

BioQule NGS System Version 1.5

Software Release Notes

Contents

1.) Introduction.....	1
2.) New Features.....	1
3.) Fixes.....	2
4.) Software Compatibility.....	2
5.) System Requirements.....	2
6.) Support Information.....	2

1.) Introduction

The BioQule software offers users an intuitive interface to execute a variety of NGS workflows and assay methods on the BioQule NGS System. The software ensures researchers and laboratory professionals can effortlessly navigate the complexities of NGS procedures on the BioQule by guiding users through setting up consumables, managing multiple instrument runs, and analyzing concentration results from each library post-run.

2.) New Features

With the latest BioQule software release version 1.5 the following features have been implemented:

- BioQule folder containing all logs, results, and assay files now located in the application folder in Program Files (x86).
- Automatic inclusion of the latest scripts for all available assays upon installation of the UI, along with removal of assays folder from any previous software version.
- Added ability to copy and paste across multiple cells in sample information and results tables by right clicking.
- Added compatibility for newly released BioQule Cartridge PN. CLS158240.
- Streamlined Assay Parameters screen to only display customizable parameters for the assay selected.
- Additional customer support contact information in homepage help screen.
- Extended editable assay parameters to include tagmentation time.

- Updated workflow animations.
- Expanded allowable range of values for editable assay parameters.
- Option to replace file when adding assay with same name as existing assay.
- Automatic confirmation that results CVS file was saved upon entering the results screen.
- Added version control to assay script files.

3.) Fixes

- Improved feedback for users when editing assay parameters.
- Optimized user and script logs for better clarity.
- Prevented on screen progress bar from filling before assay is completed.
- Resolved Home assay results displaying sample concentration data from previous assay.
- Addressed cartridge barcode disappearing when going back through workflow.
- Updated warning and alert pop ups to better communicate with users.
- Ensured device name can be reset to device serial ID or previously used names.
- Corrected glitches related to use of invalid device names.
- Fixed incorrect error logged when running the home assay with BioQule Virtual Device.

4.) Software Compatibility

Software Version 1.5 is compatible with BioQule Instruments up to Revision B.

5.) System Requirements

The BioQule NGS System requires a laptop or PC with the following recommended system requirements:

- Windows 10 Professional Full Version (64-bit)
- Processor model is Intel 11th Generation Core i5 or better.
- Storage of 256 GB or better.
- RAM specifications are 4GB DDR4 3200 MHz or better.
- Screen resolution is 1920x1080.
- Windows settings updated:
 - Windows updates are turned off
 - Windows power and sleep are set to Never

6.) Support Information

Contact Revvity Technical Support by phone, fax, or e-mail from 8:00AM to 8:00PM Eastern Standard Time.

Telephone (US Toll Free): 800-762-4000

Telephone (Worldwide): +1 203-925-4602

Fax: +1 203-944-4902

Email: L3BioQule@revvity.com (L3 Support)

Internet: <https://www.revvity.com>