

GeneChip™ Command Console™ (GCC)

Software v7.0.3

Pub. No. MAN1002339 Rev. A

Note: The following release notes provide information about GeneChip™ Command Console™ (GCC) software v7.0.3.

Included in GCC v7.0.3

- GCC v7.0.3 is a patch installer. It requires GCC 7.0.1 or GCC 7.0.2 to already be installed before updating to GCC 7.0.3.
- GCC v7.0.3 includes an improvement for Automated Axiom™ Analysis (AutoAx) customers. The GCC uploader tool which is now incorporated into the GCC Launcher is now named “Automated Axiom Configurator”.

Set up the Automated Axiom Configurator

GCC 7.0.3 is only supported on the US English version of Windows™ 10 Enterprise 2019 LTSC operating system.

1. From the GCC Launcher, doubleclick **Automated Axiom Configurator**.
The Automated Axiom Configurator window launches and is displayed.
2. Enter the IP address or port number for the AutoAx Server.
3. Click **Enable Auto Data Transfer** to start the GCC File Copier Service.
The GCC File Copier Service is started and runs with a startup type of **Automatic** in the services control panel.
The system is ready to use.

Known issues in GCC 7.0.1, GCC 7.0.2, and GCC 7.0.3

- When adding a data root to a second workstation, any projects in that root are not added to the second location. In this case, go to **Project Management** and set the project path as primary. This updates the system and allows samples to be registered.
- GeneTitan™ Instrument intermittently fails to start after a shutdown. If this occurs, restart the instrument and workstation.
- The GeneTitan™ user interface displays the full status of the next available workflows. The Thermo Fisher Cloud user interface only displays the earliest start time for various possible workflows. If you need to view these items, use the GeneTitan™ user interface.
- The recommended procedure “to add or remove a user to monitor the GeneTitan™ MC Instrument status” is to use the Thermo Fisher Cloud Administration Utility in GCC.
- In the **Details Page**, the **Error Log** Window displays the UTC time (displayed in bold) and the local time when the error occurred. To identify the time of failure, use the local instrument time displayed that is in the error message.
- The Thermo Fisher Cloud user interface displays the completion time for the current run in UTC time instead of local time. Take this difference into account when scheduling subsequent runs.
- For the instrument remote monitoring, do not use the Internet Explorer™ browser as it can sometimes not update the status properly. Use either the Chrome™ or Firefox™ browser for remote monitoring.
- Enabling and disabling cloud connectivity is currently permitted when the instrument is active. It is recommended that customers do not change this setting with the instrument control application running.
- In the v7.0.3 release, when hardware errors occur, the users can see a loss of throughput, due to automated restart/relaunch of the Scanner software.
- In the v7.0.3 release, the **Reconnector** application's Help function does not work. The alternative for this problem is to manually open the **Help** document and navigate to the section on this application.
- In the v7.0.3 release, opening a corrupt/incomplete DAT file can cause an abnormal exit of the **Image Viewer** application.

Limited product warranty

Life Technologies Corporation and its affiliates warrant their products as set forth in the Life Technologies' General Terms and Conditions of Sale at www.thermofisher.com/us/en/home/global/terms-and-conditions.html. If you have questions, contact Life Technologies at www.thermofisher.com/support.

 Affymetrix, Inc. | 6055 Sunol Blvd | Pleasanton, California 94566 USA
For descriptions of symbols on product labels or product documents, go to thermofisher.com/symbols-definition.

Revision history: Pub. No. MAN1002339 A

Revision	Date	Description
A	22 June 2026	Initial release of document. Software version 7.0.3. GCC File Uploader has been incorporated into the tool and renamed Automated Axiom Configurator.

The information in this guide is subject to change without notice.

DISCLAIMER: TO THE EXTENT ALLOWED BY LAW, THERMO FISHER SCIENTIFIC INC. AND/OR ITS AFFILIATE(S) WILL NOT BE LIABLE FOR SPECIAL, INCIDENTAL, INDIRECT, PUNITIVE, MULTIPLE, OR CONSEQUENTIAL DAMAGES IN CONNECTION WITH OR ARISING FROM THIS DOCUMENT, INCLUDING YOUR USE OF THE DOCUMENT OR THE PRODUCT.

Important Licensing Information: These products may be covered by one or more Limited Use Label Licenses. By use of these products, you accept the terms and conditions of all applicable Limited Use Label Licenses.

©2026 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified. Chrome is a trademark of Google LLC. Firefox is a trademark of the Mozilla Foundation in the U.S. and other countries. Internet Explorer and Windows are registered trademarks of Microsoft Corporation in the United States and/or other countries.