

Importing the SNaPshot™ Multiplex System Running Module for ABI PRISM® 3700 DNA Analyzer ONLY

The SNP1_1POP5 run module must be imported into the instrument database. Once it is imported, a default SNP1_1POP5 run module is created in the 3700 Data Collection Software.

To import the SNP1_1POP5 run module:

Step	Data Collection	Action
1	v 1.1	Copy "SNP1_1POP5.mtd" method file into the following directory: D: Perkin-Elmer\Abi\Support Files\Data Collection Support Files\Method Files
1	v 1.1.1	Copy "SNP1_1POP5.mtd" method file into the following directory: D: AppliedBio\Abi\Support Files\Data Collection Support Files\Method Files
2	v 1.1	Using Windows NT Explorer, navigate to the method importer utility called NewMethodImportUtility.bat in the following directory: D:\Perkin-Elmer\Abi\3700\Bin
2	v 1.1.1	Using Windows NT Explorer, navigate to the method importer utility called NewMethodImportUtility.bat in the following directory: D:\AppliedBio\Abi\3700\Bin
3	Both	Right-click NewMethodImportUtility.bat
4		Click Edit from the pop-up menu. This opens the utility's batch file in the Notepad program.
5		Look on the third line of the text for a string representing a file path to the method file. Within the string, you will see the text: INSERTNAME.mtd
6		Replace the text "INSERTNAME" with "SNP1_1POP5". IMPORTANT Do not replace or delete the .mtd file extension.
7		Select Save As from the File menu. This opens the Save As dialog box.
8		In the Save As dialog box, type "SNP1_1POP5". Click Save .
9		Select Exit from the File menu. This closes the Notepad program.
10		Double-click the file you just created to run the utility. The program window appears for approximately one minute and then will ask you to press any key to continue.
11		The run method should now appear in the Data Collection module editor as: "SNP1_1POP5DefaultModule".