

ReadMe\_IDDirect\_v1x.txt

=====  
About GeneMapper® ID-X Identifier Direct™ Software v1X  
Release Notes  
Copyright 1999-2009, Applied Biosystems. All rights reserved.  
October, 2009  
=====

TABLE OF CONTENTS:

- \* Overview
- \* Software Contents
- \* Validation Summary
- \* File Install Instructions
- \* Legal Disclaimer

-----  
OVERVIEW  
-----

GeneMapper® ID-X Identifier Direct™ Software Panel and Bin files may be imported into GeneMapper® ID-X Software v1.0.1 or later to assist in automated genotyping of data generated using the AmpFISTR® Identifier Direct™ kit chemistry. Also provided are Analysis Method and Size Standard files for import.

Note: These files have been provided to assist you in getting started with Identifier Direct™ kit data analysis. Validation of these files should be established by each individual laboratory based on the laboratory's internal validation studies.

-----  
SOFTWARE CONTENTS  
-----

This software package consists of 5 files:

- CE\_G5\_IdentifierDirect\_GS500.xml
- IdentifierDirect\_GS500\_Bins\_v1X.txt
- IdentifierDirect\_GS500\_Panels\_v1X.txt
- IdentifierDirect\_GS500\_Stutter\_v1X.txt
- IdentifierDirect\_v1X.xml

-----  
VALIDATION SUMMARY  
-----

Testing was carried out utilizing two capillary electrophoresis instruments (3730 Data Collection v3.0 and 3130xl Data Collection v3.0). Concordance testing was performed between GeneMapper® ID-X v1.1 and GeneMapper® ID-X v3.2.1 softwares on 200 blood on FTA samples, amplified using the Identifier Direct™ and Identifier® kits.

NOTE: The primer sequences for the previously validated chemistry, Identifier®, are the same primer sequences used in the Identifier Direct™ kit. The stutter values for the Identifier Direct™ kit have been determined experimentally and are different from the stutter values determined for the Identifier® Kit.

-----  
FILE INSTALL INSTRUCTIONS  
-----

Place the five files in the software package in the appropriate folder on the computer. If you used the default install configuration, this default Panels folder should be located at C:\AppliedBiosystems\GeneMapper\Panels

Installation of the Panel File:

#### ReadMe\_IDDirect\_v1x.txt

1. In the GeneMapper® ID-X software, select Tools from the menu bar
2. Select Panel Manager
3. Single Click on "Panel Manager" in the left hand window pane
4. Select File and Import Panels from the menu bar
5. A dialogue box will appear. Navigate the location of the Panel file on your computer. If you placed them in the default Panels folder, the dialogue box should open to the correct folder.
6. Select the file titled: IdentifierDirect\_GS500\_Panels\_v1x.txt
7. Select the appropriate Security Group
8. This will install the Panel in the left window pane

#### Installation of the Bin File:

1. Single click on "IdentifierDirect\_GS500\_v1x" in the left hand window pane
2. Select File and Import Bin Set from the menu bar
3. A dialogue box will appear. Navigate to the location of the Bin file on your computer. If you placed them in the default Panels folder, the dialogue box should open to the correct folder.
4. Select the file titled: IdentifierDirect\_GS500\_Bins\_v1X.txt
5. This will install the Bins in the pull down menu at the top titled "Bin Set"

#### Installation of the Stutter File:

1. Single click on "IdentifierDirect\_GS500\_v1x" in the left hand window pane
2. Select File and Import Marker Stutter from the menu bar
3. A dialogue box will appear. Navigate to the location of the Bin file on your computer. If you placed them in the default Panels folder, the dialogue box should open to the correct folder.
4. Select the file titled: IdentifierDirect\_GS500\_Stutter\_v1X.txt
5. This will install the Stutter file

#### Installation of the Analysis Methods File:

1. Navigate to Tools/Genemapper Manager/Analysis Methods
2. Select the import button on the bottom of the window.
3. A dialogue box will appear. Navigate to the location of the Analysis Methods file on your computer.
4. Select the file titled: IdentifierDirect\_v1X.xml
5. Select the appropriate Security Group
6. This will install the method in the list of analysis methods.

#### Installation of the Size Standards File:

1. Navigate to Tools/Genemapper Manager/Size Standards
2. Select the import button on the bottom of the window.
3. A dialogue box will appear. Navigate to the location of the Size Standards file on your computer.
4. Select the file titled: CE\_G5\_IdentifierDirect\_GS500.xml
5. Select the appropriate Security Group
6. This will install the standard in the list of size standards.

#### ----- LEGAL DISCLAIMER -----

For Research, Forensic or Paternity Use Only. Not for use in diagnostic procedures.

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

Notice to Purchaser: License Disclaimer.

Purchase of this software product alone does not imply any license under any process, instrument or other apparatus, system, composition, reagent or kit rights

ReadMe\_IDDirect\_v1x.txt

under patent claims owned or otherwise controlled by Life Technologies Corporation, either expressly, or by estoppel.

Applied Biosystems, AmpFISTR® and GeneMapper® are registered trademarks, and the AB Design and Identifier Direct™ are trademarks of Life Technologies Corporation or its subsidiaries in the U.S. and/or certain other countries.

GeneMapper® ID-X Software v1.0 has undergone a verification process defined by Applied Biosystems. However, human identification laboratories analyzing forensic, paternity, databasing, and single-source samples that choose to use GeneMapper® ID-X Software for data analysis should perform their own appropriate validation studies.

(C) 2009 Applied Biosystems. All rights reserved.