

Communication Patch 1

Early Expiration of Internal Password Affecting:

3130-Series Data Collection Software v3.1

3730-Series Data Collection Software v3.1

GeneMapper® Software v4.1

These instructions describe the installation of the GM_UDC_Communication_Patch1 which remedies an issue with an internal communication connection between AB software and embedded software components. The following AB products and product configurations are affected (Full and Upgrade installations):

- 3130 Data Collection Software v3.1
- 3130xl Data Collection Software v3.1
- 3730/3730xl Data Collection Software v3.1
- GeneMapper Software v4.1
 - Installed on a standalone computer
 - Co-installed with any of the above Data Collection Software v3.1 on the data collection computer of 3130/3130xl and 3730/3730xl Genetic Analyzers
 - Installed on the data collection computer of a 3500-Series Genetic Analyzer
 - Installed on the data collection computer of a 310 Genetic Analyzer with Data Collection Software v3.1

The following warning will be displayed when the error is experienced:

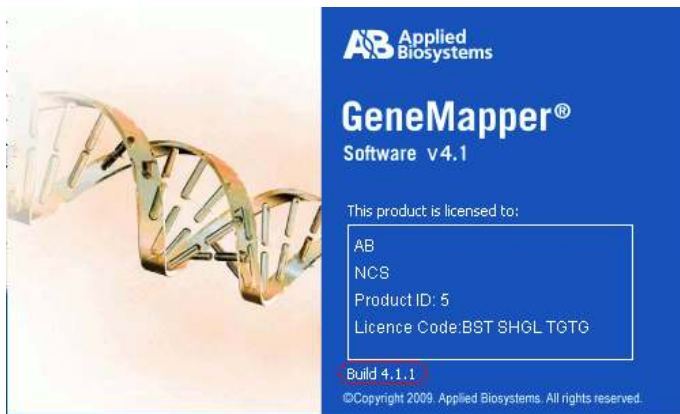
"The connection to the default database has failed"

GM_UDC_Communication_Patch1.exe can be run at any time with no need to wait for the error message to display. Close the software before running the Patch1 executable file. Once the Patch is installed, there is no need to re-install any of the software.

GM_UDC_Communication_Patch1 installation instructions:

1. Log in as Administrator
2. Close software:
 - Close GeneMapper® Software v4.1
 - Close 3130- or 3730-Series Data Collection Software v3.1
 - Close 3500-Series Data Collection Software
 - Close 310 Data Collection Software v3.1
3. Take the following steps to ensure that all Oracle Corporation software services are running:
 - a. Go to **Start > Control Panel > Administrative Tools > Services**
 - b. Make sure that the following services have the status 'Started'
 - OracleIFATNSListener or OracleUDCTNSListener
 - OracleServiceIFA or OracleServiceUDC
 - c. If the above services do not have the status 'Started', select the service, right click and select **Start** to start the services.
4. Double click on GM_UDC_Communication_Patch1.exe to start the installation process. Follow the instructions provided by the installer until the installation is complete.
5. After installing the Patch, please read the release notes.

- For GeneMapper4.1, the build number 4.1.1 will display in the About dialog box. The GeneMapper® Software v4.1.1 Release Notes.doc is located under x:\AppliedBiosystems\GeneMapper\Docs



- For Data Collection Software (3130- and 3730-Series) v3.1, the build number 3.1.1 will display on the Viewer Panel. The Data Collection Software v3.1.1 Release Notes.doc is located under x:\AppliedBiosystems\UDC\Documents



For Research Use Only. Not intended for any animal or human therapeutic or diagnostic use.