed as part of the Method Start options just before the test begins.

CAUTION

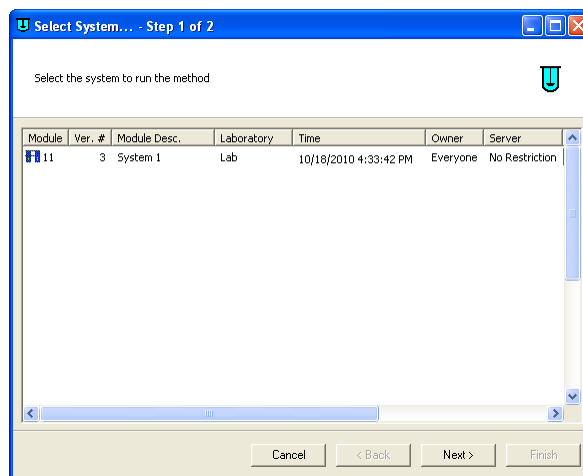
400-DS sample cells are numbered left to right. To prevent damage to the heater jackets, close the valves from right to left when less than 13 active channels are used.

CAUTION

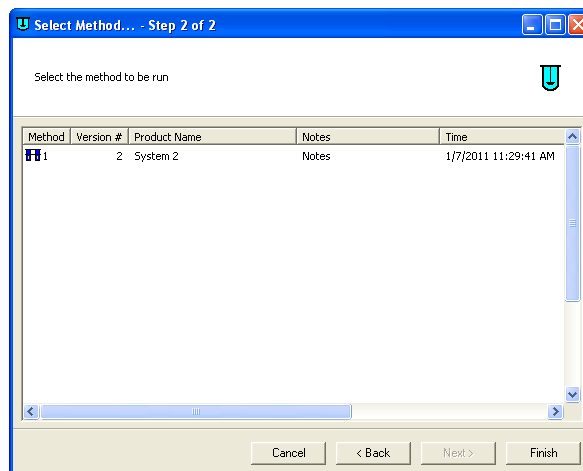
Ensure sample vials are present in positions 1 and 13 of each sample tray row if less than 13 active channels are used. This provides proper sensing, alignment, and balance for each sample tray.

Running a Method

- 1 From the navigation bar, click **Run Method**. The Select System Step 1 of 2 screen displays.



- 2 Select the desired **system** to run the method and click **Next**. The Select Method Step 2 of 2 screen displays.



- 3 Select the desired **method** to run and click **Finish**. The system status screen displays.

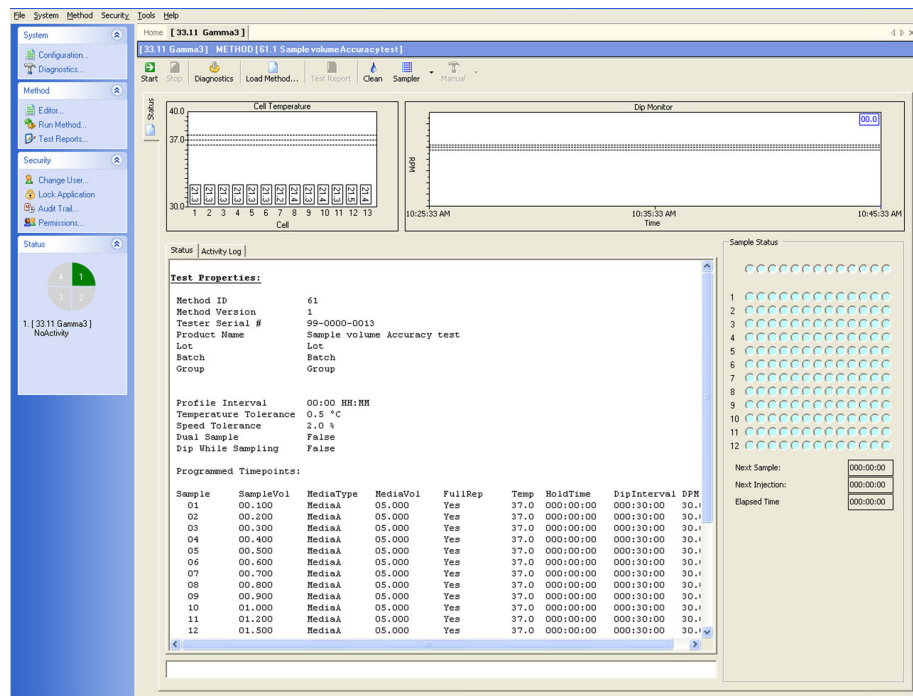


Figure 20. System status screen

- 4 If necessary, remove the evaporation cover.
- 5 Place the samples on or in the sample holder.
- 6 Insert the sample holders into the sample cells.
- 7 Reinstall the evaporation cover and tighten the two thumb screws.
- 8 Place the capped sample vials in the applicable sample rows.

CAUTION

Ensure sample vials are present in positions 1 and 13 of each sample tray row if less than 13 active channels are used. This provides proper sensing, alignment, and balance for each sample tray.

- 9 Click **Sampler > Load** to load the tray.
- 10 Fill applicable media containers.

- 11 Click **Start**. The Method Start Options screen displays (see **Figure 21**, “Method Start Options screen,” on page 75).

The screenshot shows the 'Method Start Options' dialog box. It has a title bar and a main area with several sections. The 'Sample Information' section at the top has fields for Product Name (filled with '12 Hour Test / 5mL'), Notes, Lot, Batch, and Group. Below this is the 'Active Channels' section with a numeric field set to '13'. To the right is the 'Injection Control' section with two checkboxes: 'Enable Injections' and 'Inject Only (No Sampling)', with a syringe icon. Below that is the 'Time Delayed Start' section with a date field set to '1/26/2011' and a time field set to '10:40:06 AM', accompanied by a clock icon. To the right is the 'Temperature Delayed Start' section with two checkboxes: 'Vessel Temperature Start' and 'Bath Temperature Start', with a thermometer icon. At the bottom right are 'OK' and 'Cancel' buttons.

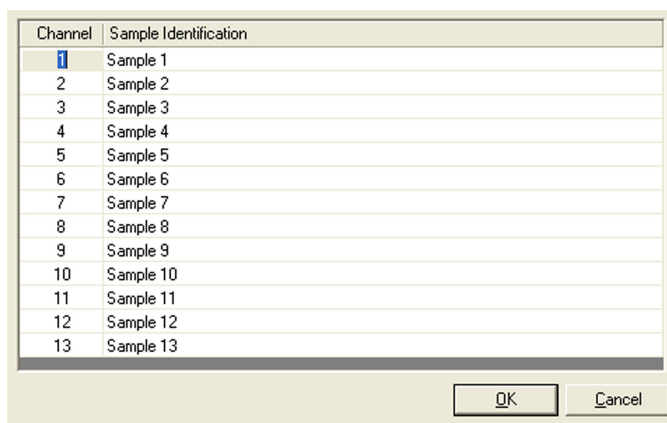
Figure 21. Method Start Options screen

Following is a description of the Method Start Options screen options:

Table 8 Method Start Options

Screen Section	Parameter	Description
Sample Information	Product Name	Enter the product name.
	Notes	Enter any appropriate notation.
	Label 1	Enter the appropriate information based on the user-defined labels (see “ User Defined Labels ” on page 62).
	Label 2	
	Label 3	
Active Channels	Enter the number of active heating channels. Note: Heating channels are counted from left to right. See “Active Channel Selection Valves” on page 72.	
Injection Control	Enable Injections	N/A
Temperature Delayed Start	Vessel Temperature Start	N/A
	Bath Temperature Start	N/A
Time Delayed Start	Use this option to program a delayed start. Enter the desired date and time to start the method.	

12 Click **OK**. The Method Sample Information screen displays.



The Method Sample Information screen is a dialog box with a table and two buttons. The table has two columns: 'Channel' and 'Sample Identification'. It contains 13 rows, numbered 1 through 13 in the 'Channel' column, with corresponding sample names in the 'Sample Identification' column. The first row is highlighted. At the bottom right, there are 'OK' and 'Cancel' buttons.

Channel	Sample Identification
1	Sample 1
2	Sample 2
3	Sample 3
4	Sample 4
5	Sample 5
6	Sample 6
7	Sample 7
8	Sample 8
9	Sample 9
10	Sample 10
11	Sample 11
12	Sample 12
13	Sample 13

Figure 22. Method Sample Information Screen

13 Enter the sample information, if applicable.

NOTE

If the method requires more than 12 sample rows, complete the following steps:

1. Upon completion of filling the 12th sample row, click **Sampler > Eject Tray**.
2. Replace the sample rows with the required amount of fresh sample vials.
3. Load the new sample tray.
4. Click **Sampler > Load New Tray**.

NOTE

Vials may be removed or replaced at any time after a sample row has been completed, provided a sample is not occurring.

14 When the method is complete, click **Test Report**. The test report displays. The report can be printed, exported, searched, verified, and/or signed.

Test Reports

To display a report of the completed method, complete the following steps:

- 1 From the navigation bar, click Test Reports. The Test Report Selection screen displays.

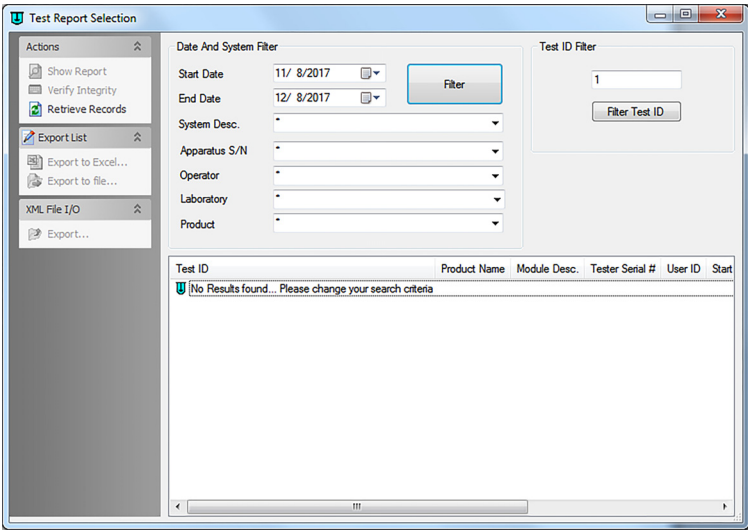


Figure 23. Test Report Selection screen

NOTE

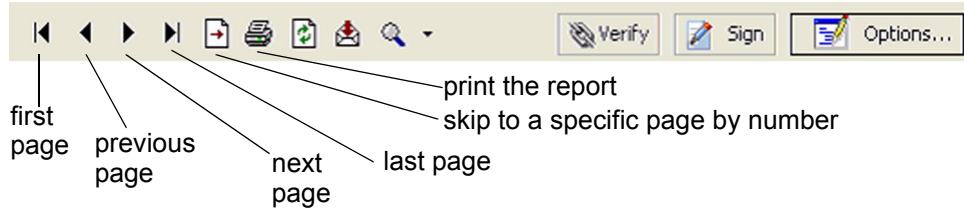
By default, the test report dialog only shows tests done in the previous week. To view tests completed earlier than the previous week, click the **Start Date** and/or **End Date** drop-down arrows to indicate date range for a test report selection and click **Filter Date and System**.

Additionally, you can select a **system name** from the System combo box to restrict the selection to a particular system.

To restrict the selection to a particular test identification, enter the value in the Test ID Filter box and click **Filter Test ID**.

- 2 Click a **test** in the list to select it.

- 3 From the navigation bar in the Test Report Selection dialog, click **Show Report**. The report for the selected test displays.
- 4 Using the buttons on the report toolbar, you can do the following:




Electronic Signatures

When the user is satisfied with the results, the results can be electronically signed. The software allows multiple electronic signings of a set of results. Each signing is accomplished using the signature dialog box shown. The user authenticity is determined by testing the user identification and password against the Windows security database.

Electronic signatures are permanently linked to the results. The software always requires the signature to be executed using all the signature components. Any attempts to sign a set of results using an invalid user identification, password, or any combination thereof that is incorrect is automatically recorded to the system audit trail.

Clean System

From the system status screen, click . All instruments initialize. Media is pulled from the Rinse Media Container and cells are filled and expelled through the valves. The cells are aspirated and purged.

NOTE

The 400-DS should be flushed / cleaned as soon as possible after a test. Multiple clean cycles may be necessary to properly rinse the system.

This page was intentionally left blank, except for this message.



5 Troubleshooting and Maintenance

Preventive Maintenance 82

Troubleshooting 84

Preventive Maintenance

WARNING

Disconnect power from the 400-DS before servicing.

WARNING

The 400-DS contains electrical circuits, devices, and components operating at dangerous voltages. Contact with these circuits, devices, and components can cause death, serious injury, or painful electric shock.

Preventive maintenance needs may vary depending on frequency of instrument usage.

Calibration

Agilent recommends the 400-DS be calibrated at six-month intervals. However, based on your own SOPs, this can be extended for up to a year. Contact the Agilent Service Department to schedule preventive maintenance and qualification. The Agilent Service Department can be reached at 1.800.227.9770.

Routine Maintenance (to be Performed Between Each Dissolution Method)

- Execute the Clean System cycle using the 400-DS Workstation software (“**Clean System**” on page 79). Ensure an appropriate cleaning solution is in place to thoroughly rinse the media lines and sample cells.

Sample Holders

- Clean sample holders thoroughly after each use and place them in an ultrasonic bath for a few minutes, if necessary.
- When using corrosive materials such as hydrochloric acid, surfactants, or medium containing salts, be sure to rinse the holders with deionized water immediately after use. Dry completely with a soft towel or cloth (preferably lint-free).
- Do not clean the sample holders with abrasive cleansers or cloths. Use deionized water whenever possible. If you must use cleanser or solvent, be sure that it is as mild as possible, non-abrasive, and fully compatible with fluorocarbons and stainless steel before use.
- It is suggested that the sample holders are cleaned by hand. Use of a laboratory dishwasher is NOT recommended.
- Handle with care at all times.
- Store the sample holders properly between each use.

Troubleshooting

Obtaining Warranty and Other Services

To place a service order (warranty or other services), please contact your local Customer Care Center. Contact information can be found at www.agilent.com under your country using the Contact Us link. Place your service request using the displayed phone number or E-mail address.

In This Book

- Chapter 1 Safety
- Chapter 2 Introduction
- Chapter 3 Setup
- Chapter 4 Operation
- Chapter 5 Troubleshooting

www.agilent.com

© Agilent Technologies, Inc. 2018

70-9064 Rev D March 2018

