



# LabChip® GX/GXII Touch™ Version 1.16 LabChip® GX Reviewer Version 5.14 Software Release Notes

## Contents

Contents.....	1
1. Introduction .....	1
2. New Features .....	1
3. Changes .....	2
4. Fixes .....	2
5. Known issues.....	2
6. System Requirements .....	3
7. Installation Notes.....	3
8. Support Information .....	4

## 1. Introduction

The LabChip® GX Touch™ software controls the LabChip® GX Touch™ and LabChip® GXII Touch™ instruments. The LabChip® GX Reviewer software is used to analyze and compare data in the form of virtual gels or electropherograms across LabChip® GX/GXII Touch™ systems.

## 2. New Features

The LabChip® GX/GXII Touch™ 1.16 and LabChip® GX Reviewer Version 5.14 release features

1. New Assay Support: HT pDNA in HT pDNA chip

### 3. Changes

LabChip® GX Touch™ Version 1.16 and LabChip® GX Reviewer 5.14 have the following major changes:

1. Operation optimization for subsequent run if Chip Door not ejected (US 21898)

### 4. Fixes

LabChip® GX Touch™ Version 1.16 and LabChip® GX Reviewer 5.14 have fixed the following major issues:

1. NA

### 5. Known issues

This list describes known issues with the LabChip® GX Touch™ software version 1.16 and Reviewer version 5.14. Please read this before reporting new bugs.

Known Issues	Recommended Workaround
<b>LabChip® GX Touch™</b>	
Display setting issues for secondary monitor. Display mode may change to Extend instead of Clone.	For instruments with LabChip® GX Touch™ ver 1.7 or older: execute the command “displayswitch /clone” at startup to force the display to duplicate on the secondary monitor. For instruments with LabChip® GX Touch™ ver 1.8 or higher: in the Task Manager Startup tab, enable the “Display Switch” task to start Windows in duplicate display mode at startup. Restart instrument.
Disconnect Error during Diagnostics.	To minimize occurrences, do not run any CPU-intensive applications/process including Windows Update or Windows Defender (full scan) while LabChip® GX Touch™ is running. Power down the instrument daily.
System Diagnostics: Random communication error from the Devices leads to failure of the Diagnostic Tests.	Relaunch LabChip® GX Touch™ software and start the Diagnostics Tests again. This issue is very rare.
<b>LabChip® GX Reviewer</b>	
Windows Installer Package error displays during the GX Reviewer GxP mode installation if installed	Only occurs on certain installation order. Install in the following order to avoid this issue:

immediately after uninstalling GX Touch from the control panel.	<ol style="list-style-type: none"> <li>1. Uninstall LabChip® GX Reviewer from the control panel</li> <li>2. Uninstall LabChip® GX Touch™ from the control panel</li> <li>3. Install LabChip® GX Touch™ with GxP mode on the system</li> </ol>
Print Dialog: Footer settings are not retained after relaunch the Print dialog.	User needs to enable or disable the footer settings manually.
Version History: Unwanted details are populated in Version History for Protein Express data files when user enable\disable exclude system peaks option.	None.
Connection failed error when uninstalling Reviewer from a PC in which GX Touch application mapped with remote server.	Restart the DatafileService, or restart the instrument.
Version History inappropriately logs Bracket Ladder Alignment changes for Glycan Extended Rage data file when user makes any Analysis Settings changes.	None.

## 6. System Requirements

LabChip® GX Touch™ requires a LabChip® GX/GXII Touch instrument.

LabChip® GX Reviewer software can be used on a personal computer with following Microsoft Windows Operating Systems:

- Windows® 10 Professional Full Version (64-bit)
- Windows® 10 Enterprise 2019 LTSC (64-bit)
- Windows 11 IoT Enterprise LTSC 24H2

Minimum RAM: 1GB (2GB recommended).

Minimum free hard drive space: 2 GB.

## 7. Installation Notes

The LabChip® GX Touch™ software 1.16 is pre-installed on new instruments manufactured after Jan 2025. The LabChip® GX Reviewer software can be installed on additional computers in reviewer mode to analyze data files.

Customers with Windows 10 on the instrument can upgrade from LabChip® GX Touch™ software version 1.9 and higher and from LabChip® GX Reviewer software version 5.7 and higher.

To upgrade non-GxP software in the instrument to GxP version, only the Reviewer need to be reinstalled with the GxP option. GxP mode in the Touch application is enabled by the USB dongle.

Both Touch and Reviewer software must be upgraded before using the software; either upgrading software revisions or upgrading to GxP mode software.

The LabChip® GX Touch – User shall login to the Instrument PC using credentials GXTUser\GXTUser

## **8. 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: +1 203-925-4602

Email: L3LabChip@Revvity.com

Internet Address: <http://www.Revvity.com>