

# Use of the MSI Analysis System with the Applied Biosystems® 3500 Genetic Analyzer and POP-7® Polymer

## MSI Analysis System Application Note

Promega Corporation

### Notes:

The MSI Analysis System is  
For Research Use Only.  
Not For Use In Diagnostic Procedures.

*Guidelines for detecting  
microsatellite instability  
using the MSI Analysis  
System, Applied Biosystems®  
3500 Genetic Analyzer and  
POP-7® Polymer*

### 1. Introduction

The MSI Analysis System was developed and optimized for use with the Applied Biosystems® Genetic Analyzers using POP-4® polymer and a 36cm capillary array. However, amplification products can be detected using other instrument configurations. This Application Note provides general guidelines for use of the MSI Analysis System with the Applied Biosystems® 3500 and 3500xL Genetic Analyzers and POP-7® polymer. Optimization may be required for individual research needs.

### 2. General Considerations for Amplification

The MSI Analysis System is optimized to amplify 1–2ng of genomic DNA in a 10µl reaction volume using the GeneAmp® PCR System 9600 and 9700 thermal cyclers. See the *MSI Analysis System, Version 1.2, Technical Manual #TM255* for amplification conditions.


For thermal cyclers other than the GeneAmp® PCR System 9600 and 9700, we recommend that you replicate as closely as possible the thermal cycling conditions described in the *MSI Analysis System, Version 1.2, Technical Manual #TM255*. Suboptimal thermal cycling conditions can result in imbalanced profiles. Contact Promega Technical Services at [genetic@promega.com](mailto:genetic@promega.com) for assistance with other thermal cyclers.


### 3. Materials to Be Supplied by the User

- MSI Analysis System, Version 1.2 (Cat.# MD1641)
- PowerPlex® 4C Matrix Standard (Cat.# DG4800)

### 4. Spectral Calibration Using the PowerPlex® 4C Matrix Standard

The PowerPlex® 4C Matrix Standard (Cat.# DG4800) contains matrix fragments labeled with four fluorescent dyes: Fluorescein, JOE, TMR and CXR. Once generated, the spectral calibration file is applied during collection of MSI Analysis System data to calculate and compensate for spectral overlap between different fluorescent dye colors. These matrix standards are compatible with the Applied Biosystems® 3500 and 3500xL Genetic Analyzers. For more information, see the *PowerPlex® 4C Matrix Standard Technical Manual #TMD048* or the *Applied Biosystems 3500/3500xL Genetic Analyzer User Guide*.

 The quality of formamide is critical. Use Hi-Di™ formamide. Freeze formamide in aliquots at –20°C. Multiple freeze-thaw cycles or long-term storage at 4°C may cause breakdown of formamide. Poor-quality formamide may contain ions that compete with DNA during injection, which results in lower peak heights and reduced sensitivity. A longer injection time may not increase the signal.

 Formamide is an irritant and a teratogen; avoid inhalation and contact with skin. Read the warning label, and take the necessary precautions when handling this substance. Always wear gloves and safety glasses when working with formamide.

#### 4.A. Matrix Sample Preparation

There may be instrument-to-instrument variation in the sensitivity of detection. The dilutions described here may need to be optimized in individual laboratories depending on the sensitivity of each Applied Biosystems® 3500 or 3500xL Genetic Analyzer.

**Note:** We have found that the use of fresh polymer and a new capillary array results in an optimal spectral calibration.


1. Thaw the matrix standards. Mix each matrix standard by vortexing for 5–10 seconds prior to use. Do not centrifuge the matrix standards as this may cause the DNA to be concentrated at the bottom of the tube.

**Note:** While the matrix dyes thaw, preheat the oven of the capillary electrophoresis instrument to 60°C as described in Section 4.B, Step 1.


2. Vortex the 4C Matrix Mix for 10–15 seconds prior to use. Add 10µl of 4C Matrix Mix to one tube of Matrix Dilution Buffer. Vortex for 10–15 seconds. Note the date of dilution on the tube.

**Note:** The diluted 4C Matrix Mix can be stored for up to 1 week at 2–10°C.

3. Add 10µl of the diluted 4C Matrix Mix prepared in Step 2 to 500µl of Hi-Di™ formamide. Vortex for 10–15 seconds.
4. For the Applied Biosystems® 3500xL Genetic Analyzer, 24 wells are used for spectral calibration on 24 capillaries (wells A1 through H3 of a 96-well plate). Add 15µl of 4C Matrix Mix with formamide prepared in Step 3 to each of the 24 wells. After placing the septa on the plate, briefly centrifuge the plate to remove any air bubbles.

 Do not heat denature.

For the Applied Biosystems® 3500 Genetic Analyzer, 8 wells are used for spectral calibration on 8 capillaries (wells A1 through H1 of a 96-well plate). Add 15µl of 4C Matrix Mix with formamide prepared in Step 3 to each of the eight wells. After placing the septa on the plate, briefly centrifuge the plate to remove any air bubbles.

 Do not heat denature.

5. Place the plate in the 3500 series 96-well standard plate base, and clamp the plate retainer squarely in place over the plate and base until it clicks twice. Place the plate assembly in Position A on the autosampler with the labels facing you.

## 4.B. Instrument Preparation

1. If you have not already preheated the oven to 60°C, select “Start Pre-Heat” (Figure 1). Applied Biosystems recommends that you preheat the oven for at least 30 minutes before you start a run if the instrument is cold.

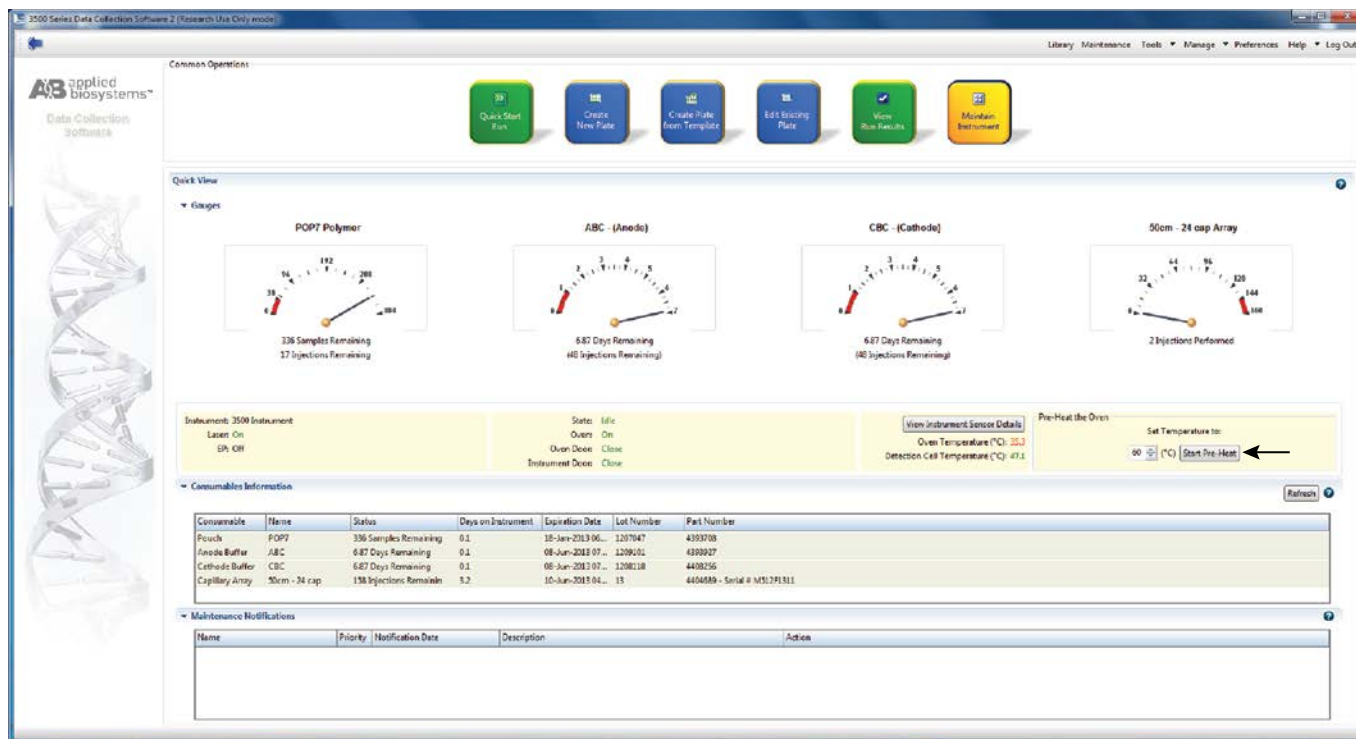


Figure 1. The Dashboard.

2. To perform a spectral calibration with the Promega 4-dye chemistry for the first time, create a new dye set.
  - a. To create the new dye set, navigate to the Library, highlight “Dye Sets” and select “Create” (Figure 2).
  - b. The Create New Dye Set window will appear. Name the Dye Set (e.g., Promega 4Dye), select “Matrix Standard” for the Chemistry and select “F Template” for the Dye Set Template. Leave all other settings as the default values. You may lock this dye set to avoid making unintended changes. Select “Save” and “Close”.

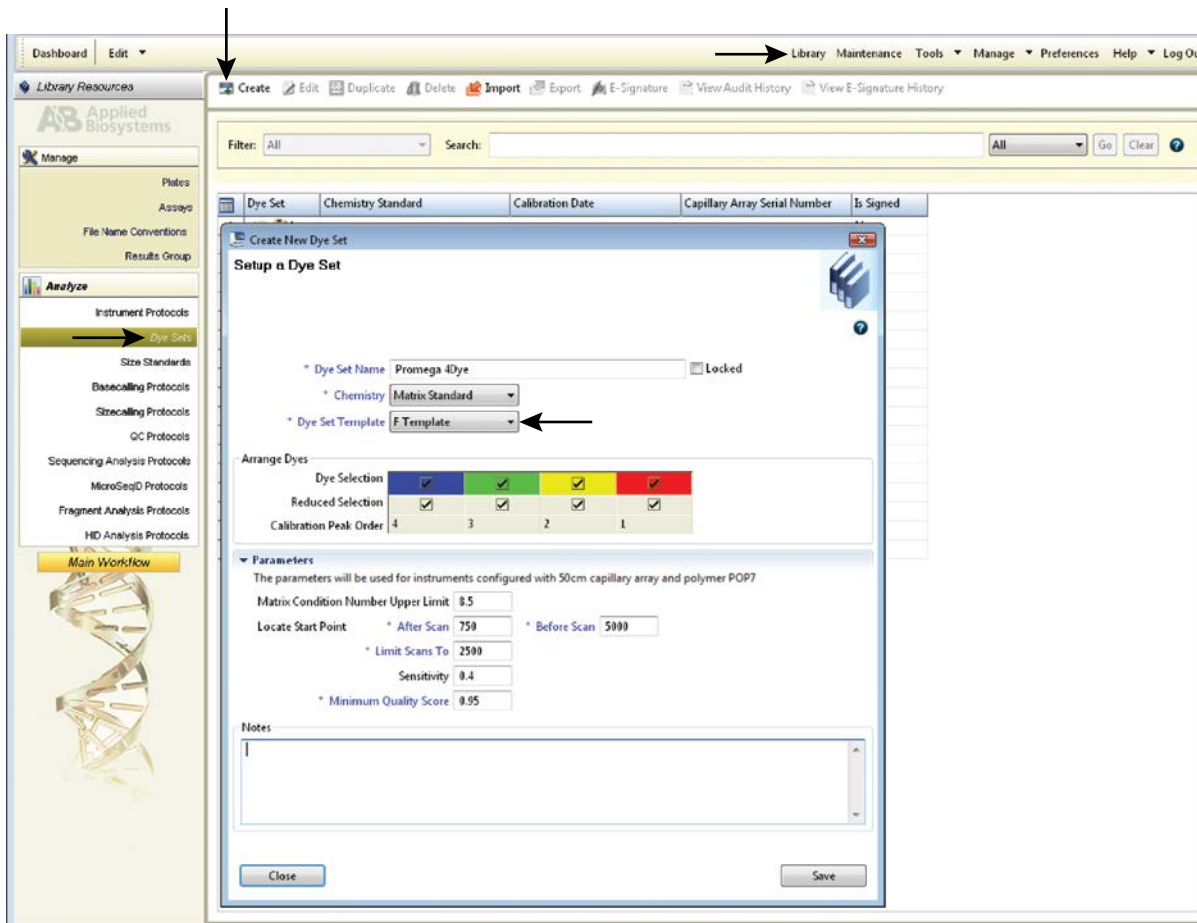


Figure 2. The Create New Dye Set window.

- c. To perform the spectral calibration with the Promega 4-dye chemistry, choose “Maintenance” from the Menu bar, then choose “Spectral” in the navigation pane. Under the Calibration Settings choose “Matrix Standard” as the Chemistry Standard, and select the name of the dye set you created in Step 2.b. Check the Allow Borrowing checkbox. Select other options as shown in Figure 3.
- d. Select “Start Run”.

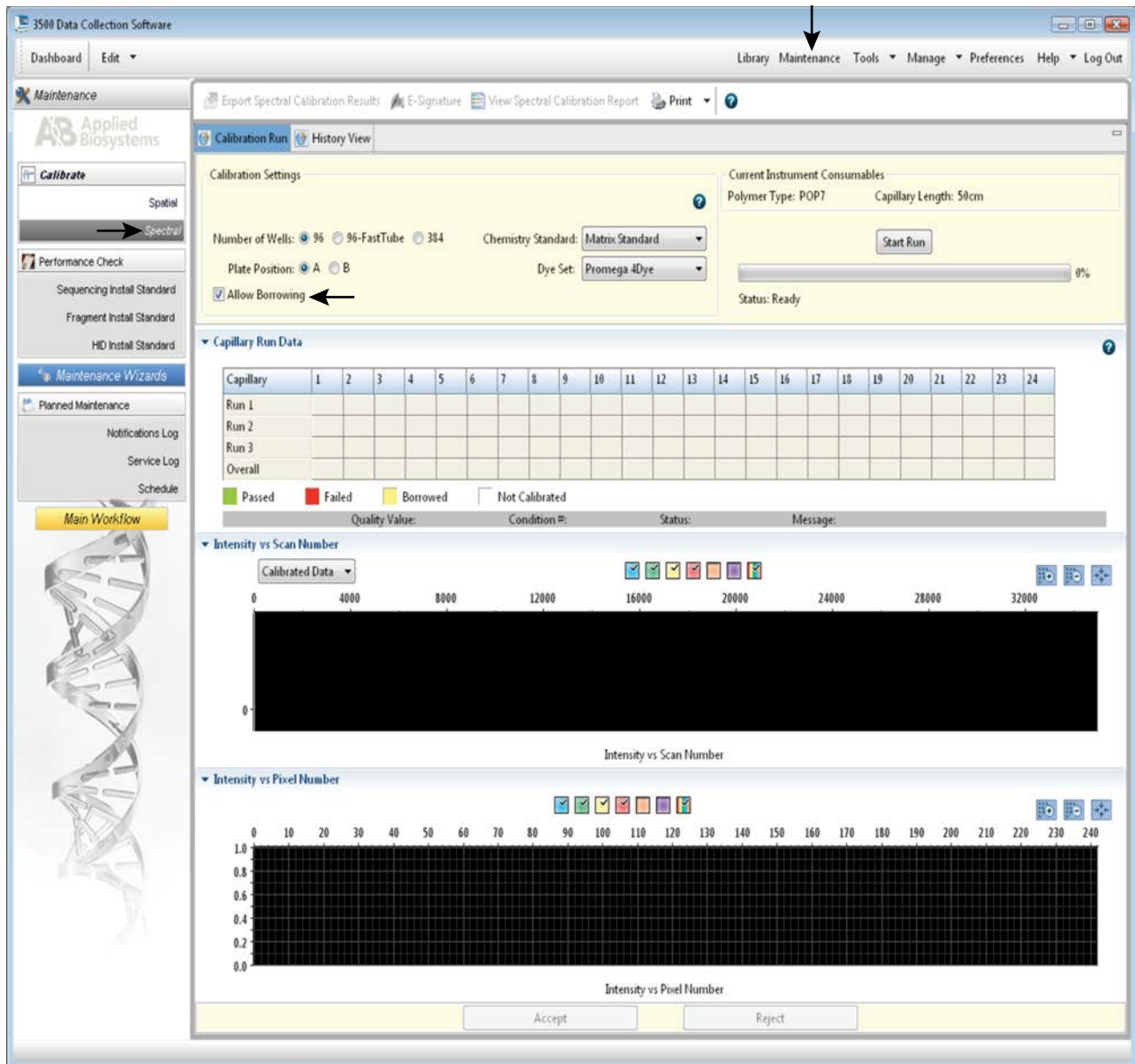


Figure 3. Calibration Run.

- If fewer than the recommended number of capillaries pass, the spectral calibration run will be repeated automatically up to two more times. Upon completion of the spectral calibration, check the quality of the spectral calibration in the Capillary Run Data display (Figure 4). Choose either “Accept” or “Reject” (not shown).

**Note:** Refer to the *3500 Series Data Collection Software User Manual* for the criteria recommended by Applied Biosystems when accepting or rejecting a spectral calibration. The data collection software will display the results for passing and failed capillaries. If the spectral calibration fails, see the Troubleshooting section of the *PowerPlex® 4C Matrix Standards Technical Manual* #TMD048.

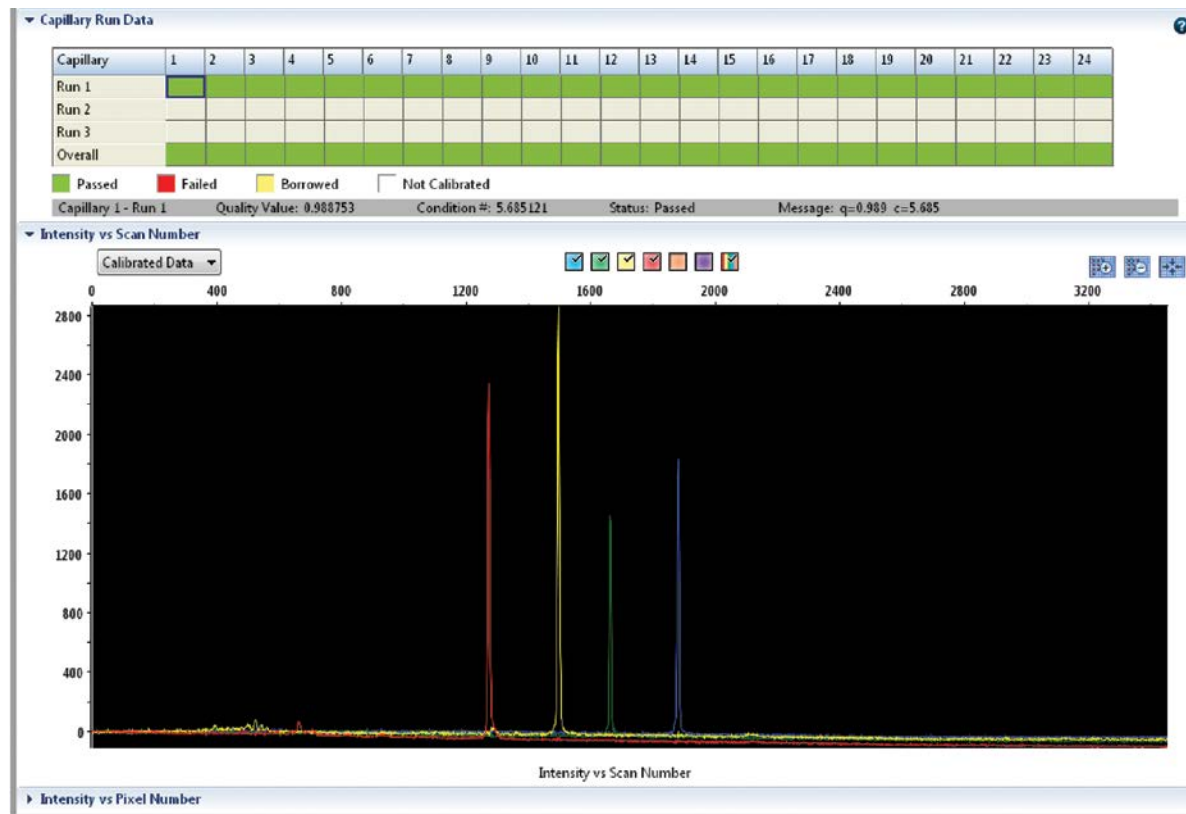




Figure 4. The Capillary Run Data display.

## 5. Detection of Amplified Fragments Using the Applied Biosystems® 3500 or 3500xL Genetic Analyzer

-  The quality of formamide is critical. Use Hi-Di™ formamide. Freeze formamide in aliquots at –20°C. Multiple freeze-thaw cycles or long-term storage at 4°C may cause breakdown of formamide. Poor-quality formamide may contain ions that compete with DNA during injection, which results in lower peak heights and reduced sensitivity. A longer injection time may not increase the signal.
-  Formamide is an irritant and a teratogen; avoid inhalation and contact with skin. Read the warning label, and take appropriate precautions when handling this substance. Always wear gloves and safety glasses when working with formamide.

### 5.A. Sample Preparation

Before preparing the samples, preheat the oven of the capillary electrophoresis instrument to 60°C as described in Section 5.B, Step 1, if desired.

1. Determine the number of samples to be analyzed, and add additional reactions to this number to compensate for pipetting error. Prepare a loading cocktail by combining and mixing Internal Lane Standard 600 (ILS 600) and Hi-Di™ formamide as follows:  
[(0.5µl ILS 600) × (# samples)] + [(9.5µl Hi-Di™ formamide) × (# samples)]

**Note:** The volume of internal lane standard used in the loading cocktail can be increased or decreased to adjust the intensity of the size standard peaks.

2. Vortex for 10–15 seconds to mix.
3. Pipet 10µl of loading cocktail into each well.
4. Add 1µl of amplified sample. Cover wells with appropriate septa.

**Note:** Instrument detection limits vary; therefore, injection time or injection voltage may need to be increased or decreased. To modify the injection time or injection voltage in the run module, select “Instrument Protocol” from the Library menu in the data collection software.

5. Centrifuge plate briefly to remove air bubbles from the wells.
6. Denature samples at 95°C for 3 minutes, then immediately chill on crushed ice or in an ice-water bath for 3 minutes. Denature samples just prior to loading the instrument.

**Note:** If using a thermal cycler to denature samples, do not close the heated lid as this may melt the plate septa.

### 5.B. Instrument Preparation

Refer to the *Applied Biosystems 3500/3500xL Genetic Analyzer User Guide* for the instrument maintenance schedule and instructions to install the capillary array, buffers and polymer pouch and perform a spatial calibration. This Application Note provides general guidelines for use of the MSI Analysis System with the Applied Biosystems® 3500 Genetic Analyzers and POP-7® polymer. Some optimization may be required for individual research needs.

1. Open the 3500 Data Collection Software. The Dashboard screen will launch (Figure 1). Ensure that the Consumables Information and Maintenance Notifications are acceptable. Set the oven temperature to 60°C, then select “Start Pre-Heat”. Applied Biosystems recommends that you preheat the oven for at least 30 minutes before you start a run if the instrument is cold.



- To set up an Instrument Protocol, navigate to the Library, select “Instrument Protocols” from the navigation pane, and choose “Fragment” as the Filter. Choose the default protocol that is the most similar to the one you will be using from the list of protocols, and duplicate it (Figure 5). Do not choose “Create”. For example, if you are using an 8-capillary 3500 Genetic Analyzer with a 50cm array and POP-7<sup>®</sup> polymer, choose “FragmentAnalysis50\_POP7\_1.”, and select “Duplicate”. Edit the copy of this Instrument Protocol, and fill in the drop-down menus as appropriate for your configuration. Make sure you choose “Fragment” as the Application Type, and select the dye set that you created in Section 4.B. The Run Module will be “FragmentAnaysis50\_POP7”. Assign the Instrument Protocol a meaningful name so that it will be clear what the Instrument Protocol can be used for in future runs. Leave all run options and Advanced Options as the default values. It is possible to change default injection parameters later if necessary later. Select “Save” and “Close”.

When you save this protocol with the Promega 4Dye dye set, the Instrument Protocol name will appear in the Library with the Dye Set Name associated with it.

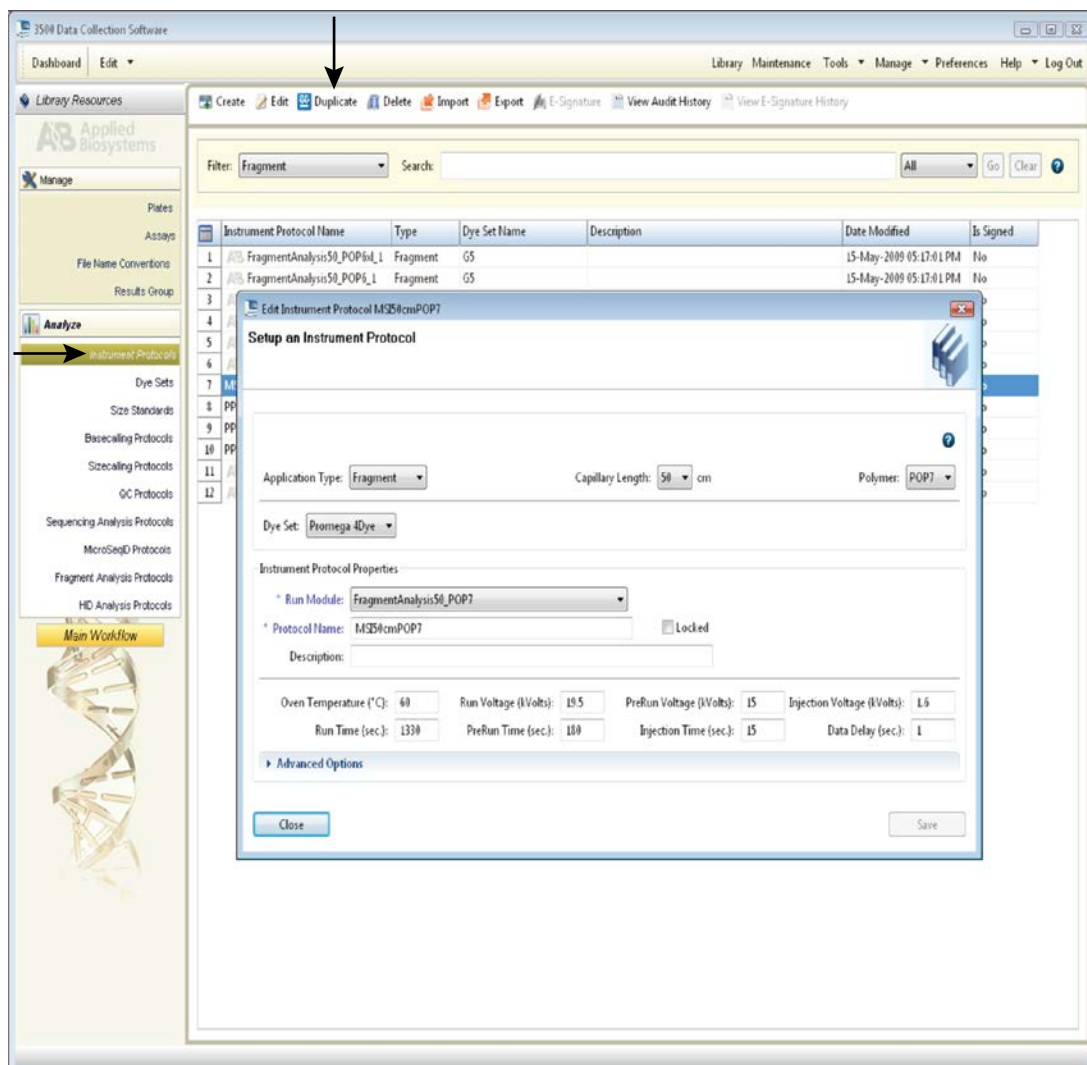


Figure 5. The Setup an Instrument Protocol window.



3. To create a new Size Standard for the Sizecalling protocol, navigate to the Library. Select “Size Standards”, then select “Create”. Assign the size standard an appropriate name (Figure 6). Choose “Red” for the Dye Color. Enter the fragment sizes for the size standard: 60, 80, 100, 120, 140, 160, 180, 200, 225, 250, 275, 300, 325, 350, 375, 400, 425, 450, 475, 500, 550 and 600 bases. Select “Add Sizes”, “Save” then “Close”.

**Note:** Definition and detection of the 600bp fragment is optional.

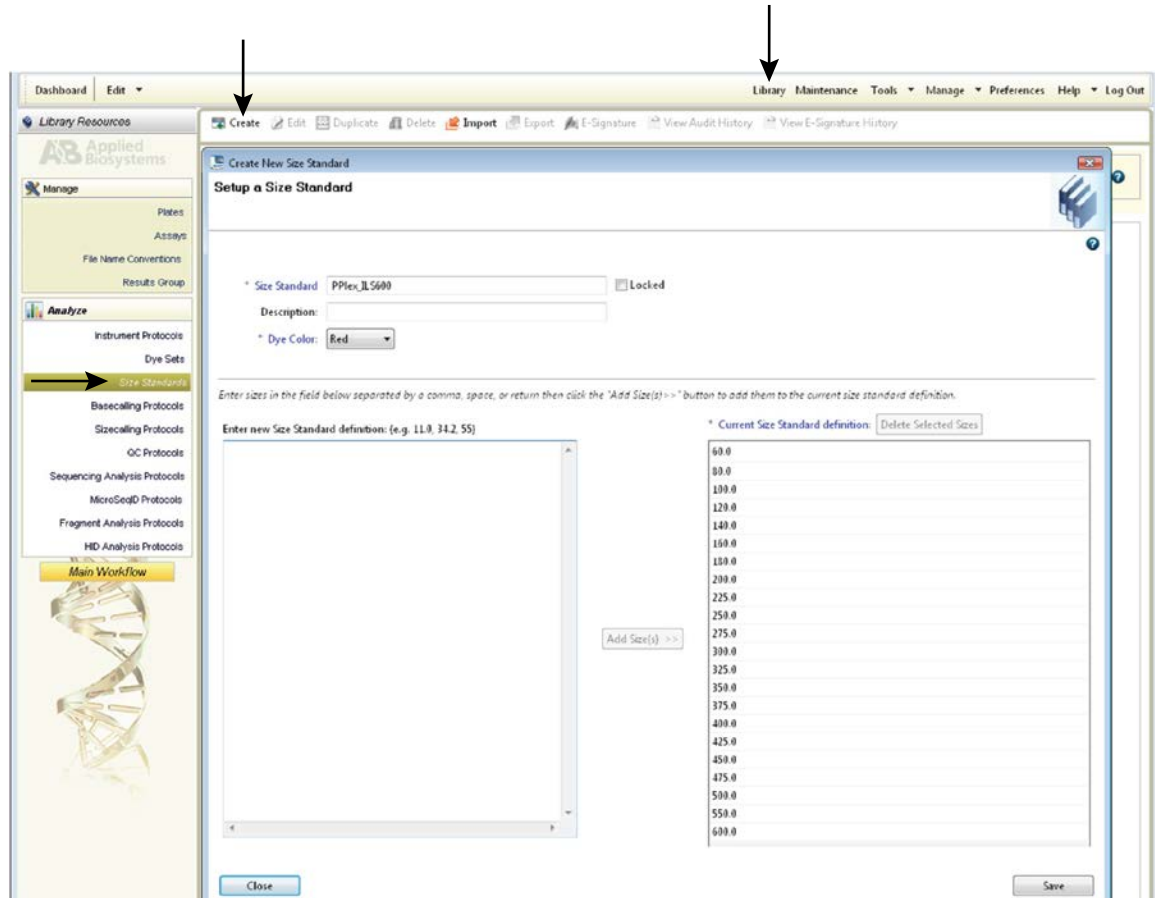


Figure 6. The Create New Size Standard window.

- To create a new Sizecalling Protocol, navigate to the Library. Select “Sizecalling Protocols”, then select “Create”. Assign a descriptive protocol name (Figure 7). Select the size standard created in Step 3. Leave all other settings as default. Select “Save” then “Close”.

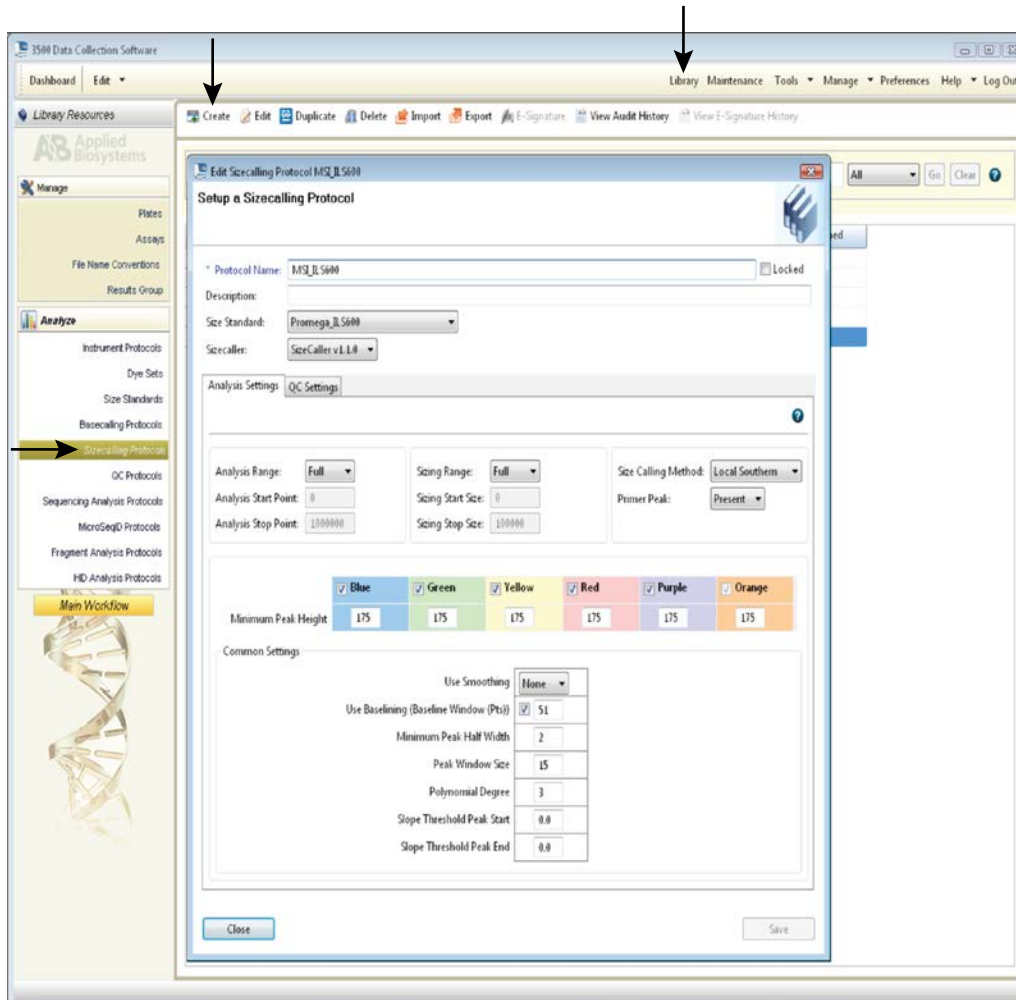


Figure 7. The Create New Sizecalling Protocol window.

5. To create a new Assay, navigate to the Library. Select “Assays”, then select “Create”. In the Assay window (Figure 8), select the application type “Fragment” and the appropriate Instrument Protocol and Sizecalling Protocol. Assign a descriptive assay name, then select “Save” and “Close”.

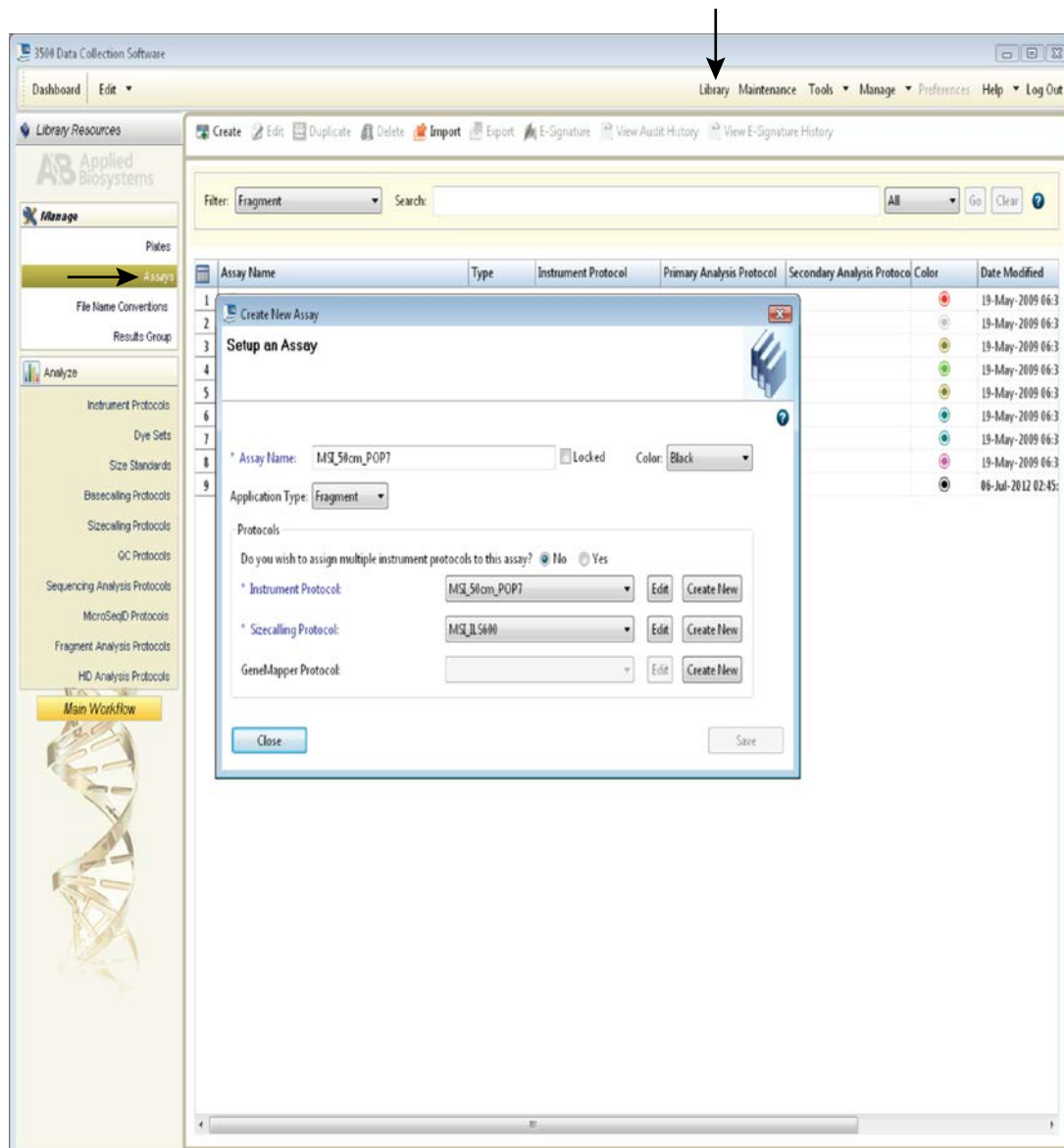


Figure 8. The Assay window.

- To create a new File Name Convention, navigate to the Library. Select “File Name Conventions”, then select “Create” (Figure 9). Select the File Name Attributes according to laboratory practices, and save with a descriptive name. Under “Select File Location” choose “Default File Location”, or choose “Custom File Location” and navigate to your custom location. Select “Save” then “Close”.

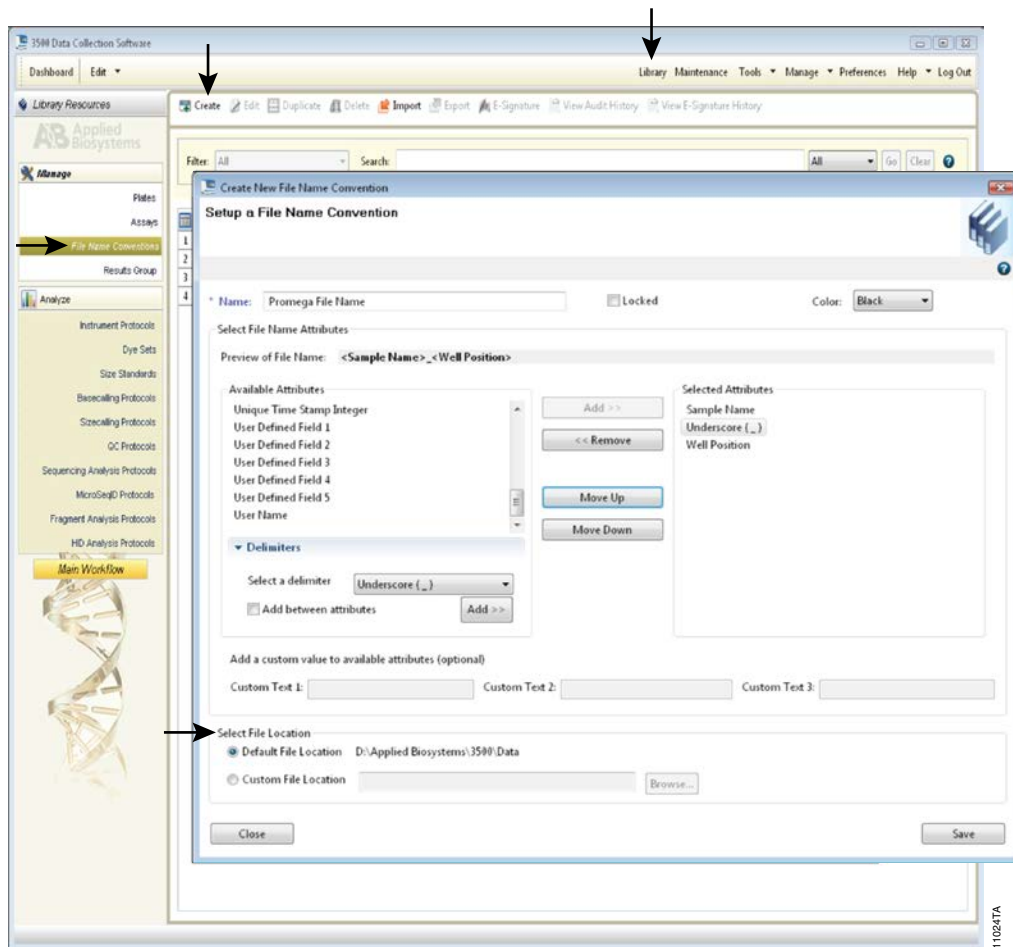


Figure 9. The Create New File Name Convention window.

7. To create a new Results Group, navigate to the Library. Select “Results Group”, then select “Create” (Figure 10). Select Results Group Attributes according to laboratory practices. Save with a descriptive name, and select “Close”.

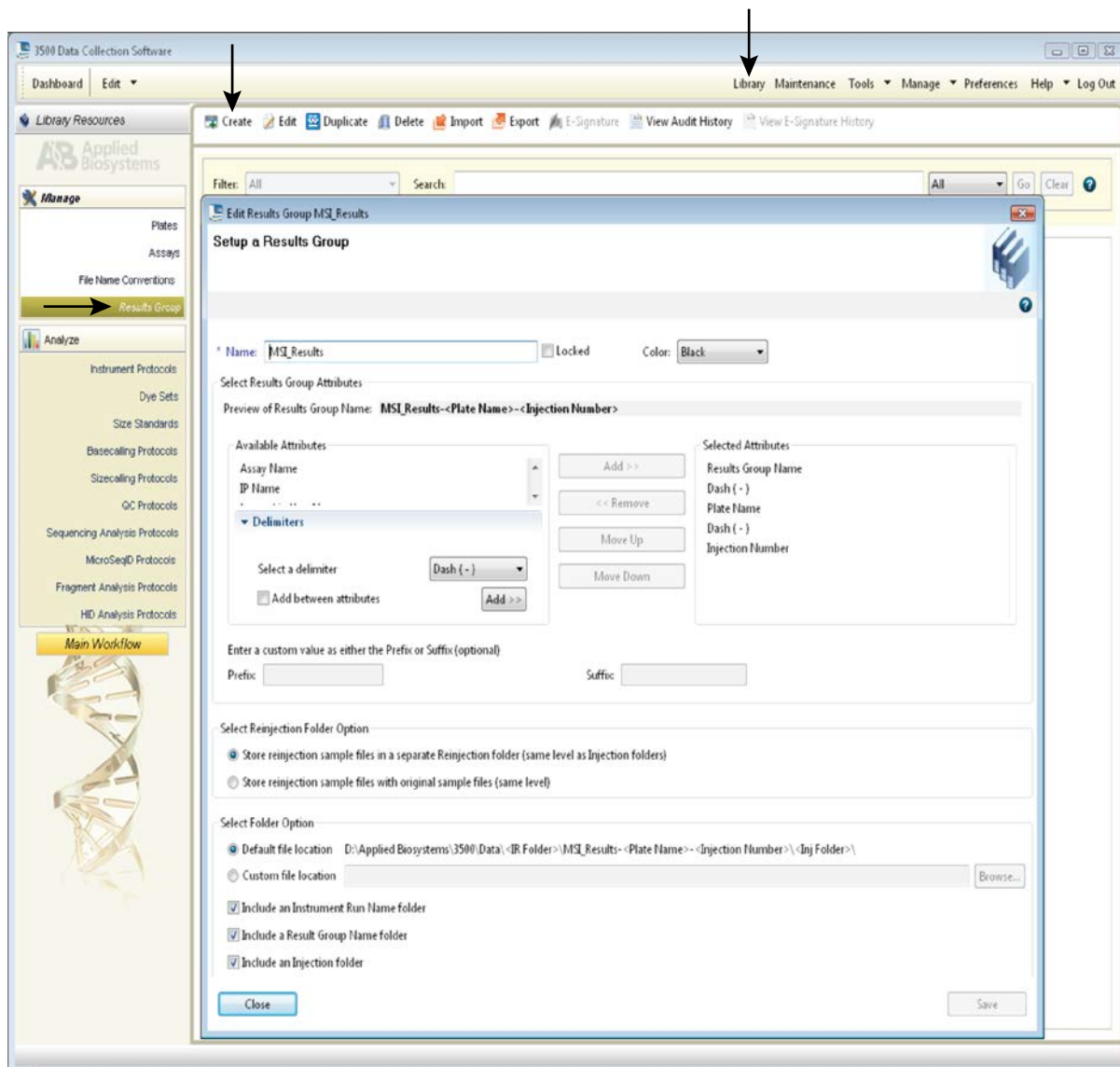
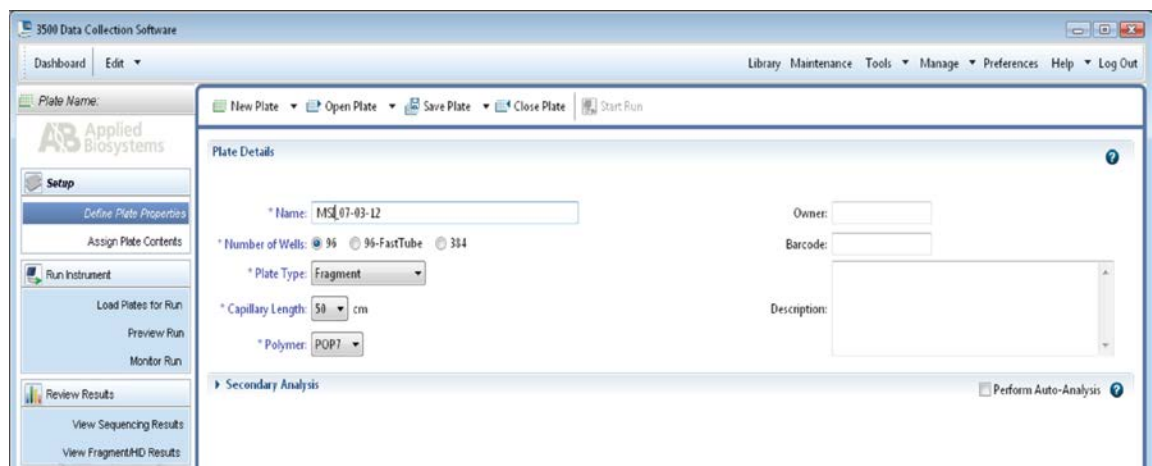


Figure 10. The Create New Results Group window.

- To create a New Plate, navigate to the Library, and from the Manage menu, select “Plates”, then “Create”.
- Assign a descriptive plate name. Select the Plate Type “Fragment” from the drop-down menu (Figure 11). Select the appropriate capillary length and polymer.
- Select “Assign Plate Contents”.



**Figure 11. Defining plate properties.**

- Assign sample names to wells.

- In the lower left portion of the screen (Figure 12), under “Assays”, use the Add from Library option (not shown) to select the Assay created in Step 5. Click on the Add to Plate button, and close the window.

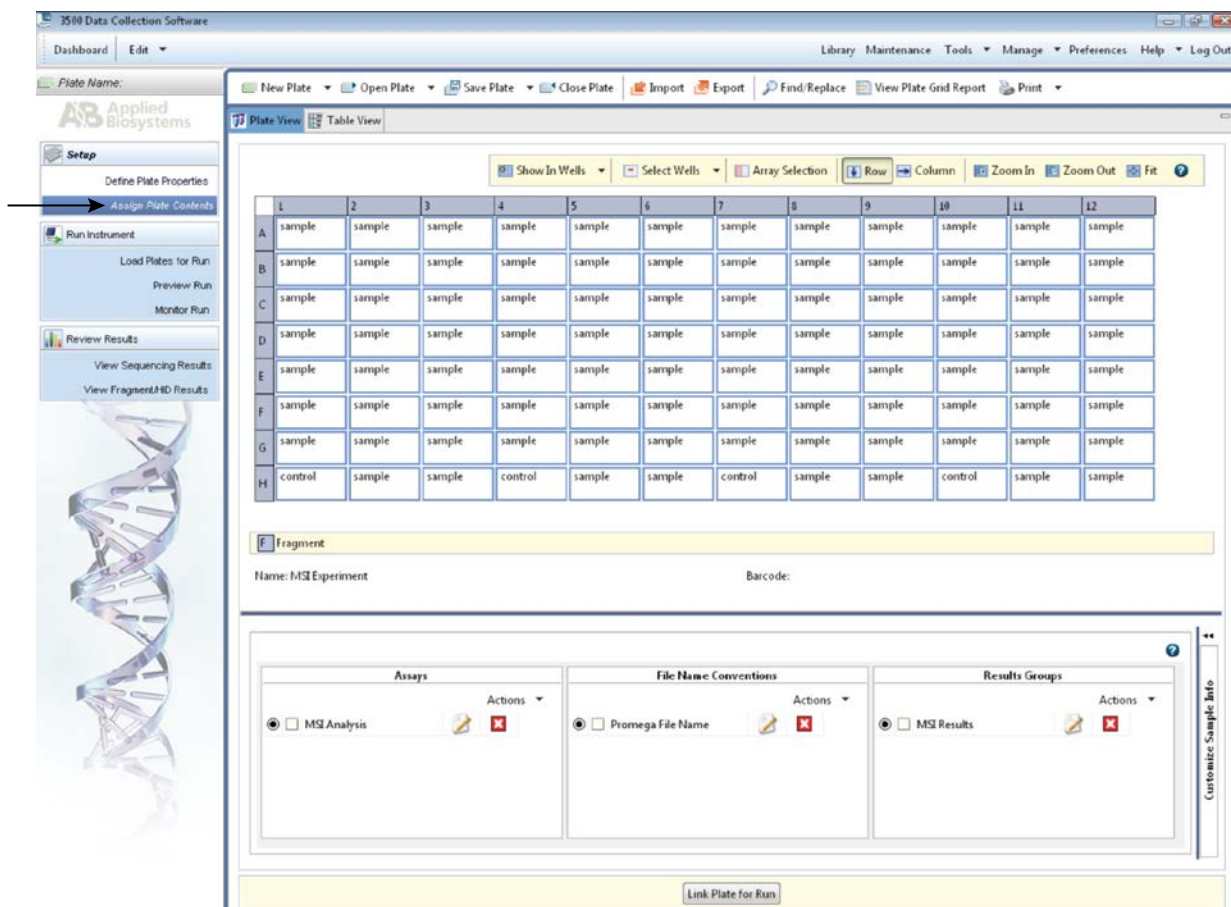


Figure 12. Assigning plate contents.

- Under “File Name Convention”, use the Add from Library option to select the File Name Convention created in Step 6. Click on the Add to Plate button, and close the window.
- Under “Results Groups”, use the Add from Library option to select the Results Group created in Step 7. Click on the Add to Plate button, and close the window.
- Highlight the sample wells, then select the boxes in the Assays, File Name Conventions and Results Groups that pertain to those samples. To assign sample types, select the Custom Sample Info tab at the bottom left side of the window, assign sample types, then save the plate.
- Select “Link Plate for Run”.
- The Load Plate window will appear. Select “Yes”.
- In the Run Information window, assign a Run Name. Select “Start Run”.



## 5.C. Data Analysis

This section describes how to obtain information for the analysis of MSI Analysis System data generated using the Applied Biosystems® 3500 series of genetic analyzers with GeneMapper® software, version 4.1 or 5.0. Note that Promega scientists have not verified the performance of GeneMapper® software, version 4.1 or 5.0, with MSI Analysis System data generated using a 50cm capillary and POP-7® polymer. Some optimization may be required.

1. To analyze .fsa files collected on the Applied Biosystems® 3500 series of genetic analyzers using GeneMapper® software, version 4.1, download the panels and bins text files at:

**[www.promega.com/resources/software-firmware/msi-panels-and-bins-for-gemapper-software/](http://www.promega.com/resources/software-firmware/msi-panels-and-bins-for-gemapper-software/)**

Install the panels and bins text files as described in the *MSI Analysis System, Version 1.2, Technical Manual #TM255*.

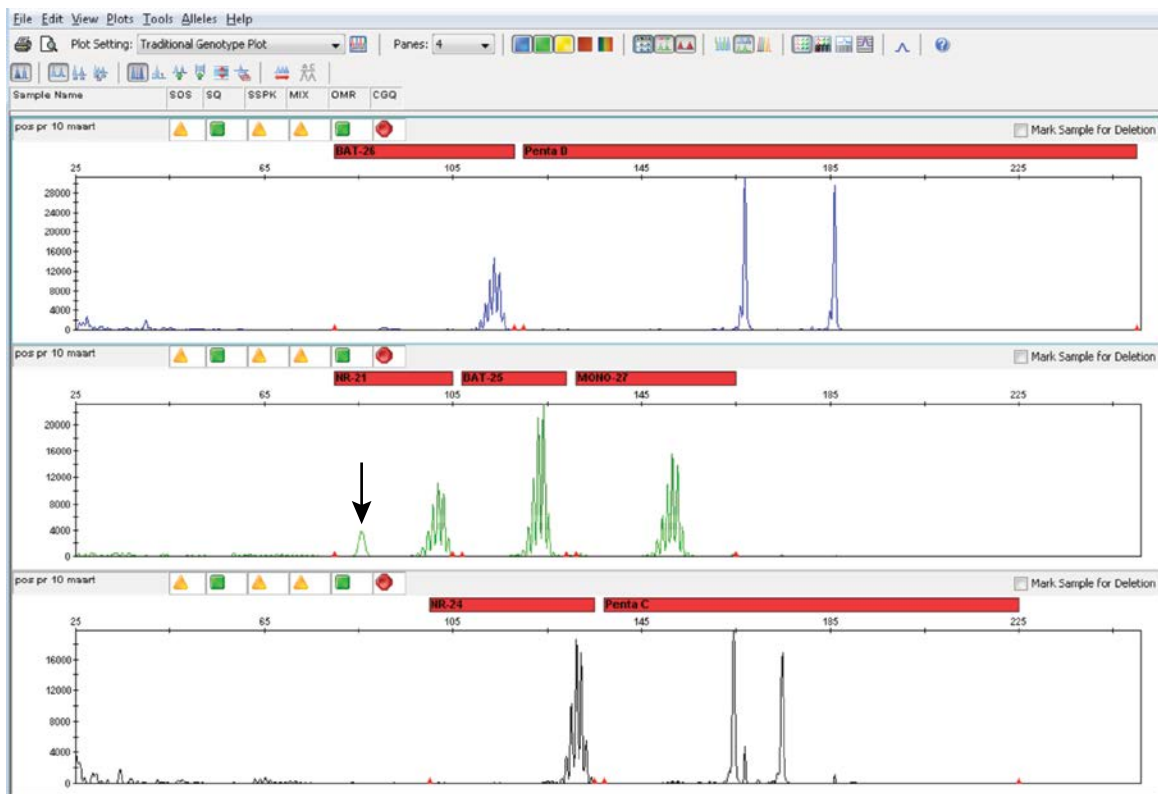
For data import and data processing, follow the instructions provided in the *MSI Analysis System, Version 1.2, Technical Manual #TM255*.

For detailed data analysis instructions refer to the *Applied Biosystems GeneMapper® 4.1 User Guide*.

2. To analyze .fsa files collected on the Applied Biosystems® 3500 series of genetic analyzers using GeneMapper® ID or GeneMapper® ID-X software, contact Promega Technical Services by e-mail at: [genetic@promega.com](mailto:genetic@promega.com) to obtain the appropriate panels and bins text files.

## 5.D. Known Artifacts

We frequently observe an artifact at 87–90bp in the JOE channel when using POP-7® polymer with the Applied Biosystems® 3500 series Genetic Analyzers (Figure 13) and Applied Biosystems® 3130 series Genetic Analyzers. However, the size of this artifact is considerably different than the sizes of the repeat markers and, as a result, should not interfere with data interpretation. Other nonspecific artifacts may occur depending on the detection conditions. See the troubleshooting section of the *MSI Analysis System, Version 1.2, Technical Manual #TM255* for more information.



**Figure 13.** An example of an artifact observed when analyzing MSI Analysis System data with an Applied Biosystems® 3500 Genetic Analyzer and POP-7® polymer.

# MSI Analysis System

---

## Ordering Information

Product	Size	Cat.#
MSI Analysis System, Version 1.2*	100 reactions	MD1641
PowerPlex® 4C Matrix Standard**	5 preps	DG4800

\*For Research Use Only. Not for Use in Diagnostic Procedures.

\*\*Not For Medical Diagnostic Use.

PowerPlex is a registered trademark of Promega Corporation.

Applied Biosystems, GeneAmp and GeneMapper are registered trademarks of Applied Biosystems. Hi-Di is a trademark of Applied Biosystems. POP-4 and POP-7 are registered trademarks of Thermo Fisher Scientific.

Products may be covered by pending or issued patents or may have certain limitations. Please visit our Web site for more information.

