


BigDye XTerminator™ Purification Kit

BDX Utility and run modules for 3730/3730x/ and 3130/3130x/ instruments with Data Collection v4 Software or earlier

Catalog Numbers 4376484, 4376485, 4376486, and 4376487

Pub. No. 4483510 Rev. C

 **WARNING!** Read the Safety Data Sheets (SDSs) and follow the handling instructions. Wear appropriate protective eyewear, clothing, and gloves. Safety Data Sheets (SDSs) are available from thermofisher.com/support.

This user bulletin provides:

- A list of BigDye XTerminator™ (BDx) run modules
- Instructions for downloading and installing the BDX Utility
- Instructions for managing BDx run modules

Overview

BDx run modules adjust the injection height of the instrument to allow direct injection of the sample supernatant following purification with the BigDye XTerminator™ Purification Kit. In addition, BDx run module injection times and voltages are optimized for the high signals obtained from samples purified with the kit.

The BDX Utility installs and calibrates BDx run modules for the Data Collection Software of the instrument. Run the BDX Utility after Data Collection Software installation and every time the instrument autosampler is recalibrated.

Note: The BDX Utility is included in the Data Collection v5 Software installation. No separate download or installation is necessary.

IMPORTANT! Do not use a BDx run module if the supernatant is transferred to a new plate after purification. See the instrument user guide for the appropriate standard run module.

BDx run modules

Table 1 BDx run modules for the 3730/3730x/ instrument

Capillary length	Polymer	BDx run module name
36-cm	POP-7™	BDx_StdSeq36_POP7
		BDx_RapidSeq36_POP7
50-cm	POP-7™	BDx_XLRSeq50_POP7
		BDx_LongSeq50_POP7
		BDx_FastSeq50_POP7

Table 2 BDx run modules for the 3130/3130x/ instrument

Capillary length	Polymer	BDx run module name
36-cm	POP-7™	BDx_RapidSeq36_POP7
	POP-6™	BDx_RapidSeq36_POP6
	POP-7™	BDx_UltraSeq36_POP7
	POP-4™	BDx_UltraSeq36_POP4
50-cm	POP-7™	BDx_StdSeq50_POP7

Capillary length	Polymer	BDx run module name
	POP-6™	BDx_StdSeq50_POP6
	POP-4™	BDx_StdSeq50_POP4
	POP-7™	BDx_FastSeq50_POP7
80-cm	POP-7™	BDx_LongSeq80_POP7
	POP-4™	BDx_LongSeq80_POP4

Install BDx run modules

Download, then install the BDx Utility

1. Power on the instrument and the computer with your Data Collection Software.
2. Exit the Data Collection Software if it is running.
3. Go to thermofisher.com/us/en/home/technical-resources/software-downloads/bigdye-xterminator-purification-kit.html.
4. Click **Download**.
5. Double-click the **bdx1.0.1updater.exe** file.
6. Follow the prompts in the **Support Files Installer** wizard.
 - Data Collection v3 Software—The run modules are installed and calibrated based on the instrument’s autosampler calibration.
 - Data Collection v3.1, v3.1.x, or v4 Software on 3730/3730xl instruments—The run modules are installed and calibrated based on the instrument’s autosampler calibration.
 - Data Collection v3.1, v3.1.x, or v4 Software on 3130/3130xl instruments—Proceed to “Install, then calibrate BDx run modules (for 3130/3130xl instruments only)” on page 2.

Install, then calibrate BDx run modules (for 3130/3130xl instruments only)

Follow this procedure if you are using the 3130/3130xl instrument with Data Collection v3.1, v3.1.x, or v4 Software.

In this procedure, the BDx Utility is uninstalled, then reinstalled to update BDx run modules.

IMPORTANT! Uninstalling the BDx Utility removes all existing BDx objects from the Data Collection Software database.

1. Power on the instrument and the computer with your Data Collection Software.
2. Exit the Data Collection Software if it is running.
3. Select **Start ▶ Programs ▶ Applied Biosystems™ ▶ BDx Utility ▶ Uninstall BDx**.
4. Follow the prompts in the **Uninstaller** wizard.
All existing BDx run modules are removed.
5. Double-click the **bdx1.0.1updater.exe** file to run the BDx Utility.
6. Follow the prompts in the **Support Files Installer** wizard.
The BDx Utility installs new BDx run modules and calibrates the run modules based on the instrument’s autosampler calibration.

Update BDx run modules

Run the BDx Utility to update BDx run modules after Data Collection Software installation and every time the instrument autosampler is recalibrated.

1. Exit the Data Collection Software if it is running. Do not shut down the data-collection computer.
2. Power on the instrument.
3. Select **Start** ▶ **Programs** ▶ **Applied Biosystems™** ▶ **BDx Utility** ▶ **Update BDx Utility**.

Uninstall BDx run modules

IMPORTANT! Uninstalling a BDx run module also removes any instrument protocols and plate records that use the run module. Run folders, containing AB1 files that have been extracted to the hard drive, are not removed.

1. Power on the instrument and the computer with the Data Collection Software.
2. Exit the Data Collection Software if it is running.
3. Select **Start** ▶ **Programs** ▶ **Applied Biosystems™** ▶ **BDx Utility** ▶ **Uninstall BDx**.
4. Follow the prompts in the **Uninstaller** wizard.



Life Technologies Corporation | 2 Preston Court | Bedford, MA 01730

For descriptions of symbols on product labels or product documents, go to [thermofisher.com/symbols-definition](https://www.thermofisher.com/symbols-definition).

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. 4483510

Revision	Date	Description
C	2 June 2020	<ul style="list-style-type: none">• Updated compatible instruments.• Updated list of BDx run modules.• Updated instructions for using the BDx Utility.• Made minor wording and formatting updates.
B	12 November 2018	Rebranded. Updated download instructions and information on Data Collection Software v5.0.
A	February 2013	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.