

User Bulletin

3730/3730xl DNA Analyzer Data Collection Software

May 2009

SUBJECT: Data Collection Software version 3.1

In this user bulletin

This user bulletin covers:

■ New features	1
■ Compatibility matrix	2
■ System requirements	2
■ Important operational notes	3

New features

- **Windows® Vista Support** - 3730/3730xl Data Collection v3.1 Software now supports running on the Windows® Vista (32 bit, Business Edition) operating system.
- **Important Results - group capacity warning.** When the 3730/3730xl Data Collection v3.1 software assigns the Results Group for Autoanalysis with Secondary applications (GeneMapper), the software now displays a warning prompt when the number of samples assigned to the result group reaches 80% of the maximum limit (maximum = 7,000 samples).
- **Support for the POP6™ Polymer** - 3730/3730xl Data Collection v3.1 Software supports the POP6™ Polymer.
- **New run modules** - Several new run modules and dye sets are included in this version:
 - BDX Run Module
 - TargetSeq Run Module
 - SNPlex Dye set and Run Module
 - GS 1200 Run Module
- **Updated wizards** support the changing of polymers.
- **Database upgrade** - The embedded database has been upgraded from Oracle 8i to Oracle 11g. This provides improved performance, and enhances database security with transparent data encryption.
- **New license scheme** - Data Collection v3.1 software now uses a new license scheme requiring registration on-line.

Note: This User Bulletin is for users of Data Collection software v3.1 and **not** for users of Data Collection software v3.0 who use GeneMapper ID v3.2 or v3.2.1, Genemapper IDX v1.0, v1.0.1 or v1.1 or MicroSeq ID v2.0.

Compatibility matrix

This release supports the Applied Biosystems® 3130/3130xl instruments. You can view and analyze sample files generated by this software with the following software products:

- Sequencing Analysis v5.4
- SeqScape® v2.7
- GeneMapper® v4.1
- GeneMapper® IDX v1.2
- Variant Reporter™ v1.1
- MicroSeq ID® 2.2

System requirements

The minimum supported computer hardware and system requirements for 3730/3730xl Data Collection v3.1 are as follows:

- **Minimum hardware requirements** - Intel® Pentium® IV computer with 1GB RAM, 3.0 GHz processor, 120 GB of free hard disk space.
Display - Recommend 17+ inches, resolution 1024 x 768 and TCP/IP protocol configured.
- Recommended hardware requirements -Intel® Dual Core 755 OptiPlex computer with 2 GB RAM, Intel® Core™ 2 Duo CPU E8400 3.0 GHz processor, 320 GB of free hard disk space.
Display - Recommend 17+ inches, resolution 1024 x 768 and TCP/IP protocol configured.
- The software has been tested on a Dell OptiPlex 755 Workstation with Intel® Core™ 2 Duo CPU E8400(3.0 GHz) with 2 GB RAM.
- 17-inch Color Monitor (1024 x 768 resolution or higher)
- **Operating system options:**
 - Windows® Vista Business Edition Operating System with Service Pack 1 or
 - Windows® XP Professional operating system with Service Pack 2 or
 - Windows® XP Professional operating system with Service Pack 3

Note: Although you can install 3730/3730xl Data Collection v3.1 on a computer with less than 1 GB of RAM and 120GB hard disk space, Applied Biosystems highly recommends using 2 GB of RAM (or higher) and 320 GB of free hard disk space (or higher) for better performance. 3730/3730xl Data Collection v3.1 has not been tested on a computer with server-type hardware or server operating systems.

Important operational notes

- Applied Biosystems recommends that you run a spectral calibration each time that a new capillary array is installed. In 3730/3730xl Data Collection v3.1 software, if you install an array that is the same length as the previously installed array, the active spectral calibration still persists. For optimal data quality, perform a new spectral calibration before you perform regular runs.
- Applied Biosystems has successfully used Norton Anti-Virus™ Corporate software in our Data Collection Software testing environment with no problems.
- 3730/3730xl DNA Analyzer Data Collection v3.1 Software does not support the switch-user mode in Windows® XP and Vista operating systems. Applied Biosystems recommends that you use the 3730User user to run this version of software, or log in as Administrator.
- Ensure that the “Automatically adjust clock for daylight saving changes” checkbox is checked. If it is not, timestamps of data run folders might be wrong by an hour.
- The C: partition on the disk drive should not be used for data storage. This partition is only 8 GB in size, and most of that space is taken by the storage and running of the operating system. If the C: partition is used as a location to store run data, the software does not do any disk space checks to ensure that there is enough space to store the run. This means that you might lose data!
- The Results Group Destination tab does not recognize remote storage locations. Map the network location to a local drive letter by using the Map Network Drive feature of the operating system. Then, specify the mapped drive letter location in the Results Group Destination tab.
- In the ESig Administration, Configuration tab, the Require User ID checkbox should be checked if electronic signature is required. Otherwise, a default username might appear when ESig is activated.
- Ensure that all Data Collection Services are running before you launch the AB Navigator tool for security, audit trail and electronic signature features. You know that all services are running when the Service Console contains four green squares.
- Do not change the computer's date or time during a run. If you change the system time more than 20 times the average run time during a run, some runs will not be analyzed.
- Copying and pasting within the plate editor -The Plate Editor Copy and Paste functionality is supported only within one plate editor. Users are not allowed to copy and paste the contents of one plate to another. Use the Duplicate Plate feature instead.

If multiple cells are selected for copying, select the same number of corresponding target cells before you execute the Paste command.
- Use the manual control of the autosampler feature with care. Give attention to the autosampler position when you use Manual Control to move the autosampler. If the autosampler reaches its maximum travel position, a fatal instrument error is generated. The software needs to be re-started after this error.
- To re-use a plate after deleting the analysis protocol associated with it, either create another plate, or re-create the analysis protocol with the same name.

- When you use the Service Console, do not press Start All or Stop All when the Data Service displays the yellow triangle. If this happens, you need to re-boot the computer.
- Data Collection software might not be able to recover properly if the free disk space on the E: drive is less than 2 GB. Use of auto-analysis increases this chance. To prevent this error, perform regular file system maintenance. Move files from the E:\ drive to archival locations regularly.
- When you add another sample run to a processed plate, all the information in the processed runs should be valid. Otherwise, that data will not be validated and a new sample run cannot be created.
- If you use the duplicate plate function, all the information in the plate to be duplicated must be valid. Otherwise, an empty plate is created.
- When you use an array of a different length, all spectral calibrations are voided. You need to go to the Spectral Viewer and choose a past calibration to set an active spectral calibration. You cannot link or run a plate unless the dye set used in the plate has been set in the Spectral Viewer. Furthermore, when a plate is running, the Set Active Spectral Calibration function is inactive. Spectral calibrations can be set only during the idle or ready mode.
- In Applied Biosystems 3730/3730xl DNA Analyzer Data Collection v2.0 Software, if no active spectral calibration is set, the spectral viewer does not display a spectral profile. However in 3730/3730xl DNA Analyzer Data Collection v3.1 software, there is a case when spectral profiles are displayed even when no active spectral calibration is set: When you install a different-length array associated with historical spectral calibrations. Be sure to set an active spectral calibration before you proceed with regular runs.
- In the 3730/3730xl DNA Analyzer Data Collection v2.0 Software, in the Re-Extraction panel, color codes that indicated ranges of values were assigned to the Score column of sequencing plates. In 3730/3730xl DNA Analyzer Data Collection v3.1 software, the numeric quality scores are written directly.
- Manual changes to the instrument calibration text file are not automatically sent to the instrument during an instrument reboot, because of a change in the software and firmware design.
- Users might observe the error message listed under the Data Service console, "The requested service has already been started. More help is available by typing NET HELPMSG 2182." This message is an informative message and should be ignored.
- Changes in the wizards found in the 3730/3730xl Data Collection v3.0.6 and v3.1 result in longer wizard setup time for the Change Polymer Type wizard and the Instrument Shutdown wizard. Both wizards take approximately 20 extra minutes to complete.

For more information about any of these topics, please refer to the 3730/3730xl Data Collection Software v3.1 Release Notes.

© Copyright 2009, Applied Biosystems. All rights reserved.

Information in this document is subject to change without notice. Applied Biosystems assumes no responsibility for any errors that may appear in this document.

APPLIED BIOSYSTEMS DISCLAIMS ALL WARRANTIES WITH RESPECT TO THIS DOCUMENT, EXPRESSED OR IMPLIED, INCLUDING BUT NOT LIMITED TO THOSE OF MERCHANTABILITY OR FITNESS FOR A PARTICULAR PURPOSE. IN NO EVENT SHALL APPLIED BIOSYSTEMS 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.

TRADEMARKS:

AB(Design) and GeneMapper are trademarks of Applied Biosystems or its subsidiaries in the U.S. and/or certain other countries. ABI PRISM and SeqScape are registered trademarks of Applied Biosystems or its subsidiaries in the U.S. and/or certain other countries.

Microsoft and Windows are registered trademarks of Microsoft Corporation.

All other trademarks are the sole property of their respective owners.

Worldwide Sales and Support

Applied Biosystems vast distribution and service network, composed of highly trained support and applications personnel, reaches 150 countries on six continents. For sales office locations and technical support, please call our local office or refer to our Web site at www.appliedbiosystems.com.

Applied Biosystems is committed to providing the world's leading technology and information for life scientists.

Headquarters

850 Lincoln Centre Drive
Foster City, CA 94404 USA
Phone: +1 650.638.5800
Toll Free (In North America): +1 800.345.5224
Fax: +1 650.638.5884

Part Number 4440403 Rev A

05/2009