

Ion Chef™ Instrument v4.4 Software

Pub. No. MAN0013498 Rev. A.0

This document highlights the main changes in the Ion Chef™ Instrument software from v4.2.

IMPORTANT! When updating Torrent Suite™ Software (TSS), it is critical that users update both their instrument software and their Torrent Server. For TSS 4.0 or higher, this is especially important because the instrument will not retrieve the proper on-instrument analysis if the instrument is running an older version of software on the Torrent Server.

IMPORTANT! If your Ion Chef™ Instrument is currently configured to see more than one Torrent Server, you will need to follow the guidelines in the *Ion Chef™ and Torrent Server Network Setup User Guide* (Pub. no. MAN0013444) to reconfigure that connection. This will require choosing a primary Torrent Server as your direct Ion Chef™ connection. Then, from that primary server, you will need to enable Planned Run sharing in the Torrent Browser, which will allow you to see additional Torrent Servers.

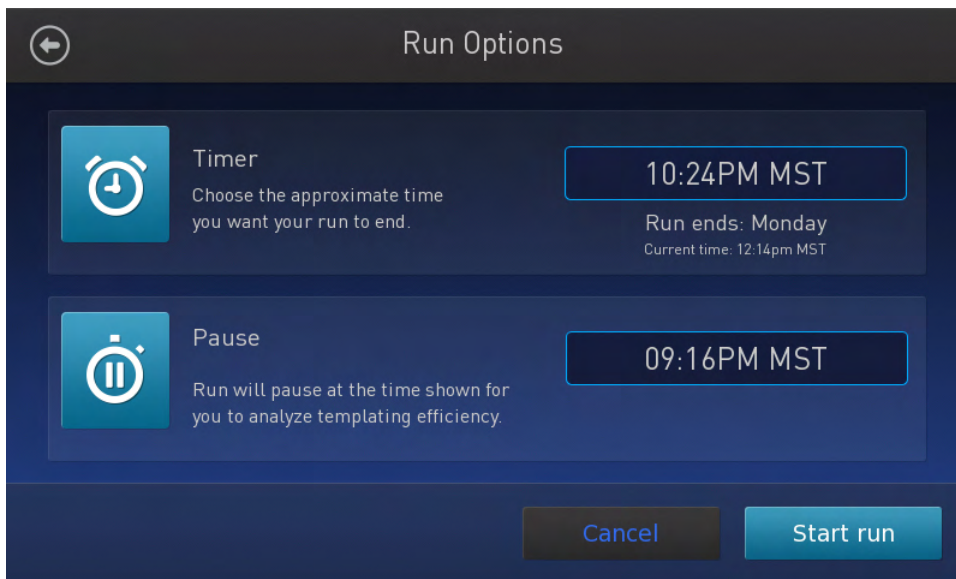
- User interface improvements 1
- Planned Run sharing 2
- Ion Chef™ System and Torrent Server connectivity 2
- Ion PGM™ Hi-Q™ Chef Kit enabled 2
- Deck Scan improvements (previously Load Check) 3
- Other improvements and bug fixes 3
- Limited product warranty 3

User interface improvements

- More intuitive user interface.



- Run Scheduler now allows selection of the desired completion time, rather than hours to completion.
- Run Progress now estimates the predicted pause time when the **Pause** option is selected for QC sample collection.



Planned Run sharing

The new Planned Run sharing function allows users to move Planned Runs from one Torrent Server to another Torrent Server. This functionality is ideal for creating chips on an Ion Chef™ Instrument and running them on multiple Ion PGM™ or Ion Proton™ Sequencing instruments.

Note: To enable this functionality, please refer to the Torrent Suite™ Software User Guide documentation and/or the *Ion Chef™ and Torrent Server Network Setup User Guide* (Pub. no. MAN0013444). The guides provide step-by-step details on how to configure your network and transfer Planned Runs between Torrent Servers.

Ion Chef™ System and Torrent Server connectivity

- An Ion Chef™ Instrument now supports only one Torrent Server connection. Confirm that your primary Torrent Server connection is correct and reconfigure if necessary. Please refer to the *Ion Chef™ and Torrent Server Network Setup User Guide* (Pub. no. MAN0013444) for detailed instructions to configure this.
- The specified Torrent Server will be the default FTP server for log transfer. All Ion Chef™ logs will stay on the designated primary server.
- Ion Chef™ status can now be monitored at the Torrent Server Monitor Page:

Ion Chef Refresh

| Date | Plan | Chef Instrument | Stage Progress | Stage Status |
|--------------------------|----------------------------------------|-----------------|---------------------------------------------------------------------------|---------------------------------------|
| Nov. 18, 2014, 1:52 p.m. | chef_monitor_IC_200_template_kit_plan | chef-alpha | <div style="width: 100%; height: 10px; background-color: #0070C0;"></div> | Starting UnlockDoor |
| Nov. 18, 2014, 4:12 p.m. | chef_monitor_IC_200_template_kit_plan2 | chef-beta-2 | <div style="width: 100%; height: 10px; background-color: #0070C0;"></div> | Starting UnlockDoor |
| Nov. 18, 2014, 7:20 p.m. | chef_monitor_IC_200_template_kit_plan3 | chef-delta-3 | <div style="width: 25%; height: 10px; background-color: #0070C0;"></div> | Enriching for Template-Positive Beads |

Ion PGM™ Hi-Q™ Chef Kit enabled

For Ion PGM™ Hi-Q™ Chef Kit users only:

- Selection of 200-bp or 400-bp template size is enabled. Ion Chef™ runs must be set up with two libraries of the same template size. Refer to the *Ion PGM™ Hi-Q™ Chef Kit User Guide* (Pub. no. MAN0010919) for details.
- Single-chip workflow is now supported. Refer to the user guide for details.

Deck Scan improvements (previously Load Check)

- Issues are detected more quickly, reducing the average instrument setup time.
- Deck Scan now spins the centrifuges to detect potential sources of run failure and common setup issues.
- Deck Scan now has improved warning messages to guide correction of deck setup issues.
- A post-run Deck Scan is now available to confirm that all consumables are removed before UV treatment.

Other improvements and bug fixes

- Better log cleanup to archive and remove old log files.
- More log files transferred to the Torrent Server to aid with troubleshooting.

Limited product warranty

Life Technologies Corporation and/or its affiliate(s) warrant their products as set forth in the Life Technologies' General Terms and Conditions of Sale found on Life Technologies' website at www.lifetechnologies.com/termsandconditions. If you have any questions, please contact Life Technologies at www.lifetechnologies.com/support.

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

DISCLAIMER

LIFE TECHNOLOGIES CORPORATION AND/OR ITS AFFILIATE(S) DISCLAIM ALL WARRANTIES WITH RESPECT TO THIS DOCUMENT, EXPRESSED OR IMPLIED, INCLUDING BUT NOT LIMITED TO THOSE OF MERCHANTABILITY, FITNESS FOR A PARTICULAR PURPOSE, OR NON-INFRINGEMENT. TO THE EXTENT ALLOWED BY LAW, IN NO EVENT SHALL LIFE TECHNOLOGIES AND/OR ITS AFFILIATE(S) BE LIABLE, WHETHER IN CONTRACT, TORT, WARRANTY, OR UNDER ANY STATUTE OR ON ANY OTHER BASIS FOR SPECIAL, INCIDENTAL, INDIRECT, PUNITIVE, MULTIPLE OR CONSEQUENTIAL DAMAGES IN CONNECTION WITH OR ARISING FROM THIS DOCUMENT, INCLUDING BUT NOT LIMITED TO THE USE THEREOF.

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

©2015 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified.

For support visit lifetechnologies.com/support or email techsupport@lifetech.com

lifetechnologies.com