

SeqStudio™ Genetic Analyzer

Pub. No. MAN0019362 Rev. A.0

This document includes a checklist to help you prepare your SeqStudio™ Genetic Analyzer for use after an extended shut down. We encourage you to complete the tasks that apply to your situation, then retain the completed checklist with the instrument service records.

Also, we recommend that you review the maintenance procedures in the *SeqStudio™ Genetic Analyzer Instrument and Software User Guide*, and ensure that you perform these tasks when and as instructed.

Resources

I need ...	Resource
Step-by-step instructions for tasks in this checklist	<i>SeqStudio™ Genetic Analyzer Instrument and Software User Guide</i> (Pub. No. MAN0016138)
A list of instrument consumables and reagents	See "Catalog numbers" on page 2.
To access more SeqStudio™ Genetic Analyzer tools, documentation, or troubleshooting content	Visit the Applied Biosystems™ SeqStudio™ Genetic Analyzer Support Center .
General product support	Visit thermofisher.com/support .
To request technical support or a service call	Visit thermofisher.com/contactus .

Instrument serial number:

Instrument location:

Instrument owner:

Restart the SeqStudio™ Genetic Analyzer after an extended shutdown

✓	Task
If the instrument has been powered off for more than 2 weeks, then perform the following tasks:	
<input type="checkbox"/>	Restart the SeqStudio™ Genetic Analyzer.
<input type="checkbox"/>	Eject the plate. Eject the autosampler tray, then: <ul style="list-style-type: none"> • Wipe down the autosampler tray with a lint-free wipe moistened with deionized water. • Install new cathode buffer container.
<input type="checkbox"/>	If the SeqStudio™ Cartridge is in the instrument, then eject it. Install a new SeqStudio™ Cartridge or a previously used SeqStudio™ Cartridge that has been properly stored. See the <i>SeqStudio™ Genetic Analyzer Instrument and Software User Guide</i> for proper cartridge storage instructions.
If the instrument has been powered on and idle with consumables installed for more than 2 weeks, then perform the following tasks:	
<input type="checkbox"/>	Eject the plate.
<input type="checkbox"/>	Remove the cathode buffer container, and note if the buffer level has fallen below the fill line.
<input type="checkbox"/>	Wipe down the autosampler tray with a lint-free wipe moistened with deionized water.
<input type="checkbox"/>	Replace the cathode buffer.
<input type="checkbox"/>	Eject the SeqStudio™ Cartridge.
<input type="checkbox"/>	Inspect the SeqStudio™ Cartridge at the capillary array, polymer delivery system, and universal polymer supply. <ul style="list-style-type: none"> • If there are white crystals, then replace the SeqStudio™ Cartridge with a new cartridge or a previously used SeqStudio™ Cartridge that has been properly stored. • If the cathode buffer level was below the fill line, then replace the SeqStudio™ Cartridge with a new cartridge or a previously used SeqStudio™ Cartridge that has been properly stored. • If neither of above, then reinstall the SeqStudio™ Cartridge.
<input type="checkbox"/>	If you did not replace the SeqStudio™ Cartridge, then touch Settings ▶ Cartridge ▶ Cartridge maintenance ▶ Refresh PDS .
<input type="checkbox"/>	Optional. Follow the instructions in the <i>SeqStudio™ Genetic Analyzer Instrument and Software User Guide</i> to perform an install run using sequencing or fragment analysis installation standards.

Catalog numbers

Part description	Cat. No.
SeqStudio™ Genetic Analyzer Cartridge v2	A41331
Cathode Buffer Container	A33401
Reservoir Septa	A35640
BigDye™ Terminator 3.1 Sequencing Standards	4404312
DS-33 GeneScan™ Installation Standards with GeneScan™ 600 LIZ™ Size Standard v2.0	4376911
Hi-Di™ Formamide (25 mL)	4311320



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 IT.

Revision history: Pub. No. MAN0019362

Revision	Date	Description
A.0	21 May 2020	New document.

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.

©2020 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified.