



**Promega**

## Technical Bulletin

---

# PowerPlex<sup>®</sup> 5-Dye Matrix Standards, 310

INSTRUCTIONS FOR USE OF PRODUCT DG4600.

### IMPORTANT

We recommend using this product once, then discarding it.



[www.promega.com](http://www.promega.com)

# PowerPlex<sup>®</sup> 5-Dye Matrix Standards, 310

All technical literature is available on the Internet at: [www.promega.com/tbs/](http://www.promega.com/tbs/)  
Please visit the web site to verify that you are using the most current version of this  
Technical Bulletin. Please contact Promega Technical Services if you have questions on use  
of this system. E-mail: [techserv@promega.com](mailto:techserv@promega.com)

1. Description.....	2
2. Product Components and Storage Conditions .....	2
3. Detection of Matrix Fragments Using the ABI PRISM <sup>®</sup> 310 Genetic Analyzer, GeneMapper <sup>®</sup> ID Software and POP-4 <sup>™</sup> Polymer .....	3
A. Instrument Preparation .....	3
B. Sample Preparation.....	4
C. Capillary Electrophoresis and Detection.....	4
D. Matrix Generation for the ABI PRISM <sup>®</sup> 310 Genetic Analyzer.....	5
4. Detection of Matrix Fragments Using the ABI PRISM <sup>®</sup> 310 Genetic Analyzer, GeneMapper <sup>®</sup> ID Software and POP-6 <sup>™</sup> Polymer .....	7
A. Instrument Preparation .....	7
B. Sample Preparation.....	9
C. Capillary Electrophoresis and Detection.....	9
D. Matrix Generation for the ABI PRISM <sup>®</sup> 310 Genetic Analyzer.....	10
5. Detection of Matrix Fragments Using the ABI PRISM <sup>®</sup> 310 Genetic Analyzer and GeneScan <sup>®</sup> Software .....	11
A. Instrument Preparation .....	11
B. Sample Preparation.....	12
C. Capillary Electrophoresis and Detection.....	12
D. Matrix Generation for the ABI PRISM <sup>®</sup> 310 Genetic Analyzer.....	12
6. Troubleshooting.....	14
7. Related Products .....	16

## 1. Description

Proper generation of a matrix file is critical to evaluate multicolor systems with the ABI PRISM® 310 Genetic Analyzer. To prepare a matrix, five standards are analyzed using the same capillary electrophoresis (CE) conditions that are used for samples and allelic ladders. The PowerPlex® 5-Dye Matrix Standards, 310<sup>(a-c)</sup>, consists of DNA fragments labeled with five different fluorescent dyes. One tube contains a DNA fragment labeled with fluorescein, one tube contains a DNA fragment labeled with JOE, one tube contains a DNA fragment labeled with TMR-ET, one tube contains a DNA fragment labeled with CXR-ET, and one tube contains a DNA fragment labeled with CC5. The PowerPlex® 5-Dye Matrix Standards, 310, was developed for use with the 5-dye PowerPlex® Systems.

Use the Fluorescein Matrix, JOE Matrix, TMR-ET Matrix, CXR-ET Matrix and CC5 Matrix for the blue, green, yellow, red and orange standards, respectively. A matrix should be generated for each individual instrument. Protocols to operate the fluorescence-detection instrumentation should be obtained from the manufacturer.

## 2. Product Components and Storage Conditions

Product	Size	Cat.#
PowerPlex® 5-Dye Matrix Standards, 310	50µl (each dye)	DG4600

Not for Medical Diagnostic Use. Includes:

- 50µl Fluorescein Matrix, 5-Dye (310)
- 50µl JOE Matrix, 5-Dye (310)
- 50µl TMR-ET Matrix, 5-Dye (310)
- 50µl CXR-ET Matrix, 5-Dye (310)
- 50µl CC5 Matrix, 5-Dye (310)

**Storage Conditions:** Store all components at -20°C in a nonfrost-free freezer. Do not store reagents in the freezer door, where the temperature can fluctuate. The fragments in the matrix standards are light-sensitive and must be stored in the dark. We strongly recommend that the matrix standards be stored with post-amplification reagents (away from pre-amplification materials) and used separately with different pipettes, tube racks, etc. We recommend using this product once, then discarding it.

Additional product information and ordering information for accessory components and related products are available upon request from Promega or at: [www.promega.com](http://www.promega.com)

### 3. Detection of Matrix Fragments Using the ABI PRISM® 310 Genetic Analyzer, GeneMapper® ID Software and POP-4™ Polymer

#### Materials to Be Supplied by the User

- 95°C dry heating block, water bath or thermal cycler
- crushed ice or ice-water bath
- 310 capillaries, 47cm × 50µm
- performance optimized polymer 4 (POP-4™ polymer; see note below)
- 10X genetic analyzer buffer
- sample tubes and septa
- aerosol-resistant pipette tips
- Hi-Di™ formamide (Applied Biosystems Cat.# 4311320)

**!** 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.

**Note:** When using the PowerPlex® ESI 16 and 17 Systems (Cat.# DC6770, DC6771, DC6780, DC6781), the use of performance optimized polymer 6 (POP-6™) might be necessary to resolve the 17.3 and 18 alleles and 18.3 and 19 alleles in the D12S391 allelic ladder and the 11.3 and 12 alleles in the D2S441 allelic ladder. See Section 4.

#### 3.A. Instrument Preparation

Refer to the *ABI PRISM® 310 Genetic Analyzer User's Manual* for instructions on cleaning the pump block, installing the capillary, calibrating the autosampler and adding polymer to the syringe.

1. Open the ABI PRISM® 310 Data Collection Software, Version 3.1.0.
2. To preheat the ABI PRISM® 310 Genetic Analyzer to 60°C, select "Manual Control" in the Window menu. In the Function menu, select "Temperature Set". Set Value to "60.0", then select "Execute". Close the Manual Control screen.
3. In the File menu, select "New" to open the Create New menu. Open a sample sheet (either "48-Tube" or "96-Tube").
4. In the upper right corner of the sample sheet, change "4 Dyes" to "5 Dyes". Enter the appropriate sample information in the Sample Name field. Matrix sample names should be descriptive; for example, add the color to the sample name. Label tubes with the corresponding sample names.
5. To save the sample sheet, select "Save As" in the File menu. Assign a name to the file, and save to the Sample Sheet folder. Close the file.

### 3.A. Instrument Preparation (continued)

6. In the File menu, select “New” to open the Create New menu.
7. Open the injection list.
8. Select the sample sheet (i.e., the .gss file) that was created in Step 5.
9. Use the **GS STR POP4 (1mL) G5v2.md5** module, which is available as a download from Applied Biosystems). Choose this module using the pull-down menu.

The settings should be:

Inj. Secs:	3
Inj. kV:	15.0
Run kV:	15.0
Run °C:	60
Run Time (minutes):	28

**Note:** The injection time may need to be increased or decreased, depending on instrument sensitivity. Peak heights of 1,000–3,000RFU are optimal for matrix generation.

10. Select “none” for the matrix file.

### 3.B. Sample Preparation

1. Thaw the matrix standards on ice. For each matrix standard, vortex the tube for 5–10 seconds to mix, then add 2µl of matrix standard to 25µl of Hi-Di™ formamide.
2. Denature each sample for 3 minutes at 95°C, and immediately chill on crushed ice or in an ice-water bath for 3 minutes. Denature samples just prior to loading.
3. Place tubes in the appropriate autosampler tray (48-tube or 96-well).
4. Place the autosampler tray in the instrument, and close the instrument doors.

### 3.C. Capillary Electrophoresis and Detection

1. After loading the sample tray and closing the doors, select “Run” to start the capillary electrophoresis system.
2. Monitor the electrophoresis by observing the raw data and status windows.

Each sample will take approximately 40 minutes for syringe pumping, sample injection and electrophoresis.

**Note:** The matrix files that are created will be .fsa files. After the run is finished, save or transfer the .fsa files to a secure location where they can be opened in a GeneMapper® project.

### 3.D. Matrix Generation for the ABI PRISM® 310 Genetic Analyzer

1. Open a new GeneMapper® project. To add matrix sample files to the new project, select "Add Samples to Project" in the File menu. Choose the appropriate run folder containing the .fsa files from Section 3.C. Highlight the run folder, and select "Add To List", then "Add".
2. To open the raw data for a specific matrix sample file, locate "Project" in the upper left corner of the screen, and double-click on the run folder to reveal the .fsa files.
3. Choose a single .fsa file to observe the raw data. While viewing the raw data, move the cursor to the region that is to the right of the primer peak and to the left of at least five peaks. Choose a region in a flat part of the baseline.
4. Record the data point value found at the lower left portion of the screen for use in Step 6. Repeat this step for each matrix standard.

<u>Dye Color</u>	<u>Corresponding Matrix</u>	<u>"Start At" Value</u>
Blue	Fluorescein Matrix	
Green	JOE Matrix	
Yellow	TMR-ET Matrix	
Red	CXR-ET Matrix	
Orange	CC5 Matrix	

5. To create a new matrix, select "GeneMapper Manager" in the Tools menu. Select the Matrices tab, then "New".
6. Define the new matrix in the Matrix Editor (Figure 1).

**Note:** The Matrix Name, "Start At" values and Matrix Result values shown in Figure 1 are instrument-specific and will change, depending on your instrument and whether you are using POP-4™ or POP-6™ polymer.

- a. Assign a matrix name in the Matrix Name field.
- b. Set Number of Dyes to "5".
- c. To select each matrix standard sample file, click on the dye color for each matrix (B for fluorescein, G for JOE, Y for TMR-ET, R for CXR-ET and O for CC5). Navigate to the .fsa sample file that corresponds to that dye, and double-click on it to add the sample file. Repeat this step for each matrix standard.

**Note:** To find the .fsa files in the default location, go to: "My Computer", "AB SW8DATA (D:)", "Applied Bio", "310", then "Runs", and locate the correct run folder.

- d. Enter the data point value recorded from Step 4 in the "Start at" field. Repeat this step for each matrix standard.

### 3.D. Matrix Generation for the ABI PRISM® 310 Genetic Analyzer (continued)

- e. Click on the Create button. The Matrix Result should give a value of 1.000 when comparing a dye to itself. Typically, all other values will be less than 1.000 except for the value indicated in red in Figure 1.

Select “OK”, and the matrix will be created in the Matrices tab of the GeneMapper Manager. Select “Done”.

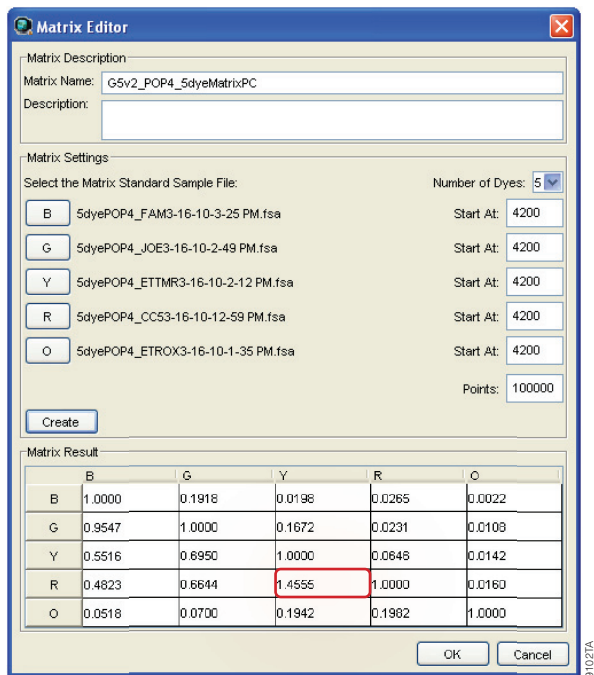




Figure 1. The Matrix Editor.

#### 4. Detection of Matrix Fragments Using the ABI PRISM® 310 Genetic Analyzer, GeneMapper® ID Software and POP-6™ Polymer

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

- 95°C dry heating block, water bath or thermal cycler
- crushed ice or ice-water bath
- 310 capillaries, 47cm × 50µm
- performance optimized polymer 6 (POP-6™ polymer; see note below)
- 10X genetic analyzer buffer
- sample tubes and septa
- aerosol-resistant pipette tips
- Hi-Di™ formamide (Applied Biosystems Cat.# 4311320)

 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.

**Note:** When using the PowerPlex® ESI 16 and 17 Systems (Cat.# DC6770, DC6771, DC6780, DC6781), the use of performance optimized polymer 6 (POP-6™) might be necessary to resolve the 17.3 and 18 alleles and 18.3 and 19 alleles in the D12S391 allelic ladder and the 11.3 and 12 alleles in the D2S441 allelic ladder.

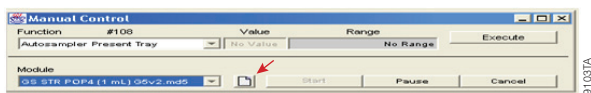
##### 4.A. Instrument Preparation

Refer to the *ABI PRISM® 310 Genetic Analyzer User's Manual* for instructions on cleaning the pump block, installing the capillary, calibrating the autosampler and adding polymer to the syringe.

1. Open the ABI PRISM® 310 Data Collection Software, Version 3.1.0.
2. To preheat the ABI PRISM® 310 Genetic Analyzer to 60°C, select "Manual Control" in the Window menu. In the Function menu, select "Temperature Set". Set Value to "60.0", then select "Execute".
3. It is necessary to create a module for use with POP-6 polymer the first time it is used. This module can be saved and used for subsequent runs. To make and save a module for use with POP-6™ polymer, choose the **GS STR POP4 (1mL) G5v2.md5** module using the pull-down menu under "Module".

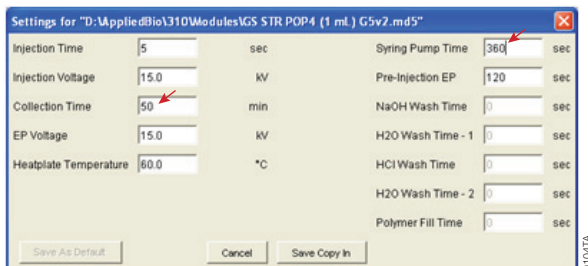
#### 4.A. Instrument Preparation (continued)

- Click on the folded page icon (Figure 2).

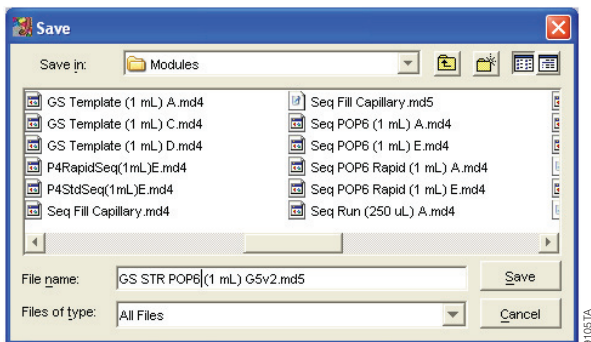


**Figure 2. The Manual Control screen.**

- Change the Collection Time to “50” and Syringe Pump Time to “360” (Figure 3). Select “Save Copy In”.
- Save the new module in the Modules folder. Change the file name to “GS STR POP6 (1mL) G5v2.md5”, and select “Save” (Figure 4).
- In the File menu, select “New” to open the Create New menu. Open a sample sheet (either “48-Tube” or “96-Tube”).
- In the upper right corner of the sample sheet, change “4 Dyes” to “5 Dyes”. Enter the appropriate sample information in the Sample Name field. Matrix sample names should be descriptive; for example, add the color to the sample name. Label tubes with the corresponding sample names.



**Figure 3. Changing the collection time and syringe pump time.**



**Figure 4. The Save screen.**

9. To save the sample sheet, select "Save As" in the File menu. Assign a name to the file, and save to the Sample Sheet folder. Close the file.
10. In the File menu, select "New" to open the Create New menu.
11. Open the injection list.
12. Select the sample sheet (i.e., the .gss file) that was created in Step 9.
13. Choose the **GS STR POP6 (1mL) G5v2.md5** module created in Step 6 using the pull-down menu.  
The settings should be:

Inj. Secs:	3
Inj. kV:	15.0
Run kV:	15.0
Run °C:	60
Run Time (minutes):	50
14. Select "none" for the matrix file.

#### 4.B. Sample Preparation

1. Thaw the matrix standards. For each matrix standard, vortex the tube for 5–10 seconds to mix, then add 2µl of matrix standard to 25µl of Hi-Di™ formamide.
2. Denature each sample for 3 minutes at 95°C, and immediately chill on crushed ice or in an ice-water bath for 3 minutes. Denature samples just prior to loading.
3. Place tubes in the appropriate autosampler tray (48-tube or 96-tube).
4. Place the autosampler tray in the instrument, and close the instrument doors.

#### 4.C. Capillary Electrophoresis and Detection

1. After loading the sample tray and closing the doors, select "Run" to start the capillary electrophoresis system.
2. Monitor the electrophoresis by observing the raw data and status windows.

Each sample will take approximately 60 minutes for syringe pumping, sample injection and electrophoresis.

**Note:** The matrix files that are created will be .fsa files. After the run is finished, save or transfer the .fsa files to a secure location where they can be opened in a GeneMapper® project.

#### 4.D. Matrix Generation for the ABI PRISM® 310 Genetic Analyzer

1. Open a new GeneMapper® project. To add matrix sample files to the new project, select "Add Samples to Project" in the File menu. Choose the appropriate run folder containing the .fsa files from Section 4.C. Highlight the run folder, and select "Add To List", then "Add".
2. To open the raw data for a specific matrix sample file, locate "Project" in the upper left corner of the screen, and double-click on the run folder to reveal the .fsa files.
3. Choose a single .fsa file to observe the raw data. While viewing the raw data, move the cursor to the region that is to the right of the primer peak and to the left of at least five peaks. Choose a region in a flat part of the baseline.
4. Record the data point value found at the lower left portion of the screen for use in Step 6. Repeat this step for each matrix standard.

Dye Color	Corresponding Matrix	"Start At" Value
Blue	Fluorescein Matrix	
Green	JOE Matrix	
Yellow	TMR-ET Matrix	
Red	CXR-ET Matrix	
Orange	CC5 Matrix	

5. To create a new matrix, select "GeneMapper Manager" in the Tools menu. Select the Matrices tab and "New".
6. Define the new matrix in the Matrix Editor (Figure 1).

**Note:** The Matrix Name, "Start At" values and Matrix Result values shown in Figure 1 are instrument-specific and will change, depending on your instrument and whether you are using POP-4™ or POP-6™ polymer.

- a. Assign a matrix name in the Matrix Name field.
- b. Set Number of Dyes to "5".
- c. To select each matrix standard sample file, click on the dye color for each matrix (B for fluorescein, G for JOE, Y for TMR-ET, R for CXR-ET and O for CC5). Navigate to the .fsa sample file that corresponds to that dye, and double-click on it to add the sample file. Repeat this step for each matrix standard.

**Note:** To find the .fsa files in the default location, go to: "My Computer", "AB SW8DATA (D:)", "Applied Bio", "310", then "Runs", and locate the correct run folder.

- d. Enter the data point value recorded from Step 4 in the "Start at" field. Repeat this step for each matrix standard.
- e. Click on the Create button. The Matrix Result should give a value of 1.000 when comparing a dye to itself. Typically, all other values will be less than 1.000 except for the value indicated in red in Figure 1.

Select "OK", and the matrix will be created in the Matrices tab of the GeneMapper Manager. Select "Done".

## 5. Detection of Matrix Fragments Using the ABI PRISM® 310 Genetic Analyzer and GeneScan® Software

### Materials to Be Supplied by the User

- 95°C dry heating block, water bath or thermal cycler
- crushed ice or ice-water bath
- 310 capillaries, 47cm × 50µm
- performance optimized polymer 4 (POP-4™ polymer)
- 10X genetic analyzer buffer
- sample tubes and septa
- aerosol-resistant pipette tips
- Hi-Di™ formamide (Applied Biosystems Cat.# 4311320)

**!** 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.

### 5.A. Instrument Preparation

Refer to the *ABI PRISM® 310 Genetic Analyzer User's Manual* for instructions on cleaning the pump block, installing the capillary, calibrating the autosampler and adding polymer to the syringe.

1. Open the ABI PRISM® 310 Data Collection Software.

**Note:** The software must be compatible with 5-dye chemistry (i.e., it must be version 2.1 for Macintosh® operating systems.)

2. Prepare a GeneScan® sample sheet for five dyes as described in the *ABI PRISM® 310 Genetic Analyzer User's Manual*. Enter the appropriate sample information in the Sample Info column. Create a new GeneScan® injection list. Select the appropriate sample sheet using the pull-down menu.

3. The recommended module is **GS STR POP4 (1mL) G5v2**; choose this module using the pull-down menu. If this module is not available, choose GS STR POP4 (1mL) G5. The settings should be:

Inj. Secs:	3
Inj. kV:	15.0
Run kV:	15.0
Run °C:	60
Run Time (minutes):	28

**Note:** The injection time may need to be increased or decreased, depending on instrument sensitivity. Peak heights of 1,000-3,000RFU are optimal for matrix generation.

4. Select "none" for the matrix file.

## 5.B. Sample Preparation

1. Thaw the matrix standards. For each matrix standard, vortex the tube for 5–10 seconds to mix, then add 2µl of matrix standard to 25µl of Hi-Di™ formamide.
2. Denature each sample for 3 minutes at 95°C, and immediately chill on crushed ice or in an ice-water bath for 3 minutes. Denature samples just prior to loading.
3. Place tubes in the appropriate autosampler tray (48-tube or 96-tube).
4. Place the autosampler tray in the instrument, and close the instrument doors.

## 5.C. Capillary Electrophoresis and Detection

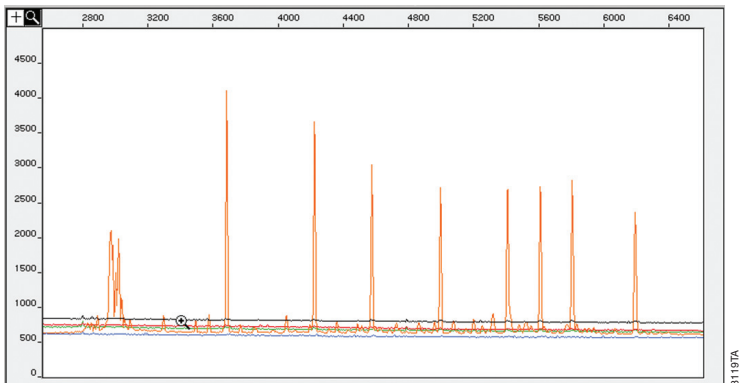
1. After loading the sample tray and closing the doors, select “Run” to start the capillary electrophoresis system.
2. Monitor the electrophoresis by observing the raw data and status windows.  
  
Each sample will take approximately 40 minutes for syringe pumping, sample injection and electrophoresis.

## 5.D. Matrix Generation for the ABI PRISM® 310 Genetic Analyzer

1. Open the GeneScan® project.
2. Review the raw data from the individual matrix samples. Highlight the sample file name, then go to the Sample menu and select “raw data”. Move the cursor to the right of the primer peak and to the left of at least 5 peaks, so the crosshair is on a flat portion of the baseline. Record the X value shown at the bottom of the window. Select an area for matrix generation. For optimal results, use as many peaks as possible. See Figure 5.
3. In the File menu, select “New,” then click on the Matrix icon. Set Number of Dyes to “5”. Click on the dye color for each matrix, and indicate the sample file that corresponds to that dye. Enter the X value recorded from Step 2 in the “Start at” field.

<u>Dye Color</u>	<u>Corresponding Matrix</u>	<u>“Start At” Value</u>
Blue	Fluorescein Matrix	
Green	JOE Matrix	
Yellow	TMR-ET Matrix	
Red	CXR-ET Matrix	
Orange	CC5 Matrix	

4. Select “OK”, and the matrix file will be generated.



**Figure 5. CC5 Matrix raw data.** The CC5 Matrix standard was analyzed using an ABI PRISM® 310 Genetic Analyzer. GeneScan® analysis software was used to view the raw data (in the Sample menu). The cursor was placed on the baseline, and the “Start at” value of 3407 was determined using the readout in the lower left corner of the window (not shown).

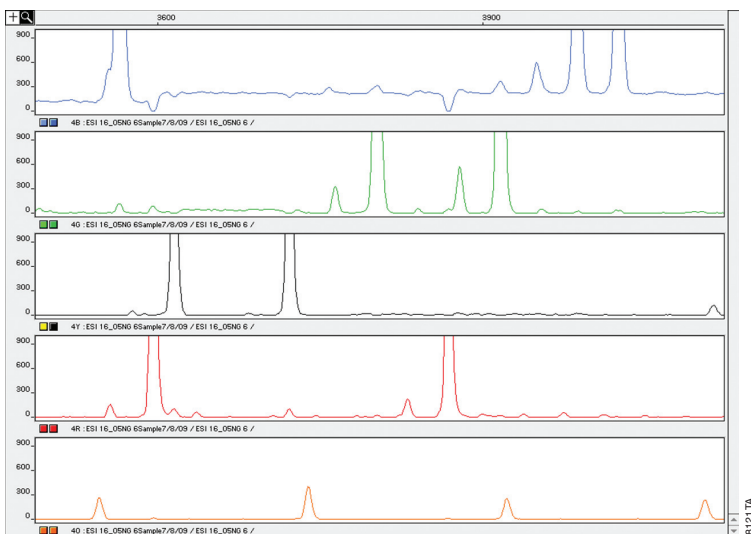
5. Save the matrix file in the Matrix Standards folder located in the GeneScan® folder. For the Macintosh® version of the software, a copy of the matrix file is automatically saved in the GS Matrix folder. For the Windows NT® version of the software, store a copy of the matrix file in the Matrix folder at: C:\appliedbio\shared\analysis\sizecaller\matrix.
6. A new matrix can be applied to previously run samples by highlighting the sample in the GeneScan® project. In the Sample menu, select “Install new matrix”, highlight the new matrix and select “Open”. The new matrix will be applied to the sample file, and the samples can be analyzed using the new matrix.

## 6. Troubleshooting

For questions not addressed here, please contact your local Promega Branch Office or Distributor. Contact information available at: [www.promega.com](http://www.promega.com). E-mail: [techserv@promega.com](mailto:techserv@promega.com)

<b>Symptoms</b>	<b>Causes and Comments</b>
Unable to generate a matrix due to faint or no peaks	Peak heights were too low. Peak heights should be 1,000-3,000RFU for the ABI PRISM® 310 Genetic Analyzer. To increase peak heights, increase the injection time or loading volume.
	Poor capillary electrophoresis (CE) injection. Re-inject the sample. Check the syringe for leakage. Check the laser power.
	Poor-quality formamide was used. Use only fresh Hi-Di™ formamide when running samples on the ABI PRISM® 310 Genetic Analyzer.
	Samples were degraded due to improper storage. Store matrix standards at -20°C protected from light. Do not store in the freezer door or in a frost-free freezer.
	Samples were not denatured. Heat-denature samples, and immediately chill on crushed ice or in an ice-water bath before loading the gel or capillary. Denature samples just prior to loading.
Peak heights too high	Peak heights are above 3,000RFU. To decrease peak heights, decrease the injection time or loading volume.
Poor-quality matrix (extra peaks visible in one or all of the color channels)	CE-related artifacts (“spikes”). Minor voltage changes or urea crystals passing by the laser can cause “spikes” or unexpected peaks. Spikes sometimes appear in one color but often are easily identified by their presence in more than one color. Re-inject the samples to confirm.
	CE-related artifacts (contaminants). Contaminants in the water used with the ABI PRISM® 310 Genetic Analyzer and for diluting the 10X genetic analyzer buffer can generate peaks in the fluorescein and JOE dye channels. Use autoclaved water to clean the pump block and prepare sample dilutions. Change vials, and wash the buffer reservoir.
Extra peaks in the CC5 and CXR-ET dye channels	Extra peaks may appear in the CC5 and CXR-ET dye channels but do not affect matrix quality.

Symptoms	Causes and Comments
Poor-quality matrix [elevated baseline (Figure 6) and/or inverted peaks in analyzed samples]	<p>Matrix used was generated on another instrument. A matrix must be generated for each instrument.</p> <p>Wrong dye was used. Generate the matrix using the same dyes as those in the samples.</p> <p>Oversubtraction of signal occurred because signal was saturated. When generating a matrix, avoid choosing samples with peak heights that are higher than the recommended RFU values, as this can result in a matrix that causes inverted peaks or elevated baseline. Analyzed sample results may be improved by diluting matrix samples in water before preparing them for use. Alternatively, decrease the injection time.</p>
Inverted peaks in matrix baseline	<p>Incorrect or no "Start At" value was entered. The "Start At" value entered in Sections 3.D, 4.D and 5.D should be chosen in a region with a flat baseline.</p>



**Figure 6. Elevated baseline.** A sample was analyzed using an ABI PRISM® 310 Genetic Analyzer and GeneScan® analysis software. The resulting electropherogram shows an elevated baseline in the blue channel (top panel).

## 6. Troubleshooting (continued)

Symptoms	Causes and Comments
Previously generated matrix no longer performs optimally	Changes to or aging of instrument components. Instrument sensitivity can change if the instrument is moved or recently serviced (replacement or realignment of the laser, CCD camera, power supply or mirrors). The sensitivity also can change over time due to aging of the instrument. These changes can result in poor matrix performance. Generate a new matrix.

## 7. Related Products

Product	Size	Cat.#
PowerPlex® ESX 16 System	100 reactions	DC6711
	400 reactions	DC6710
PowerPlex® ESX 17 System	100 reactions	DC6721
	400 reactions	DC6720
PowerPlex® ESI 16 System	100 reactions	DC6771
	400 reactions	DC6770
PowerPlex® ESI 17 System	100 reactions	DC6781
	400 reactions	DC6780

Not for Medical Diagnostic Use.

### Accessory Components

Product	Size	Cat.#
Nuclease-Free Water*	50ml	P1193
Water, Amplification Grade*	6,250µl (5 × 1,250µl)	DW0991
PowerPlex® 5-Dye Matrix Standards, 3100/3130**	25µl (each dye)	DG4700

\*For Laboratory Use.

\*\*Not for Medical Diagnostic Use.

<sup>(4)</sup>This product or portions thereof is manufactured and sold under license from GE Healthcare under Australia Pat. No. 692230, Austria Pat. No. E236994, Belgium Pat. No. 0743987, Canada Pat. No. 2231475, EP Pat. Nos. 0743987 and 0851867, France Pat. Nos. 0743987 and 0851867, Germany Pat. Nos. 19581489, 69530286.8 and 0851867, Italy Pat. Nos. 0743987 and 0851867, Japan Pat. No. 3066984, Liechtenstein Pat. Nos. 0743987 and 0851867, Netherlands Pat. Nos. 0743987 and 0851867, Spain Pat. Nos. 2197193 and 2173310, Sweden Pat. Nos. 0743987 and 0851867, Switzerland Pat. Nos. 0743987 and 0851867, United Kingdom Pat. Nos. 0743987 and 0851867, U.S. Pat. Nos. 5,654,419, 5,688,648, 5,869,255, 6,177,247, 5,707,804, 6,028,190, 6,544,744, 7,015,000 and 5,728,528 and other pending and foreign patent applications.

<sup>(5)</sup>**End User Terms and Conditions**

**Acceptance.** These terms and conditions shall govern the purchase, use, transfer and acceptance of the products described in the purchase order, quotation or invoice, which products are sold and distributed by Promega to the buyer/transferee of such products (the "End User"). The transfer/sale of products to the End User is expressly conditional upon End User's acceptance of these terms and conditions.

**Restrictions on Use.** End Users are specifically not authorized to and are forbidden from reselling, transferring or distributing any products either as a stand alone product or as a component of another product. The right to use the products does not, in and of itself, include or carry any right of the End User to any GE Healthcare Bio-Sciences Corp.'s technology or intellectual property other than expressly provided herein. End Users may not use sequence(s) in an attempt to reverse engineer parameters of any of GE Healthcare Bio-Sciences Corp. proprietary products or services.

**Disclaimer of Warranties.** GE Healthcare Bio-sciences Corp. provides no warranties to end user (statutory or implied), including without limitation, as to product quality, condition, description, merchantability or fitness for a particular purpose, and all such warranties are hereby expressly disclaimed. GE Healthcare Bio-Sciences Corp. hereby expressly disclaims any warranty regarding results obtained through the use of the products, including without limitation any claim of inaccurate, invalid or incomplete results.

**Exclusion of Liability.** GE Healthcare Bio-Sciences Corp. and its affiliates shall have no liability to an End User, including, without limitation, for any loss of use or profits, business interruption or any consequential, incidental, special or other indirect damages of any kind, regardless of how caused and regardless of whether an action in contract, tort, strict product liability or otherwise.

<sup>(6)</sup>TMR-ET, CXR-ET and CC5 dyes are proprietary.

© 2009, 2011 Promega Corporation. All Rights Reserved.

PowerPlex is a registered trademark of Promega Corporation.

ABI PRISM, GeneMapper and GeneScan are registered trademarks of Applied Biosystems Corporation. Hi-Di, POP-4 and POP-6 are trademarks of Applied Biosystems Corporation. Macintosh is a registered trademark of Apple Computer, Inc. Windows NT is a registered trademark of Microsoft Corporation.

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

All prices and specifications are subject to change without prior notice.

Product claims are subject to change. Please contact Promega Technical Services or access the Promega online catalog for the most up-to-date information on Promega products.