

iBlot 2 Gel Transfer Device Firmware Release Notes

This guide covers upgrading the Invitrogen iBlot™ 2 Gel Transfer Device to firmware version 1.5.1. This update is recommended for any iBlot 2 containing firmware version v1.4 or earlier. Update not required for system containing firmware version 1.5.0. The current firmware version installed on a iBlot 2 is displayed on the **Firmware Version** button within the **Options** screen. Download a copy of the latest firmware by visiting www.thermofisher.com/iblotfirmwareupdate.

Firmware version v1.5.1 does not add any new features; however, v1.5.1 is a required update as it eliminates a bug within the memory and file management systems that can result in a system failure during updates or when using custom transfer programs. Installation of firmware v1.5.1 can also be used to directly recover a system that has become frozen due to this specific file maintenance bug.

Product #IB21001
Revision Date: October 17, 2018
Firmware version 1.5.1

Upgrading iBlot 2 Gel Transfer Device Firmware

Note: Before upgrading your firmware, it is recommended that parameters for custom methods be recorded in a lab notebook, and any run logs that you want to keep be transferred to a USB storage device as these files will not be maintained during the update process.

Standard Update Process or recovering a unit with a frozen screen:

1. Visit www.thermofisher.com/iblotfirmwareupdate and follow the instructions to download the latest firmware version to a USB storage device.

Note: The firmware file should be saved in the main directory of the USB storage device. Please delete any previous firmware (.ib2 files) on the USB storage device.

2. Power off the iBlot2 Gel Transfer Device.
3. Insert the USB storage device containing the new firmware version (".ib2" file) into the USB port (Type A) of the iBlot 2 Gel Transfer Device.
4. Turn on the iBlot2 Gel Transfer Device.
5. Installation will start automatically. Do not remove the USB storage device until the unit beeps and the **Date & Time** screen is displayed. Total time should be approximately 90 seconds.

Note: The download is complete after **Date & Time** is displayed, after the unit beeps it will take approximately another 15-20 seconds for the software to complete the setup process before the **Date & Time** screen appears. USB should remain in the device during this process

6. After the **Date & Time** screen appears, turn off the device.
7. Remove the USB.
8. Power on the device and complete set-up.
9. After set-up is complete, go to the Options screen to verify that the new firmware version is displayed.

If you experience any issues with the iBlot 2 update process or if you have additional questions, please contact technical support (technicalsupport@thermofisher.com).

Sincerely,



Paul Haney Ph.D.
Sr. Product Manager
Thermo Scientific Pierce Protein Research
3747 N Meridian Road
Rockford, IL 611