

Magnis/MagnisDx NGS Prep System

Quick Reference Guide



K1007A MagnisDx NGS Prep system: For In Vitro Diagnostic Use

G9710A Magnis NGS Prep system: For Research Use Only. Not for use in diagnostic procedures.

Before You Begin 2

Safety Precautions 3

Environmental Requirements 6

Instrument Components 8

Starting Up the Magnis/MagnisDx NGS Prep System 11

Turn on the instrument 11

Log in to the system 11

Creating User Accounts 12

About user access levels 12

Add new user accounts 12

Decontaminating the System 13

Running Protocols 14

1) System prep 14

2) Reagent prep 14

3) Protocol set up and run 15

4) Collection and clean up 16

Before You Begin

- Intended Use

The MagnisDx NGS Prep System is an automated liquid handling system for next generation sequencing library preparation and/or target enrichment of human nucleic acid samples.

The MagnisDx NGS Prep System is to be used only by operators trained in laboratory techniques and procedures.

The customer is responsible for validation of assays and compliance with regulatory requirements that pertain to their procedures and uses of the instrument.
- Principle of the Procedure

The Agilent Magnis/MagnisDx NGS Prep System is a liquid handler that offers start-to-finish automation of library preparation and target enrichment protocols for next generation sequencing (NGS). The starting material is total RNA or fragmented genomic DNA (gDNA) purified from a cell or tissue sample, blood sample, or formalin-fixed paraffin-embedded (FFPE) sample. The end result is a target-enriched DNA library ready for sequencing.

The hardware components make up the instrument. See **“Instrument Components”** on page 8 for a list of these components.

The instrument is controlled through the software component of the system, which is displayed and operated through the LCD touchscreen.

Magnis reagents must be vortexed and centrifuged as described in the target enrichment protocol for the reagents to ensure optimal performance.

Limitations of Use

The Magnis/MagnisDx NGS Prep System has been validated for use with Agilent Magnis NGS kits.

Materials Provided

Table 1 Materials provided with the Magnis/MagnisDx NGS Prep System

Materials
Instrument with preloaded touchscreen software
Power cord
Functional test certificate

Safety Precautions

The Magnis/MagnisDx NGS Prep System is designed for safe operation when it is used in the intended manner. Using the system for purposes other than those intended may impair those safety protections.

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.

Installation

CAUTION

The system must be installed by an Agilent engineer or Agilent authorized service provider.

WARNING

Do not attempt to manually lift the instrument. To move the instrument, use an automated forklift or lift table that can lift at least 100 kg (220 lbs).

WARNING

When adjusting the placement of the instrument on the laboratory bench, take care to avoid pinching hazards on the instrument.

CAUTION

Retain the locking jigs that secure the gantry during shipment. Anytime the instrument is moved, secure the gantry using the jigs.

Electrical

Standard electrical safety precautions should be applied, including the following:

WARNING

Per North American and IEC requirements, install the instrument in a location where branch circuit protection is available in the power network.

Safety Precautions

WARNING

Install the instrument with the Agilent supplied power cord that is compatible with the electrical outlets in your region. Do not substitute with a power cord from another source.

WARNING

Install the instrument away from any flammable sources.

CAUTION

Install the instrument in a location where the power cord can be easily reached for quick disconnection of the power supply.

CAUTION

Make sure that the ventilation slots of the instrument are free of any obstruction. Maintain 10 cm (4 inches) of cleared space on each side of the instrument and 18 cm (7 inches) of cleared space at the back of the instrument.

CAUTION

Connect the power cord to a wall output that provides 100–240 VAC, 50/60 Hz, 1000 W.

CAUTION

Make sure that proper voltage is supplied before turning on the instrument for the first time.

WARNING

Connect the instrument to a grounded socket. Do not operate the instrument from a power outlet that has no ground connection.

WARNING

Do not connect the instrument to the same circuit as other devices with a high current draw (e.g., freezers, centrifuges). If possible, connect the instrument to an independent or dedicated AC circuit.

WARNING

Do not touch any switches or outlets with wet hands.

CAUTION

Before disconnecting the power cord, turn off the instrument at both the power button on the front and the power switch on the back.

WARNING

Unplug the instrument before you clean any major liquid spills and before servicing any of the electrical or internal components.

Safety Precautions

WARNING

Do not operate the instrument in a hazardous or potentially explosive environment.

CAUTION

Do not service the electrical components unless you are qualified to do so.

Fluids and Reagents

CAUTION

Magnis reagents must be vortexed and centrifuged as described in the target enrichment protocol prior to initiating a Magnis run. Never process reagents not intended for use with the Magnis/MagnisDx NGS Prep System.

CAUTION

During deck setup, make sure that all labware is positioned flat on the designated deck platform or fully seated in the appropriate labware holder. Improperly positioned labware can cause low or no final library yield for some or all samples.

CAUTION

When preparing for a Magnis run, and using critical input DNA or RNA samples with limited quantity, verify that there is a sufficient quantity of sample to complete at least two Magnis runs.

WARNING

Observe the relevant safety regulations when handling pathogenic material, radioactive substances or other substances hazardous to health.

WARNING

Do not submerge the instrument in any liquid.

Danger of Ultraviolet (UV) Light Exposure

The instrument door and side panels are not UV-transparent, therefore exposure to UV light is minimal. However, the following precautions are still needed.

WARNING

During decontamination of the instrument deck with UV light, do not look directly or indirectly at the UV light source.

WARNING

Always perform decontamination with the instrument door closed and locked. The instrument door is programmed to remain locked while the UV light is on.

Environmental Requirements

WARNING

Replacement UV tubes must be provided by Agilent and must be installed by an Agilent engineer or Agilent authorized service provider.

Danger of Burns

WARNING

During protocol runs, the thermal block and other components of the thermal cycler module quickly attain temperatures of greater than 50°C. To ensure safe operation, the instrument door must remain closed during runs. The instrument door is programmed to remain locked while protocol runs are in progress.

CAUTION

Use only Agilent materials (plates, adhesive seals, foils, mats) intended for use on the Magnis/MagnisDx NGS Prep System. These materials are sufficiently temperature-stable (up to 120°C).

Electrostatic Discharge

CAUTION

The Magnis/MagnisDx NGS Prep System is static sensitive. Electrostatic discharges greater than 8000 volts may interfere with the normal operation of the USB ports on the instrument. Handling precautions are required when working in high static environments. Wear a grounded wrist strap and take other anti-static precautions prior to making contact with the instrument in high static environments. ESD STM5.1-1998 Class 3B.

Environmental Requirements

CAUTION

Operating temperature

Keep the ambient temperature between 15°C and 25°C (53.6°F to 77°F).

CAUTION

Operating humidity

Keep the humidity levels between 30% and 70% non-condensing.

CAUTION

Altitude

Maximum altitude for instrument operation is 2000 m (6562 feet)

Installation Requirements

Review the “**Electrical**” safety precautions on [page 3](#) for power supply requirements that impact safe operation of the instrument.

Place the instrument in a post-PCR laboratory environment.

CAUTION

Instrument is for indoor use only.

CAUTION

The laboratory humidity levels must be between 30% and 70% non-condensing. Operating the system in humidity levels outside of this range can impact performance.

CAUTION

Do not place the instrument in proximity to other lab equipment that is sensitive to vibrations or that generates vibrations during use. Proximity to lab equipment that generates vibrations can impact performance.

CAUTION

Maintain the following minimum distances for cleared space around the instrument.

Instrument sides: 10 cm (4 inches) on each side to allow for proper ventilation from the side vents.

Back of the instrument: 18 cm (7 inches) at the back to allow for proper ventilation from the back vent.

Front of the instrument: 5 cm (2 inches) in front to avoid unintentional contact with the power switch.

Top of the instrument: 111 cm (44 inches) above to allow for opening the door.

CAUTION

Install the instrument in a location where the power cord can be easily reached for quick disconnection of the power supply.

WARNING

Install the instrument with the Agilent supplied power cord. Do not substitute with a power cord from another source.

WARNING

Install the instrument away from any flammable sources.

CAUTION

After installation is complete, avoid moving the instrument or making any adjustments to the placement of the instrument as it will disturb some of the settings made by the Agilent engineer or service provider, resulting in an additional service visit.

Instrument Components

Front and sides of instrument

- Instrument door
- Waste bin drawer containing disposable tip waste bin
- LCD touchscreen for display of the firmware software
- Power button
- ports (2)
- Side vents (1 on each side)



Figure 1 Instrument front, door closed

Back of instrument

- Power switch
- Ethernet port
- Inlet for power cord
- Back vent



Figure 2 Instrument back

Instrument Components

Instrument interior

- Instrument deck consisting of the following modules:
 - Thermal cycler module for incubation and PCR steps
 - Heater/shaker/magnet (HSM) module for various processing steps
 - Chiller module for reagent storage
 - Tube holders for liquid reagents (6 total)
 - Platforms to hold tip boxes (4)
 - Platform to hold Beads/Buffers Plate
 - Opening to tip waste bin
 - LED indicator lights (2)
 - Micropipettor for liquid transfer
 - Barcode scanner for labware verification and sample tracking
 - Gantry for positioning of the micropipettor and barcode scanner
 - UV tube for decontamination of surfaces of the instrument deck with UV light

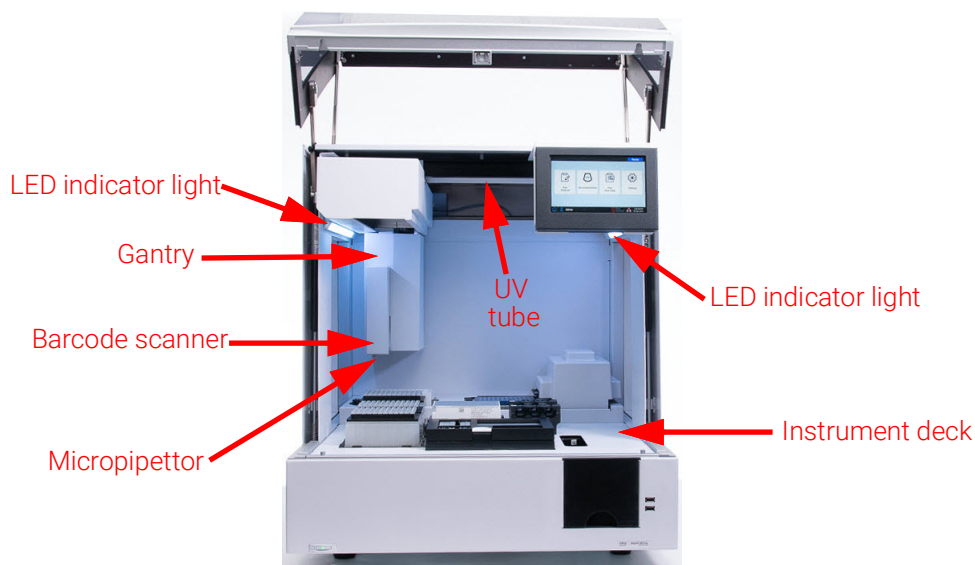


Figure 3 Instrument interior

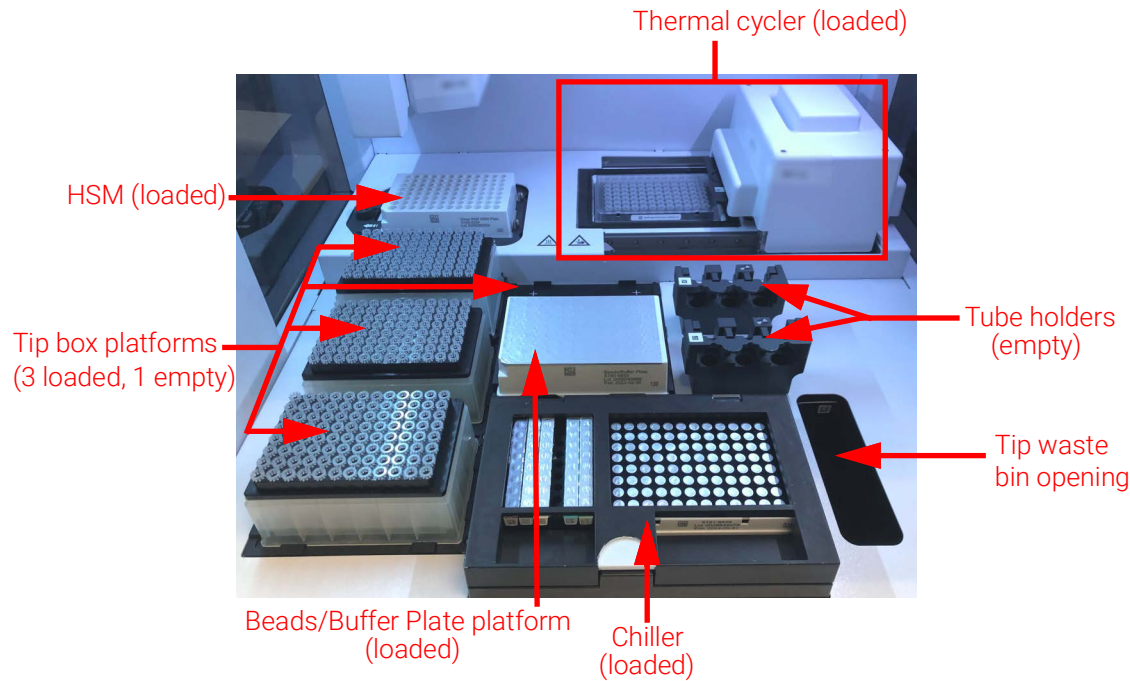


Figure 4 Instrument deck

Indicator Lights You can quickly and easily check the status of the instrument based on the color of the LED indicator lights that illuminate the entire plate filling area.

Table 2 Colors and descriptions of the LED indicator lights in the plate filling area

LED Color	Instrument Status	Description
White	Ready	The lights are illuminated white when the system is idle but the door is open, when the system is performing auto-teaching or teach-point verification, and when a user is setting up a protocol run.
Blue	Ready	The lights are illuminated blue anytime the system is idle and the door is closed, including at completion of a protocol run. The lights are also blue when the system is performing diagnostic tests.
Green	Running	The lights are illuminated green when the system is running a protocol.
Red	Error	The lights are illuminated red when the system has encountered an error. Check the touchscreen for an error message containing further details.

WARNING During UV decontamination, the indicator lights are off and the instrument deck is illuminated by the UV light. Do not look directly at the UV light.

Starting Up the Magnis/MagnisDx NGS Prep System

Turn on the instrument

The power button on the front of the instrument is to toggle the instrument on and off.

- 1 Press the power button on the front of the instrument.

The light on the power button turns green, the instrument turns on, the LED indicator lights inside the instrument illuminate, and the software launches on the touchscreen.

If pressing the power button fails to turn on the instrument, verify that the power switch on the back of the instrument is in the ON position.

Log in to the system

If you do not yet have your own personal user account, use the user name and password provided by the Agilent engineer or service provider who installed your system.

- 1 Access the Login screen of the software.

The Login screen opens automatically after the turning on the instrument.

If another user is already logged in, press the user name at the bottom of the screen then press **Log Out**. The previously logged in user is logged out and the Login screen opens.

- 2 Type the user name and password for your account into the provided fields.

The Agilent engineer or Agilent authorized service provider creates a user account during system installation that has advanced user access.

- 3 Press **Login**.

You are now logged in to the software.

Stand by as the system performs the series of preparatory activities described in [Table 3](#).

Table 3 Preparatory steps

Step	Description
Hardware initialization	During hardware initialization, the system returns all motorized parts (i.e., the gantry, HSM module, and thermal cycler) back to their home positions. Additionally, the system checks for the presence of any tips on the micropipettor and, if necessary, disposes of those tips in the tip waste bin.
Instrument Health Check (IHC)	The system performs an IHC each time it is powered on (after login) and each time a protocol run is initiated. The checks performed during the IHC ensure that the hardware is functioning within specifications.
Teach-point verification	<p>Auto-teaching is a process in which the system locates and records the positions of markers (called teach points) that are printed on the instrument deck. Auto-teaching ensures that, during a run, the micropipettor is accurately aligned with the tubes or wells at each deck position.</p> <p>You can configure the system settings to include teach-point verification as part of the initial IHC that is performed after powering up the instrument. During teach-point verification, the current teach-point locations are compared to those that were previously recorded during the most recent auto-teaching run to ensure that the values are sufficiently within range of each other. If the values do not fall within the expected range, the system prompts you to rerun auto-teaching.</p>

Creating User Accounts

About user access levels

The access level assigned to a user account – *Standard* or *Advanced* – determines the user's access to certain software settings and functions. The table below summarizes the differences in privileges between the two access levels.

Table 4 Actions permitted for Standard and Advanced user account

Action	Permitted by Standard users?	Permitted by Advanced users?
Set up and run protocols	Yes	Yes
View data from their own protocol runs	Yes	Yes
Update the number of PCR cycles during run setup	No	Yes
View data from protocol runs of other users	No	Yes
Edit user accounts of other users	No	Yes
Edit chiller temperature setting	No	Yes
Install firmware and protocol updates	No	Yes
Delete diagnostic reports	No	Yes
Change default protocol version	No	Yes
Bypass reagent expiration error	No	Yes*

* Only permissible on the G9710A Magnis NGS Prep System.

Add new user accounts

Each user who uses the system needs an account.

- 1 From the Home screen, press **Settings**.
- 2 Press **User Management**.
The User Management screen opens listing the available user names and corresponding access levels and statuses.
- 3 Press **Add**.
The Add New User screen opens.
- 4 In the User Name field, enter a user name for the new account.
User names can contain a combination of letters and numbers, but numbers are only permitted at the end of the user name (e.g., abc123). Do not include special characters in the user name.
- 5 Next to **Access Level**, select the access level (Standard or Advanced) for the new user. The default selection is *Standard*.
See **Table 4** for a summary of the differences between the two access levels.
- 6 In the Password and Confirm Password fields, enter a password for the account.
- 7 Press **OK** to save the user account.
The Add New User screen closes and you are returned to the User Management screen. The new user name appears in the list on the User Management screen.

Decontaminating the System

The Magnis/MagnisDx NGS Prep System includes a UV tube that can be used to decontaminate the surfaces of the instrument deck. The *quick cycle* is a 30-minute decontamination. Agilent recommends running the quick cycle before each protocol run. The 2-hour *extended cycle* is recommended in the event of a spill or leak on the deck surfaces.

Instructions for running a quick cycle are provided below. For instructions on running an extended cycle, see the full-length *Magnis/MagnisDx NGS Prep System User Guide* (publication K1007-90000).

WARNING Do not look directly at the UV light while decontamination is in progress.

Run a quick cycle

Agilent recommends running the decontamination quick cycle before each protocol run.

- 1 Verify that the instrument deck is cleared of all labware, then make sure that the instrument door is closed.
- 2 From the Home screen, press **Decontamination**.
The Decontamination screen opens.

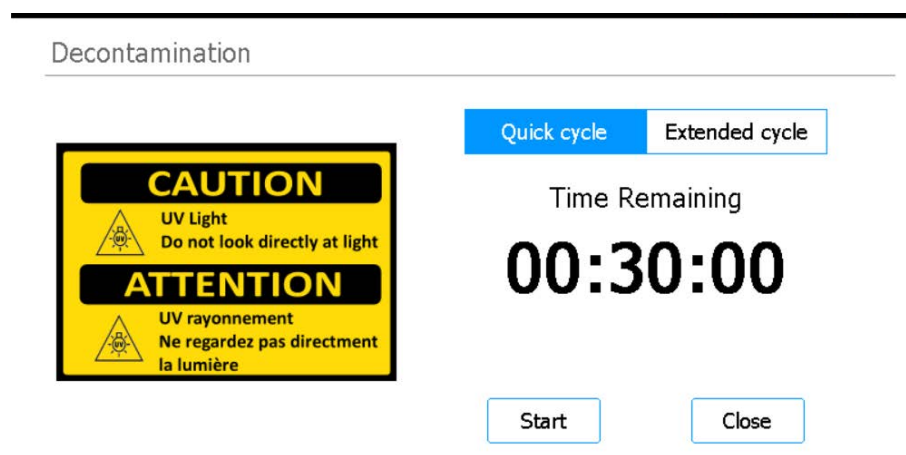


Figure 5 Decontamination screen with Quick cycle selected

- 3 At the top of the screen, select **Quick cycle**, as shown in **Figure 5**
- 4 Press **Start**.
The decontamination cycle begins and the screen displays a countdown of the time remaining. At the completion of the cycle, the UV light turns off and the instrument remains idle. If needed, press **Abort** at any time to stop the decontamination cycle.

Running Protocols

- 1) System prep** Prepare the Magnis/MagnisDx NGS Prep System so that it is ready to run a protocol.
- 1 Verify that the instrument deck is cleared of all labware from previous runs.
 - 2 Turn on the instrument and close the instrument door.
See **"Turn on the instrument"** on page 11 for instructions.
 - 3 At the Login screen, enter the credentials for your user account.
Each time the system is turned on, it performs a series of start up activities (see **Table 3** on page 11) immediately after login. These activities may take several minutes to perform. Make sure the instrument door remains closed for the duration of these activities. After completion of the start up activities, the software opens to the Home screen and the system's LED light is blue, indicating that it is ready for use.
 - 4 (Optional) Perform the 30-minute UV decontamination quick cycle to decontaminate the surfaces of the instrument deck. See **"Decontaminating the System"** on page 13.
You can begin **"2) Reagent prep"** while the decontamination is running.
For instructions on manually decontaminating the surfaces of the instrument deck, consult the full-length *Magnis/MagnisDx NGS Prep System User Guide* (publication K1007-90000).

- 2) Reagent prep** Prepare the samples, target enrichment reagents, and other materials needed for the protocol run according to the user manual for your specific Magnis Target Enrichment Kit. The user manual contains details on the materials required to run a protocol and instructions on loading your DNA samples into the Magnis Sample Input Strip.

The image below describes the orientation of the Magnis Sample Input Strips. Make sure to track the locations of your samples as you load them into the strip.

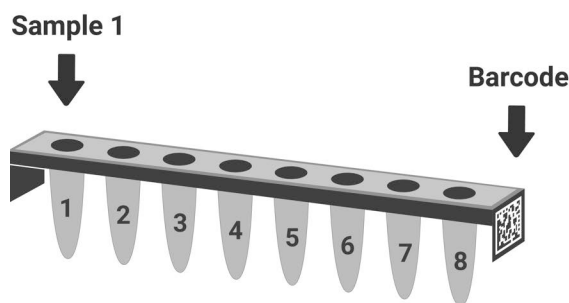


Figure 6 Orientation of the Magnis Sample Input Strips

CAUTION

Avoid touching the foil seals of sample strips, reagent strips, and reagent plates, even using gloved hands. Any contaminants deposited on strip or plate foil seals can be introduced into samples during the Magnis liquid-handling steps.

CAUTION

Magnis reagents must be vortexed and centrifuged as described in the target enrichment protocol prior to initiating a Magnis run.

CAUTION

During deck setup, make sure that all labware is positioned flat on the designated deck platform or fully seated in the appropriate labware holder.

CAUTION

Do not add any writing or labels to the strips, plates, and other labware that may obscure the barcodes.

3) Protocol set up and run

For detailed instructions on running a protocol for your particular sample type, refer to the user manual for your Magnis Target Enrichment Kit.

- 1 From the Home screen, press **Run Protocol**.

The system locks the instrument door and performs an IHC. At completion of the IHC, the Enter Run Info step opens.

- 2 Complete the Run Setup steps and start the protocol run, as instructed in the user manual for your Magnis Target Enrichment Kit.

The software walks you through the individual Run Setup steps (**Figure 7**), starting with the Enter Run Info step. Press the forward and back arrow buttons to navigate through the steps. Because the steps vary depending on the type or target enrichment you are running, you will need to refer to the instructions and screen images provided in the user manual for your Magnis Target Enrichment Kit.

After you start the run, the system's LED light turns green, indicating that a protocol run is in progress. The LED changes from green to blue upon completion of a the run.

If the system encounters an error during the protocol, the LED light changes to red.

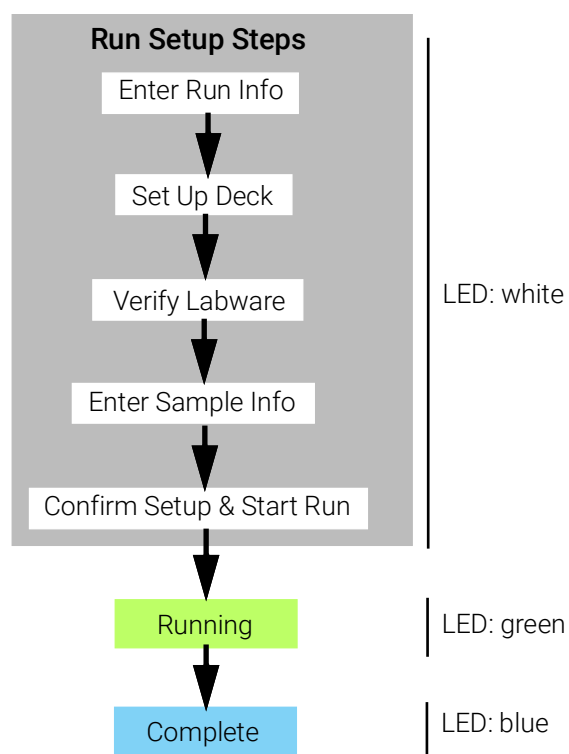


Figure 7 Protocol workflow, with corresponding color of LED status indicators lights

4) Collection and clean up

At the completion of the run, the system keeps the prepared library solutions in the PCR plate on the thermal cycler module, which is held at 12°C for up to 72 hours. Collect the samples within that 72-hour period.

- 1 Wait until the touchscreen indicates that the run is complete. Press **OK** when you are ready to collect the library samples from the instrument.

The system transfers the libraries from the PCR plate to the green library strip tube in the chiller. Allow the instrument-mediated transfer process to complete before continuing.

- 2 At the completion of the transfer process, open the instrument door and collect the final library samples (i.e., the green library strip tube) from the chiller module.
- 3 Re-seal the wells of the strip tube using a fresh foil seal strip, then store at the temperature recommended in the user manual for your Magnis Target Enrichment Kit.
- 4 If the optional pre-capture library QC samples were collected for the run, remove the blue QC sample strip from the chiller module. Process and store the samples as recommended in the user manual for your Magnis Target Enrichment Kit.
- 5 Remove all remaining consumables from the instrument deck and dispose of them according to local guidelines. Close the instrument door.
- 6 Log out of the software or turn off the instrument.

To log out, press the user name at the bottom of the screen then press **Log Out**.

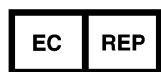
To turn off the instrument, press the power button on the front of the instrument.

Legal Manufacturer



Agilent Technologies Singapore (International) Pte Ltd.
No. 1 Yishun Avenue 7, Singapore 768923
Manufactured at:
Agilent Technologies LDA Malaysia Sdn. Bhd.
Bayan Lepas, Free Industrial Zone 11900 Penang, Malaysia
www.agilent.com

Authorized Representative for the European Union



Agilent Technologies Denmark ApS
Produktionsvej 42
2600 Glostrup, Denmark

Authorized Representative for the United Kingdom



Agilent Technologies LDA UK Limited
5500 Lakeside, Cheadle Royal Business Park
Cheadle, Cheshire, SK8 3GR, UK

Authorized Representative for Switzerland



Agilent Technologies (Schweiz) AG,
Lautengartenstrasse 6
4052 Basel, Switzerland

European Union and Switzerland Importer



Agilent Technologies Deutschland GmbH
Hewlett-Packard-Str. 8
76337 Waldbronn, Germany



Agilent Technologies (Schweiz) AG,
Lautengartenstrasse 6
4052 Basel, Switzerland

Agilent Worldwide Technical Support

Telephone support for US and Canada

Call 800-227-9770

Telephone and email support for all regions

Agilent's world-wide Sales and Support Center contact details for your location can be obtained at
www.agilent.com/en/contact-us/page.

When contacting Agilent Worldwide Technical Support with a support issue, please be prepared to provide the following information:

- Instrument serial number
- Description of the issue
- Photos of the deck and barrels
- Log files
- TapeStation QC files

Any serious incident that has occurred in relation to the device shall be reported to the manufacturer and the competent authority of the country in which the user and/or the patient is established.

© Agilent Technologies, Inc. 2019, 2022, 2024

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.

The material contained in this document is provided "as is," and is subject to being changed, without notice, in future editions.

Revision D0, December 2024



K1007-90010