

## Files for High Throughput SNP and Microsatellite Analysis with the 22cm Capillary Array for the ABI PRISM® 3100 Genetic Analyzer

November 11, 2001

The installer CD contains all the necessary files for performing high throughput SNP or microsatellite analysis with the 22cm capillary array. In order to install the necessary files the system must be running the following versions of Data Collection and GeneScan® Analysis Software:

1. Data Collection Software version 1.0.1
2. GeneScan® Analysis Software version 3.7.

What does the installer contain?

- 1) Contains all files needed to collect and analyze data generated with ABI PRISM® SNaPshot® Multiplex Kit, Linkage Mapping Set version 2.5 or custom oligos.
- 2) Contains new module files for running both applications with ABI PRISM® 3100 POP-4(TM) Polymer.
- 3) Contains a Spectral Calibration module to generate a matrix calibration with the 22 cm array.

-----

For installation, what is required before starting?

- 1) Before running the updater, reboot the computer. After the computer has finished restarting, close all open windows and shut down running applications.

Following installation:

- 2) Restart the ABI PRISM® 3100 instrument. (You will be prompted to do this at the end of installation procedure.)

-----

Installation Instructions:

Insert the installation CD and follow the directions provided in the installation procedure that appears.

Note: If you un-install the Data Collection Software V1.0.1, you will need to re-install the software support files using this installer CD.

- 1) Check to see all files have been successfully installed in the following locations:

GeneScan22\_POP4.mtd D:\AppliedBio\Support Files\Data Collection Support Files\Method Files

Spect22\_POP4.mtd D:\AppliedBio\Support Files\Data Collection Support Files\Method Files

SNP22\_POP4.mtd      D:\AppliedBio\Support Files\Data Collection Support  
Files\Method Files

For more product information, please visit [www.appliedbiosystems.com/ga](http://www.appliedbiosystems.com/ga)

ABI PRISM, GeneScan, and SNaPshot are registered trademarks and POP-4 is a  
trademark of Applied  
Corporation or its subsidiaries in the US and certain other countries.