



PowerPlex® Matrix Standards for Use on the Spectrum Compact CE System Technical Manual

Instructions for Use of Products DG4800, DG4850, DG4900 and DG5010



Table of Contents

1	Description.....	4
2	PowerPlex® 8C Spectral Calibration	5
	2.1 Product Components and Storage Conditions	5
	2.2 Matrix Sample Preparation	6
	2.3 Assembling Sample Cartridge.....	7
	2.4 Instrument Preparation and Spectral Calibration	8
	2.5 Results	13
3	PowerPlex® 6C Spectral Calibration	16
	3.1 Product Components and Storage Conditions	16
	3.2 Matrix Sample Preparation	17
	3.3 Assembling Sample Cartridge.....	18
	3.4 Instrument Preparation and Spectral Calibration	19
	3.5 Results	24
4	PowerPlex® 5C Spectral Calibration	27
	4.1 Product Components and Storage Conditions	27
	4.2 Matrix Sample Preparation	28
	4.3 Assembling Sample Cartridge.....	29
	4.4 Instrument Preparation and Spectral Calibration	30
	4.5 Results	35

5	PowerPlex® 4C Spectral Calibration	38
5.1	Product Components and Storage Conditions	38
5.2	Matrix Sample Preparation	39
5.3	Assembling Sample Cartridge.....	40
5.4	Instrument Preparation and Spectral Calibration	41
5.5	Results	46
6	Troubleshooting.....	49
7	Summary of Changes.....	51

All technical literature is available at: www.promega.com/protocols/

Visit the web site to verify that you are using the most current version of this Technical Manual.

E-mail Promega Technical Services if you have questions on use of this system: genetic@promega.com

Description

Proper generation of spatial and spectral calibration files is critical to evaluate multicolor systems with the Spectrum Compact CE System. Refer to the *Spectrum Compact CE System Operating Manual #TMD058* for the instrument maintenance schedule and instructions for installation of the capillary array, buffers and polymer cartridge and performing spatial calibration.

The PowerPlex® Matrix Standards (see Table 1) consist of DNA fragments labeled with different fluorescent dyes in one tube. Once generated, the spectral calibration file is applied during sample detection to calculate the spectral overlap and separate the raw fluorescent signals into individual color signals. A spectral calibration must be generated for each individual instrument, and must be performed after the installation of a capillary array following a spatial calibration. A spectral calibration should also be performed after any major maintenance on the system, such as changing the laser, replacing the camera or if a decrease in spectral separation is observed in the STR results. We also recommend generating a new matrix any time the instrument is moved to a new location.

Note: The instructions contained in this manual can be used with the HITACHI DS3000 Compact CE Sequencer.

Table 1. PowerPlex® Matrix Standards

Matrix Standard	Dyes	Cat.#	STR Systems
PowerPlex® 8C Matrix Standard ^(a,b)	FL-8C, JOE-8C, AQA-8C, TMR-8C, CXR-8C, TOM-8C, WEN-8C and CCO-8C	DG5010	Powerplex® 35GY PowerPlex® 18E
PowerPlex® 6C Matrix Standard ^(a,c)	FL-6C, JOE-6C, TMR-6C, CXR-6C, TOM-6C and WEN	DG4900	PowerPlex® Fusion 6C
PowerPlex® 5C Matrix Standard ^(a,d)	Fluorescein, JOE, TMR-ET, CXR-ET and WEN	DG4850	PowerPlex® Fusion PowerPlex® ESX/ESI 17 Fast PowerPlex® ESX/ESI 16 Fast PowerPlex® ESX 17 PowerPlex® ESI 17 Pro PowerPlex® 21 PowerPlex® 18D PowerPlex® Y23
PowerPlex® 4C Matrix Standard	Fluorescein, JOE, TMR and CXR	DG4800	PowerPlex® 16 HS PowerPlex® CS7

PowerPlex® 8C Spectral Calibration

2.1 Product Components and Storage Conditions

PRODUCT	SIZE	CAT.#
PowerPlex® 8C Matrix Standard	5 preps	DG5010

Not For Medical Diagnostic Use. Includes:

- 150µl 8C Matrix Mix
- 5 × 200µl Matrix Dilution Buffer

Storage Conditions

Upon receipt, store all components at +2°C to +10°C, protected from light. We strongly recommend that the PowerPlex® 8C Matrix Standard be stored with the post-amplification reagents. The PowerPlex® 8C Matrix Standard is light-sensitive; dilute the 8C Matrix Mix in the Matrix Dilution Buffer in the provided amber tube. Store the diluted 8C Matrix Mix at 2–10°C for up to 1 week.



Do not freeze the PowerPlex® 8C Matrix Standard components.

Materials to be Supplied by the User

- centrifuge compatible with 8-tube strips
- aerosol-resistant pipette tips
- Spectrum Compact Capillary Cartridge, 4-Capillary 36cm (Cat.# CE2340)
- Spectrum Compact Polymer4 (Cat.# CE2304, CE2404)
- Spectrum Compact Buffer (Cat.# CE2300)
- Spectrum Compact Cathode Septa Mat (Cat.# CE2301)
- Spectrum Compact Cathode Retainer (Cat.# CE2302)
- Spectrum Compact Strip Base & Retainer, 32-Well (Cat.# CE2332)
- MicroAmp® Optical 8-Tube Strip, 0.2ml (Applied Biosystems, Cat.# 4316567)
- Strip Septa Mat, 8-Well (Cat.# CE2308)
- Hi-Di™ formamide (Applied Biosystems, Cat.# 4311320)

For additional information on performing spectral calibration, refer to the *Spectrum Compact CE System Operating Manual #TMD058*.



The quality of formamide is critical. Use Hi-Di™ formamide. Freeze the formamide in aliquots at –20°C. Multiple freeze-thaw cycles or long-term storage at 4°C can cause breakdown of formamide. Poor-quality formamide can contain ions that compete with DNA during injection, which results in lower peak heights.



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.

Notes:

- a. Only use MicroAmp® Optical 8-Tube Strips, 0.2ml (Applied Biosystems, Cat.# 4316567) as a source of 8-well strip tubes. Use of other 8-well strip tubes may affect performance or damage the Spectrum Compact CE System.
- b. Wear gloves when handling consumables and sample cartridges.

2.2 Matrix Sample Preparation

1. Vortex the 8C Matrix Mix for 10–15 seconds prior to use. Add 10µl of the 8C Matrix Mix to one tube of the Matrix Dilution Buffer. Vortex for 10–15 seconds. Label the tube with the date of dilution. The diluted 8C Matrix Mix can be stored for up to 1 week at 2–10°C.
2. Vortex the diluted 8C Matrix Mix prepared in Step 1 for 10–15 seconds, then add 10µl to 500µl of Hi-Di™ formamide.
3. Vortex the diluted 8C Matrix Mix with formamide prepared in Step 2 for 10–15 seconds, then add 15µl to each of the first four wells of an 8-well strip tube. After placing the septa on the 8-well strip tube, briefly centrifuge the plate to remove bubbles. Do not heat denature.

		Wells							
		1	2	3	4	5	6	7	8
Lane	A	Matrix Mix							
	B								
	C								
	D								

2.3 Assembling Sample Cartridge

1. Place the 8-well strip tube into the strip base in Lane A with the samples in positions 1–4 (Figure 1).

Note: Lane names A to D and well numbers 1 to 8 are embossed on the strip base. Be sure to check the lane name and well numbers when placing the 8-well strip tube into the base.

2. To complete the assembly, place the retainer over the strip in the strip base, aligning the lane names A to D and well numbers 1 to 8 on the retainer to those on the strip base and pressing until the retainer clicks into the strip base (Figure 2).

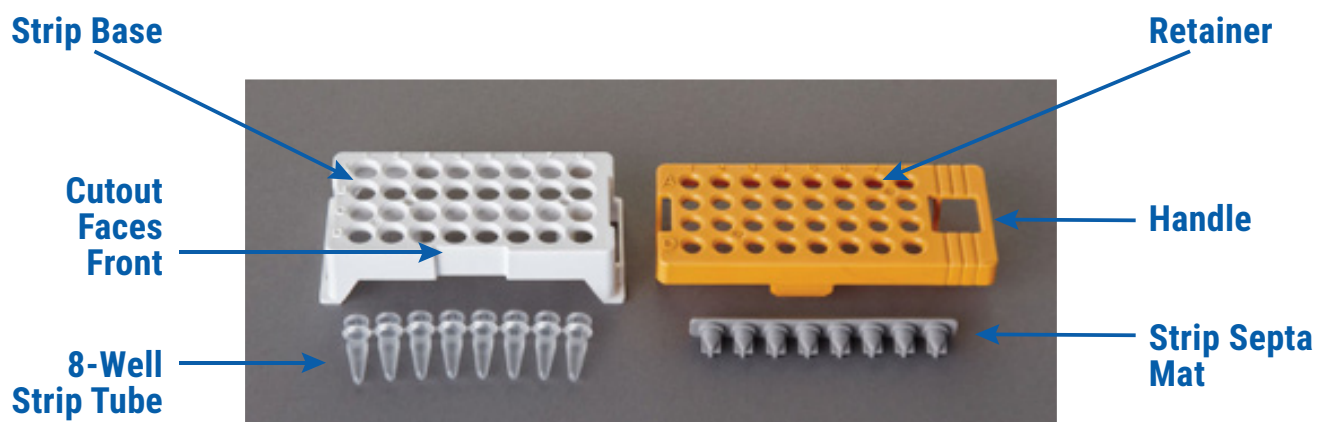


Figure 1. Assembling the Spectrum Compact Strip Base and Retainer.

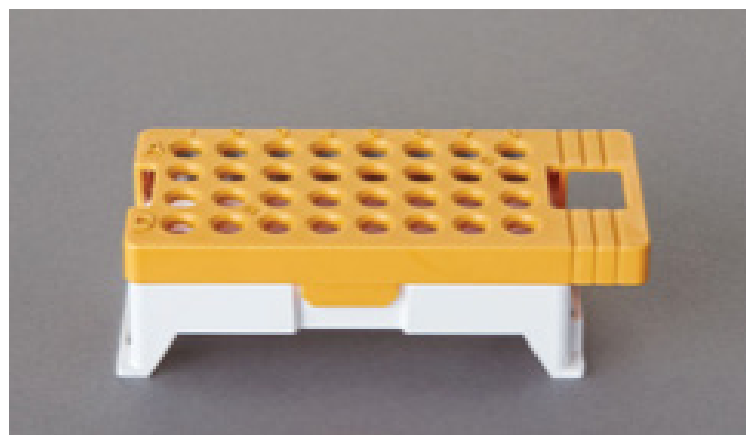


Figure 2. Assembled Spectrum Compact Sample Cartridge.

2.4 Instrument Preparation and Spectral Calibration

These instructions are intended as a guide for running PowerPlex® Matrix Standards on the Spectrum Compact CE System. They are not intended as comprehensive instructions for using the Spectrum Compact CE System. Refer to the *Spectrum Compact CE System Operating Manual #TMD058* for more details on performing spectral calibration.

Notes:

- We have found that the use of fresh polymer and new capillary array results in an optimal spectral calibration.
 - We do not recommend performing spectral calibration with expired reagents. Expired reagents should be replaced before performing a spectral calibration.
 - Refer to the *Spectrum Compact CE System Operating Manual #TMD058* for more details on installation of consumables, instrument maintenance and spatial calibration.
- Select the **Consumables** icon in the Header on the 'Main Menu' screen (Figure 3). Ensure that the consumables are not expired and that adequate injections remain for consumables installed.

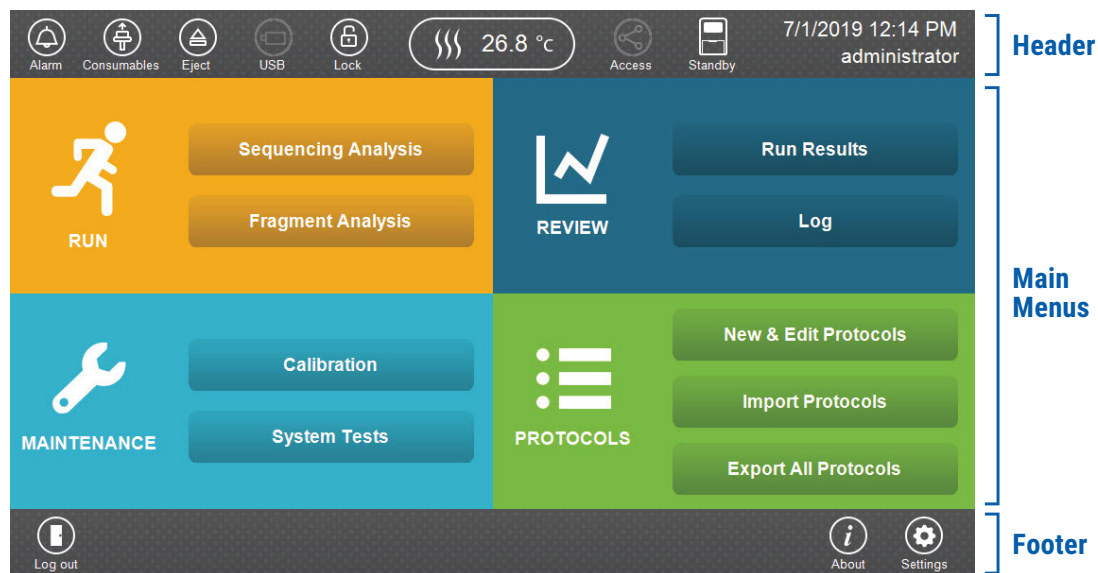


Figure 3. Spectrum Compact CE System Software 'Main Menu' screen.

2. Select the **Oven Temperature** icon in the Header on the 'Main Menu' screen, as shown in Figure 4, to start preheating the oven temperature to 60°C. The temperature displayed will change as the temperature of the oven increases. When 60°C is reached, a check mark will appear adjacent to the temperature.

Note: We recommend you preheat the oven for at least 30 minutes prior to starting a run. The oven will automatically turn off after 2 hours if a run is not started.

Oven Temperature Icon

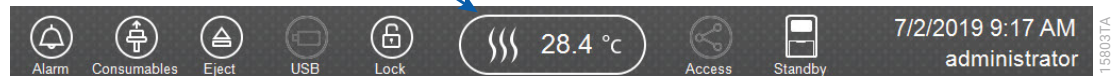


Figure 4. Preheating Oven.

3. Select **Calibration** on the maintenance portion of the 'Main Menu' screen (Figure 3) then select **Spectral Calibration** on the 'Maintenance Calibration' screen (Figure 5).

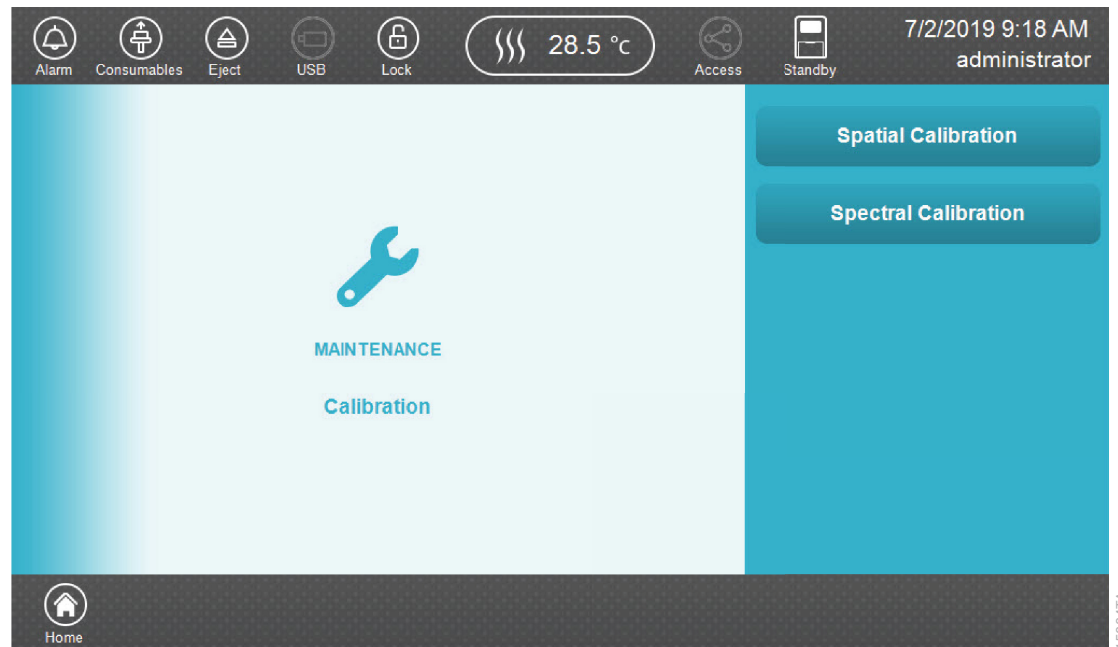


Figure 5. 'Maintenance Calibration' screen.

4. Use the scroll arrows on the right-hand side of the 'Dye Set List' screen (Figure 6) to find the correct Dye Set/Application Type/Polymer combination from the displayed list. To perform a spectral calibration using the PowerPlex® 8C Matrix Standard on Polymer4, select **Promega 8-dye** with "Fragment" and "Polymer4" as application and polymer types, respectively, then select **Calibration**. The 'Assemble the Cartridge' screen will open (Figure 7).

No	Calibrated Date / Dye Set	Application	Polymer	Capillary
001	08/09/2023 13:05:09 Promega 8-dye	Fragment	Polymer4	0000557870P323B0102
002	08/07/2023 12:43:33 Promega 8-dye	Fragment	Polymer7	0000557870P323B0102
003	06/28/2023 16:06:25 Promega 5-dye	Fragment	Polymer4	0000557870P323B0102
004	05/09/2023 12:44:50 T 4-dye sequencing	Sequencing	Polymer7	0000543661P322I0411
005	03/02/2023 10:17:32 Promega 4-dye sequencing	Sequencing	Polymer7	0000546543P322J0001

Figure 6. 'Dye Set List' screen.

5. Select **Next** on the 'Assemble the Cartridge' screen (Figure 7). A message window will open indicating that the autosampler is moving and telling the user to not open the door. In addition, the status indicator flashes green while the autosampler is moving. After autosampler movement is complete, the message window closes and the status indicator returns to a steady green.

Note: Do not open the front door of the Spectrum Compact CE System while the autosampler is in motion.

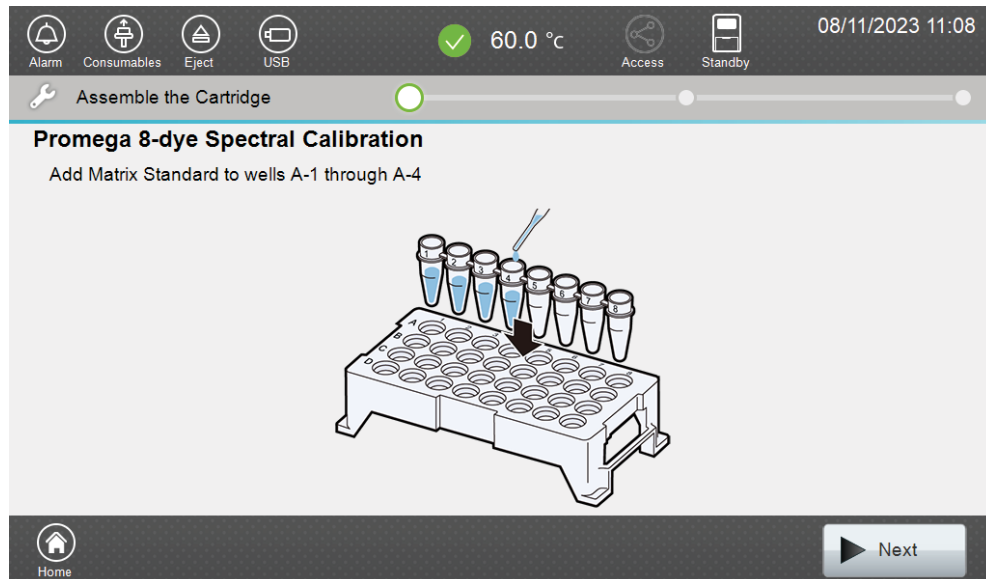


Figure 7. 'Assemble the Cartridge' screen.

6. Open the front door of the Spectrum Compact CE System and mount the sample cartridge on the autosampler by following the instructions displayed on the 'Install the Cartridge' screen (Figure 8).

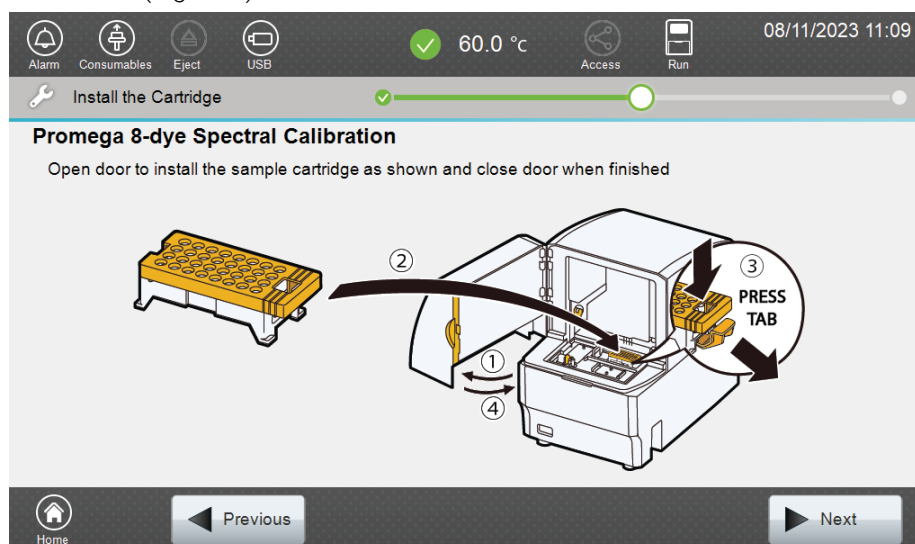


Figure 8. 'Install the Cartridge' screen.

- After mounting the sample cartridge on the autosampler, close the front door of the Spectrum Compact CE System and wait for the status indicator to stop flashing amber and turn steady green.

Note: Do not open the front door of the Spectrum Compact CE System while the autosampler is in motion.

- After the autosampler has returned to its home position, the 'Spectral Calibration' screen will automatically be displayed (Figure 9). Select **Run** to start the spectral calibration.

Note: The 'Raw Data' tab can be used to monitor the run.

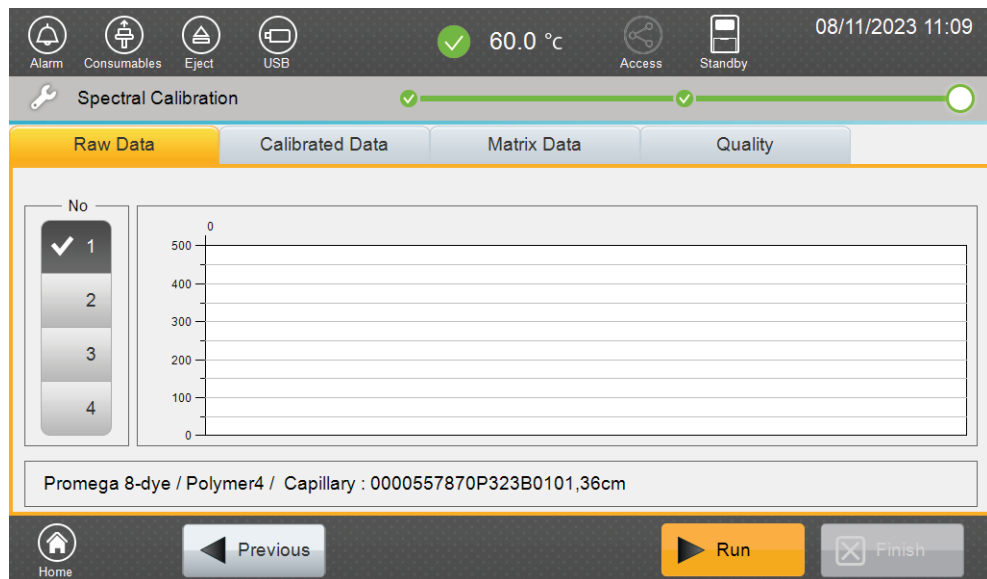


Figure 9. 'Spectral Calibration' screen.

2.5 Results

- Following the run, the 'Raw Data' tab (Figure 10) will be displayed. The minimum peak height for spectral calibration is 500 relative fluorescent units (RFU) and the maximum peak height is 32,767RFU.

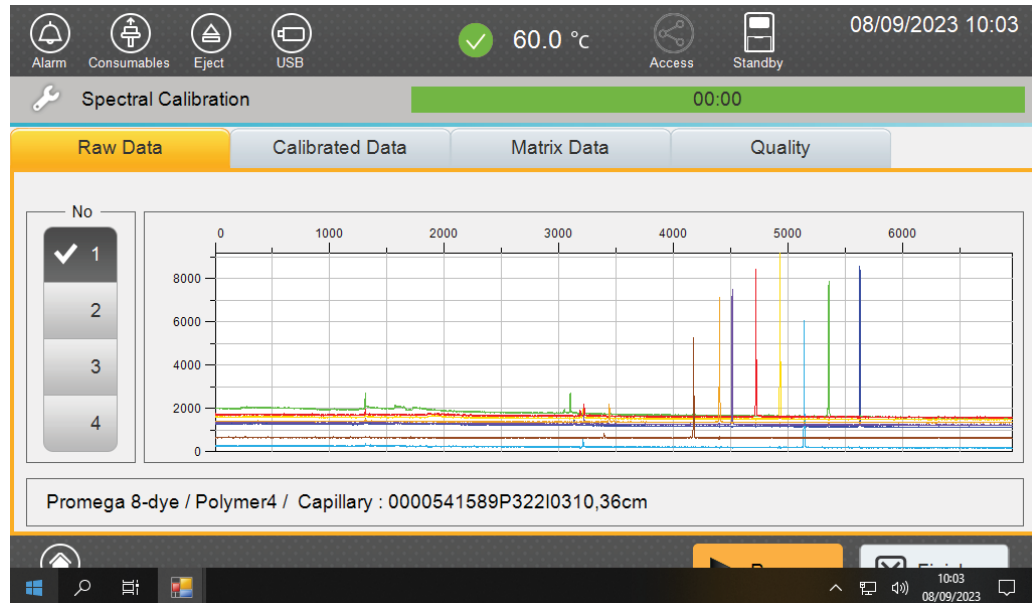


Figure 10. Spectral Calibration 'Raw Data' tab screen.

- The 'Calibrated Data' tab can be used to view the matrix peaks with both baseline and spectral applied for each capillary (Figure 11).

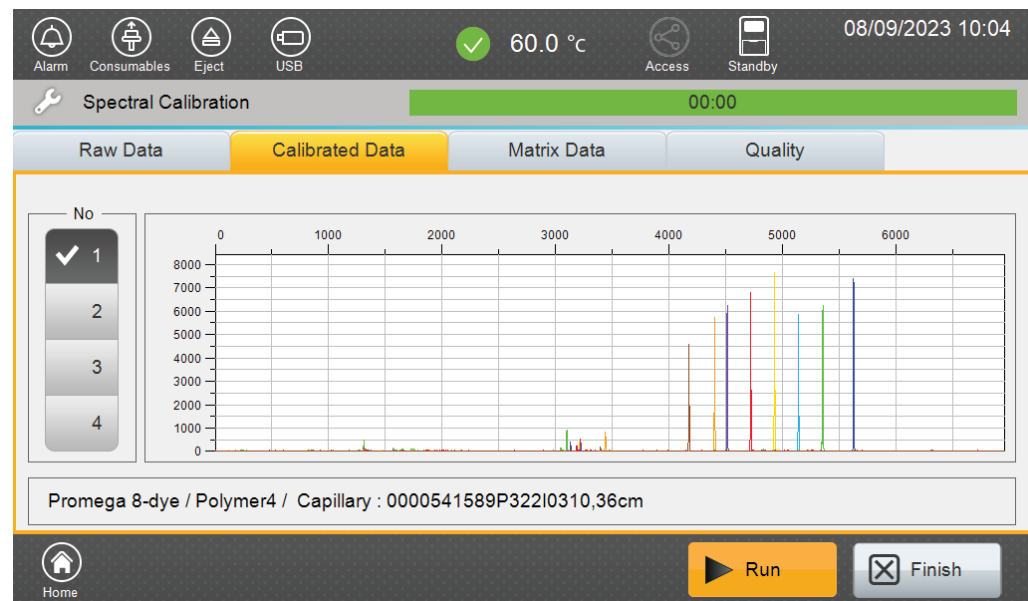


Figure 11. Spectral Calibration 'Calibrated Data' tab screen.

- The 'Matrix Data' tab can be used to view emission spectra for each capillary (Figure 12).

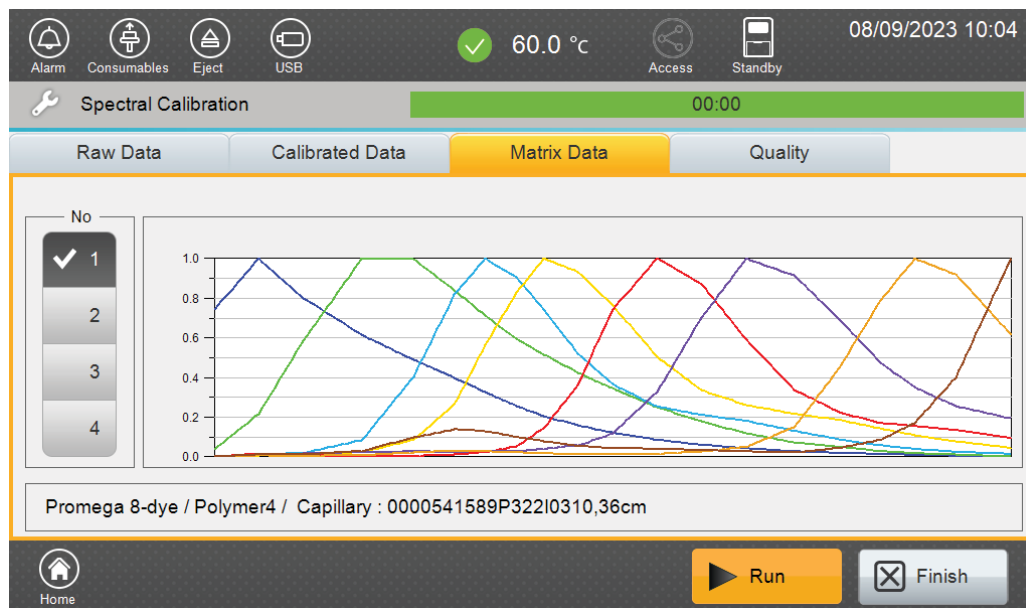


Figure 12. Spectral Calibration 'Matrix Data' tab screen.

- Review the quality of the spectral calibration by selecting the 'Quality' tab (Figure 13).

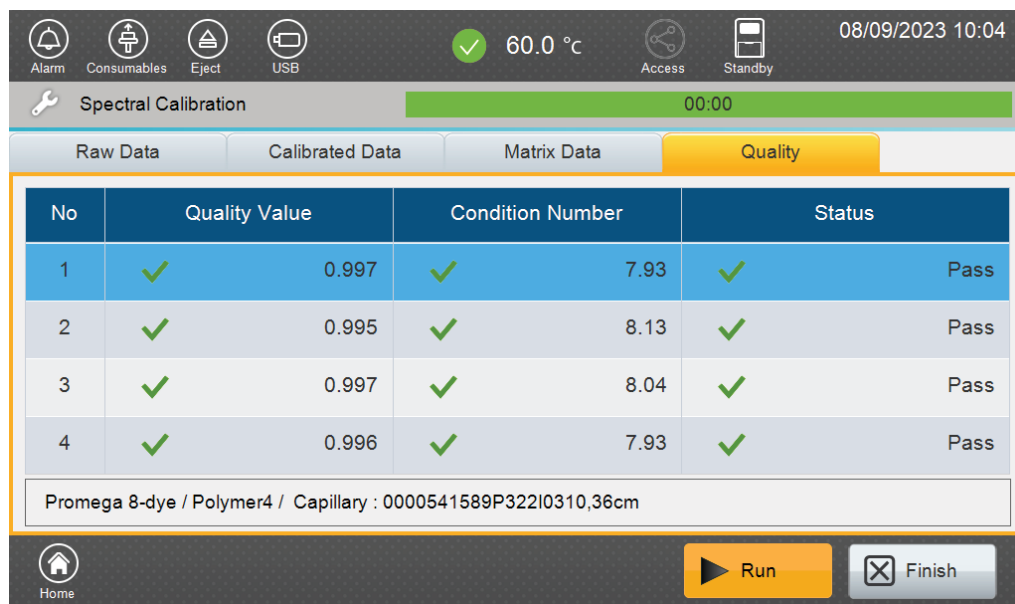


Figure 13. Spectral Calibration 'Quality' tab screen.

- Each capillary must meet the passing criteria of ≥ 0.95 for the Quality Value and < 8.5 for the Condition Number.

6. If one capillary fails to meet the criteria, it is possible to borrow spectral data from an adjacent capillary. Refer to the *Spectrum Compact CE System Operating Manual #TMD058* for details. If more than one capillary fails, the spectral must be rerun.

Notes:

- a. Selecting **Run** will rerun the spectral calibration.
 - b. Refer to Section 5 for troubleshooting if more than one capillary fails to meet the criteria.
7. After reviewing the results, select **Finish**. This will open a confirmation window. Select **Yes** to apply the spectral calibration results (Figure 14). The spectral calibration result will not be saved unless you select **Yes** on this window.

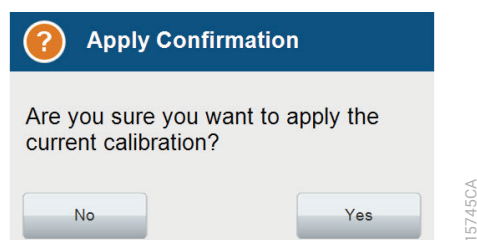


Figure 14. Apply spectral calibration confirmation window.

PowerPlex® 6C Spectral Calibration

3.1 Product Components and Storage Conditions

PRODUCT	SIZE	CAT.#
PowerPlex® 6C Matrix Standard	5 preps	DG4900

Not For Medical Diagnostic Use. Includes:

- 150µl 6C Matrix Mix
- 5 × 200µl Matrix Dilution Buffer

Storage Conditions

Upon receipt, store all components at -30°C to -10°C in a nonfrost-free freezer, protected from light. Do not store reagents in the freezer door, where the temperature can fluctuate. After the first use, store the PowerPlex® 6C Matrix Standard components at $2-10^{\circ}\text{C}$, protected from light. We strongly recommend that the PowerPlex® 6C Matrix Standard be stored with post-amplification reagents. The PowerPlex® 6C Matrix Standard is light-sensitive; dilute the 6C Matrix Mix in the Matrix Dilution Buffer in the provided amber tube. Store the diluted 6C Matrix Mix at $2-10^{\circ}\text{C}$ for up to 1 week.



Do not refreeze the PowerPlex® 6C Matrix Standard components.

Materials to be Supplied by the User

- centrifuge compatible with 8-tube strips
- aerosol-resistant pipette tips
- Spectrum Compact Capillary Cartridge, 4-Capillary 36cm (Cat.# CE2340)
- Spectrum Compact Polymer4 (Cat.# CE2304, CE2404)
- Spectrum Compact Buffer (Cat.# CE2300)
- Spectrum Compact Cathode Septa Mat (Cat.# CE2301)
- Spectrum Compact Cathode Retainer (Cat.# CE2302)
- Spectrum Compact Strip Base & Retainer, 32-Well (Cat.# CE2332)
- MicroAmp® Optical 8-Tube Strip, 0.2ml (Applied Biosystems, Cat.# 4316567)
- Strip Septa Mat, 8-Well (Cat.# CE2308)
- Hi-Di™ formamide (Applied Biosystems, Cat.# 4311320)

For additional information on performing spectral calibration, refer to the *Spectrum Compact CE System Operating Manual #TMD058*.



The quality of formamide is critical. Use Hi-Di™ formamide. Freeze the formamide in aliquots at –20°C. Multiple freeze-thaw cycles or long-term storage at 4°C can cause breakdown of formamide. Poor-quality formamide can contain ions that compete with DNA during injection, which results in lower peak heights.



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.

Notes:

- a. Only use MicroAmp® Optical 8-Tube Strips, 0.2ml (Applied Biosystems, Cat.# 4316567) as a source of 8-well strip tubes. Use of other 8-well strip tubes may affect performance or damage the Spectrum Compact CE System.
- b. Wear gloves when handling consumables and sample cartridges.

3.2 Matrix Sample Preparation

1. At the first use, thaw the 6C Matrix Mix and Matrix Dilution Buffer completely. After the first use, store the reagents at 2–10°C, protected from light.
2. Vortex the 6C Matrix Mix for 10–15 seconds prior to use. Add 10µl of the 6C Matrix Mix to one tube of the Matrix Dilution Buffer. Vortex for 10–15 seconds. Label the tube with the date of dilution. The diluted 6C Matrix Mix can be stored for up to 1 week at 2–10°C.
3. Vortex the diluted 6C Matrix Mix prepared in Step 2 for 10–15 seconds, then add 10µl to 500µl of Hi-Di™ formamide.

- Vortex the diluted 6C Matrix Mix with formamide prepared in Step 3 for 10–15 seconds, then add 15µl to each of the first four wells of an 8-well strip tube. After placing the septa on the 8-well strip tube, briefly centrifuge the plate to remove bubbles. Do not heat denature.

		Wells							
		1	2	3	4	5	6	7	8
Lane	A	Matrix Mix							
	B								
	C								
	D								

3.3 Assembling Sample Cartridge

- Place the 8-well strip tube into the strip base in Lane A with the samples in positions 1–4 (Figure 15).

Note: Lane names A to D and well numbers 1 to 8 are embossed on the strip base. Be sure to check the lane name and well numbers when placing the 8-well strip tube into the base.

- To complete the assembly, place the retainer over the strip in the strip base, aligning the lane names A to D and well numbers 1 to 8 on the retainer to those on the strip base and pressing until the retainer clicks into the strip base (Figure 16).

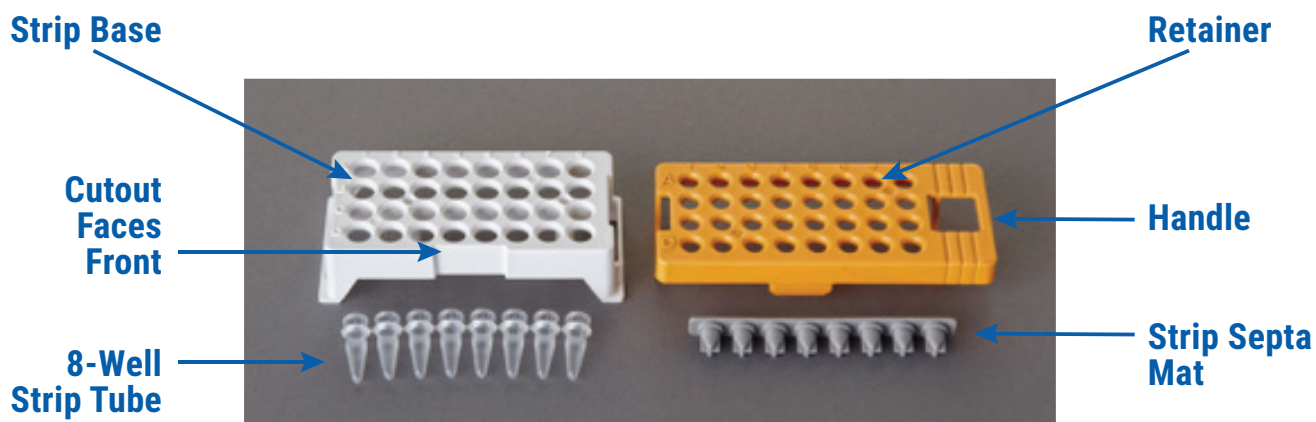


Figure 15. Assembling the Spectrum Compact Strip Base and Retainer.

15773TA

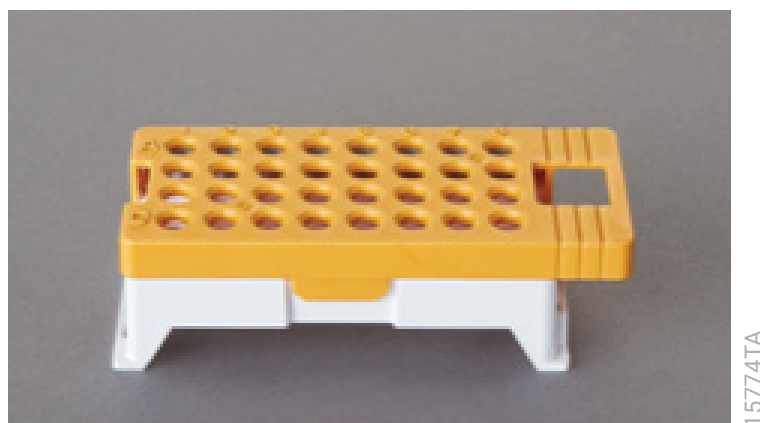


Figure 16. Assembled Spectrum Compact Sample Cartridge.

3.4 Instrument Preparation and Spectral Calibration

These instructions are intended as a guide for running PowerPlex® Matrix Standards on the Spectrum Compact CE System. They are not intended as comprehensive instructions for using the Spectrum Compact CE System. Refer to the *Spectrum Compact CE System Operating Manual* #TMD058 for more details on performing spectral calibration.

Notes:

- a. We have found that the use of fresh polymer and new capillary array results in an optimal spectral calibration.
- b. We do not recommend performing spectral calibration with expired reagents. Expired reagents should be replaced before performing a spectral calibration.
- c. Refer to the *Spectrum Compact CE System Operating Manual* #TMD058 for more details on installation of consumables, instrument maintenance and spatial calibration.

1. Select the **Consumables** icon in the Header on the 'Main Menu' screen (Figure 17). Ensure that the consumables are not expired and that adequate injections remain for consumables installed.



Figure 17. Spectrum Compact CE System Software 'Main Menu' screen.

2. Select the **Oven Temperature** icon in the Header on the 'Main Menu' screen as shown in Figure 18 to start preheating the oven temperature to 60°C. The temperature displayed will change as the temperature of the oven increases. When 60°C is reached, a check mark will appear adjacent to the temperature.

Note: We recommend you preheat the oven for at least 30 minutes prior to starting a run. The oven will automatically turn off after 2 hours if a run is not started.

Oven Temperature Icon

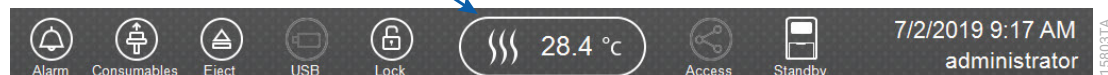


Figure 18. Preheating Oven.

3. Select **Calibration** on the maintenance portion of the 'Main Menu' screen (Figure 17) then select **Spectral Calibration** on the 'Maintenance Calibration' screen (Figure 19).

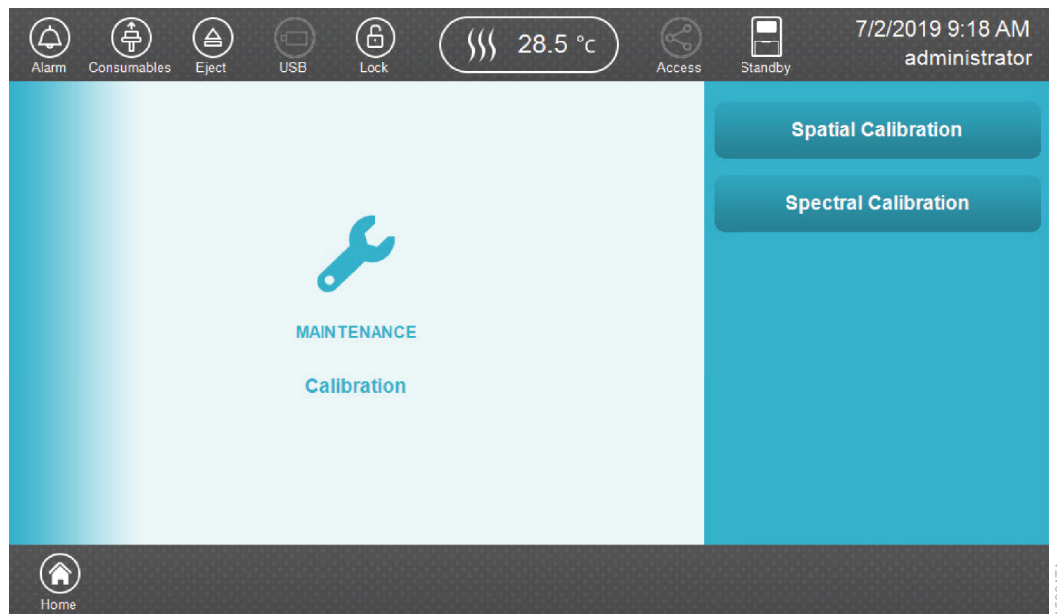


Figure 19. 'Maintenance Calibration' screen.

4. Use the scroll arrows on the right-hand side of the 'Dye Set List' screen (Figure 20) to find the correct Dye Set/Application Type/Polymer combination from the displayed list. To perform a spectral calibration using the PowerPlex® 6C Matrix Standard on Polymer4, select **Promega 6-dye** with "Fragment" and "Polymer4" as application and polymer types, respectively, then select **Calibration**. The 'Assemble the Cartridge' screen will open (Figure 21).

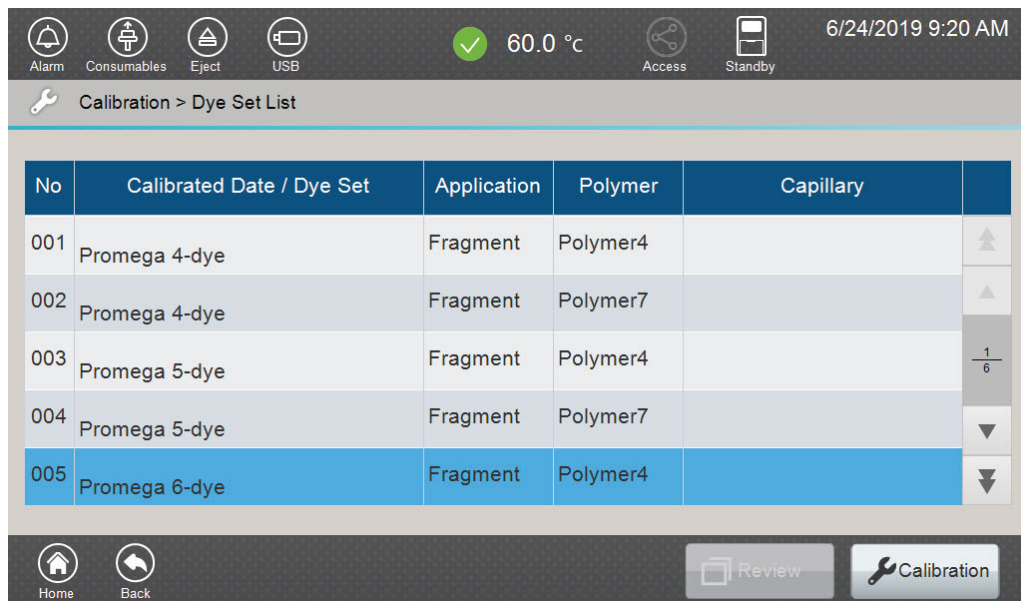


Figure 20. 'Dye Set List' screen.

5. Select **Next** on the 'Assemble the Cartridge' screen (Figure 21). A message window will open indicating that the autosampler is moving and telling the user to not open the door. In addition, the status indicator flashes green while the autosampler is moving. After autosampler movement is complete, the message window closes and the status indicator returns to a steady green.

Note: Do not open the front door of the Spectrum Compact CE System while the autosampler is in motion.

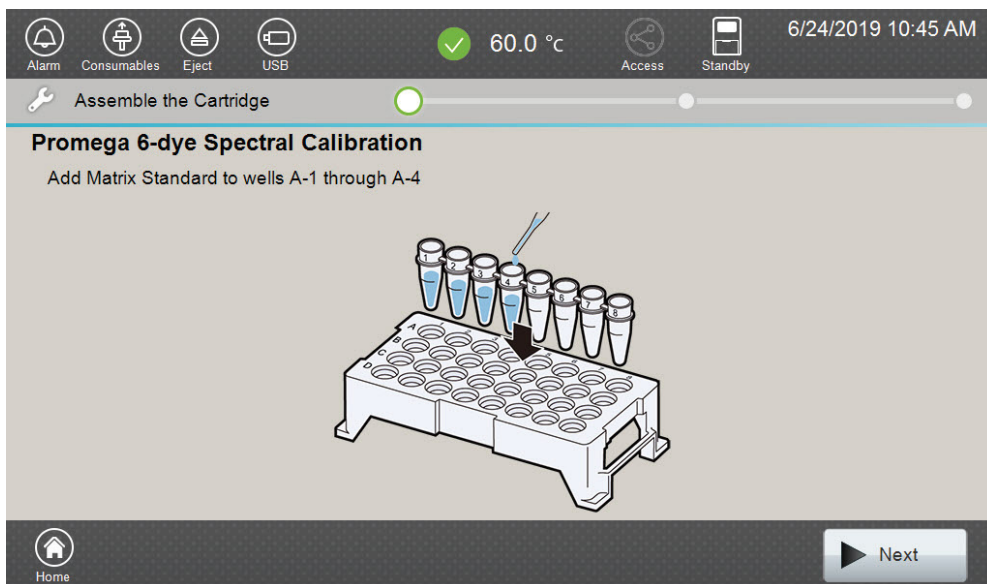


Figure 21. 'Assemble the Cartridge' screen.

6. Open the front door of the Spectrum Compact CE System and mount the sample cartridge on the autosampler following the instructions displayed on the 'Install the Cartridge' screen (Figure 22).

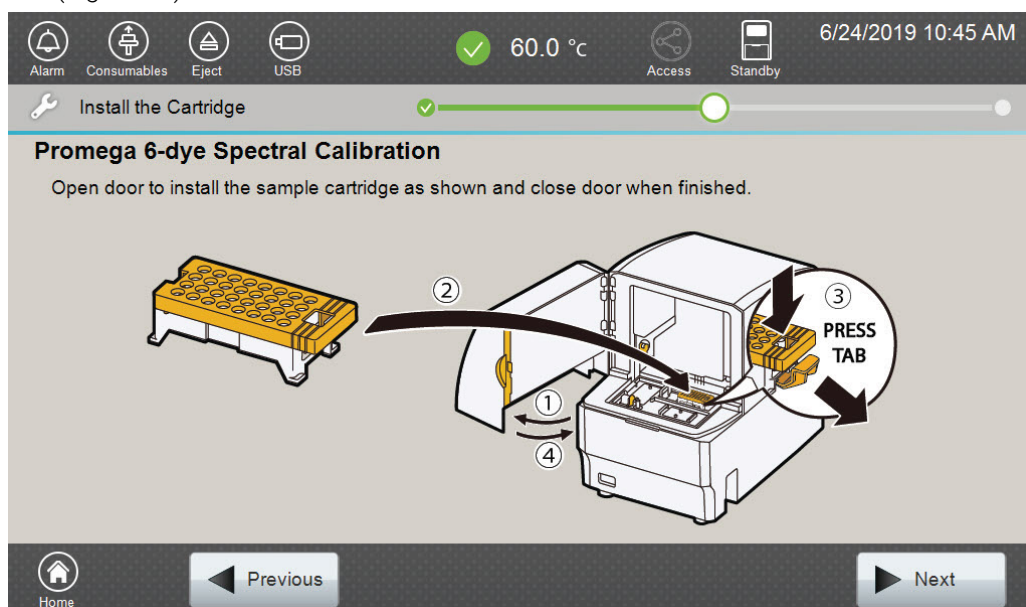


Figure 22. 'Install the Cartridge' screen.

- After mounting the sample cartridge on the autosampler, close the front door of the Spectrum Compact CE System and wait for the status indicator to stop flashing amber and turn steady green.

Note: Do not open the front door of the Spectrum Compact CE System while the autosampler is in motion.

- After the autosampler has returned to its home position, the 'Spectral Calibration' screen will automatically be displayed (Figure 23). Select **Run** to start the spectral calibration.

Note: The 'Raw Data' tab can be used to monitor the run.

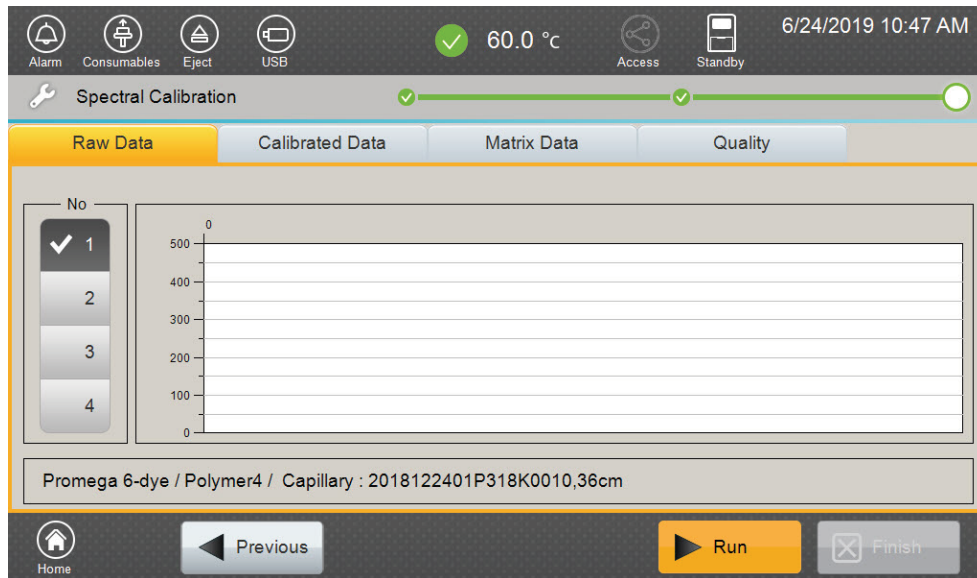


Figure 23. 'Spectral Calibration' screen.

3.5 Results

1. Following the run, the 'Raw Data' tab (Figure 24) will be displayed. The minimum peak height for spectral calibration is 500 relative fluorescent units (RFU) and the maximum peak height is 32,767RFU.

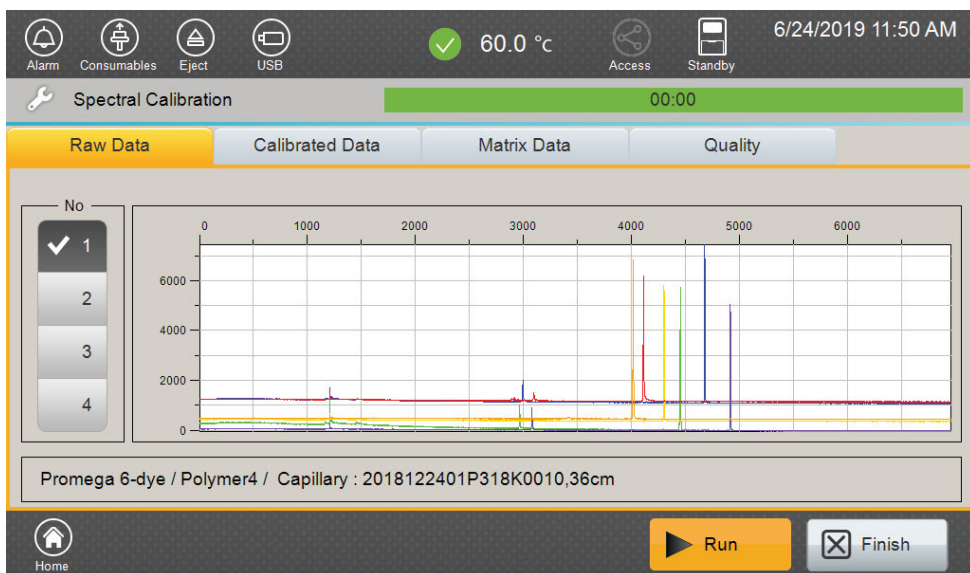


Figure 24. Spectral Calibration 'Raw Data' tab screen.

2. The 'Calibrated Data' tab can be used to view the matrix peaks with both baseline and spectral applied for each capillary (Figure 25).

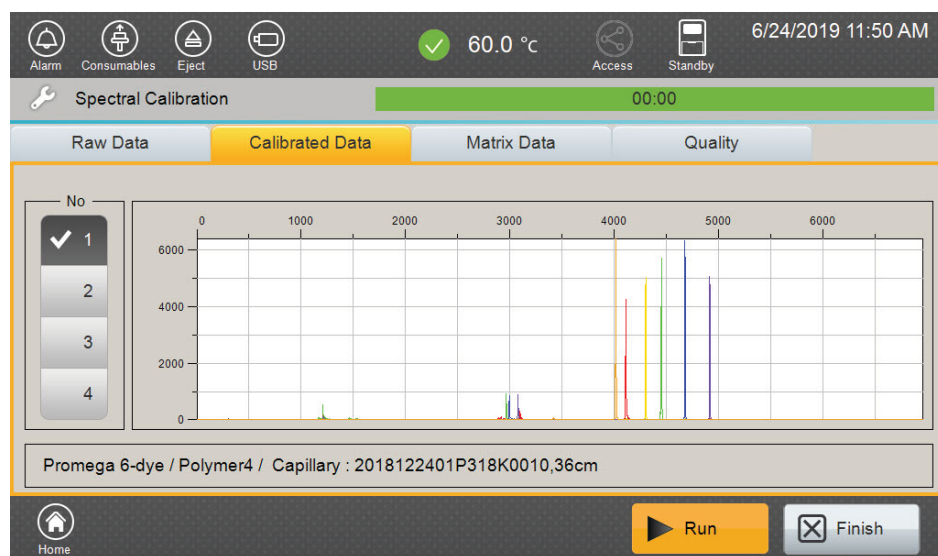


Figure 25. Spectral Calibration 'Calibrated Data' tab screen.

3. The 'Matrix Data' tab can be used to view emission spectra for each capillary (Figure 26).

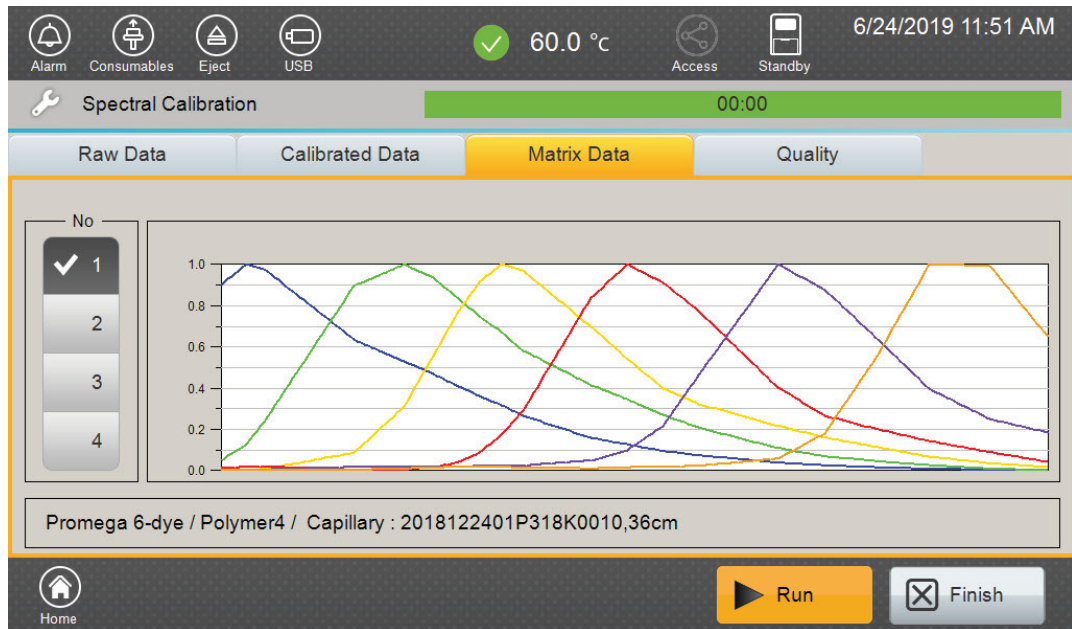


Figure 26. Spectral Calibration 'Matrix Data' tab screen.

4. Review the quality of the spectral calibration by selecting the 'Quality' tab (Figure 27).

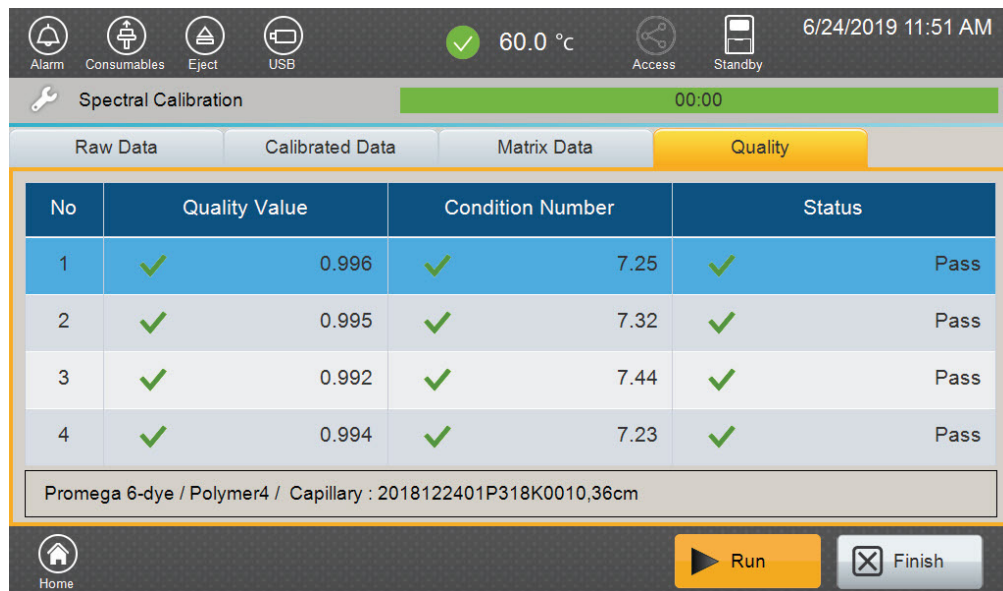


Figure 27. Spectral Calibration 'Quality' tab' screen.

5. Each capillary must meet the passing criteria of ≥ 0.95 for the Quality Value and < 8.5 for the Condition Number.

6. If one capillary fails to meet the criteria, it is possible to borrow spectral data from an adjacent capillary. Refer to the *Spectrum Compact CE System Operating Manual #TMD058* for details. If more than one capillary fails, the spectral must be rerun.

Notes:

- a. Selecting **Run** will rerun the spectral calibration.
- a. Refer to Section 5 for troubleshooting if more than one capillary fails to meet the criteria.
7. After reviewing the results, select **Finish**. This will open a confirmation window. Select **Yes** to apply the spectral calibration results (Figure 28). The spectral calibration results will not be saved unless you select **Yes** on this window.

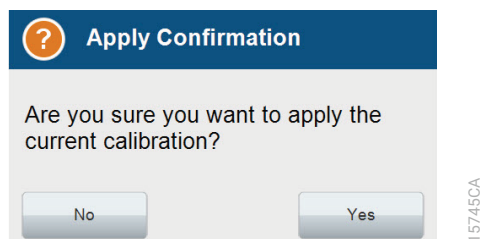


Figure 28. Apply spectral calibration confirmation window.

PowerPlex® 5C Spectral Calibration

4.1 Product Components and Storage Conditions

PRODUCT	SIZE	CAT.#
PowerPlex® 5C Matrix Standard	5 preps	DG4850

Not For Medical Diagnostic Use. Includes:

- 150µl 5C Matrix Mix
- 5 × 200µl Matrix Dilution Buffer

Storage Conditions

Upon receipt, store all components at -30°C to -10°C in a nonfrost-free freezer, protected from light. Do not store reagents in the freezer door, where the temperature can fluctuate. After the first use, store the PowerPlex® 5C Matrix Standard components at $2-10^{\circ}\text{C}$, protected from light. We strongly recommend that the PowerPlex® 5C Matrix Standard be stored with post-amplification reagents. The PowerPlex® 5C Matrix Standard is light-sensitive; dilute the 5C Matrix Mix in the Matrix Dilution Buffer in the provided amber tube. Store the diluted 5C Matrix Mix at $2-10^{\circ}\text{C}$ for up to 1 week.



Do not refreeze the PowerPlex® 5C Matrix Standard components.

Materials to be Supplied by the User

- centrifuge compatible with 8-tube strips
- aerosol-resistant pipette tips
- Spectrum Compact Capillary Cartridge, 4-Capillary 36cm (Cat.# CE2340)
- Spectrum Compact Polymer4 (Cat.# CE2304, CE2404)
- Spectrum Compact Buffer (Cat.# CE2300)
- Spectrum Compact Cathode Septa Mat (Cat.# CE2301)
- Spectrum Compact Cathode Retainer (Cat.# CE2302)
- Spectrum Compact Strip Base & Retainer, 32-Well (Cat.# CE2332)
- MicroAmp® Optical 8-Tube Strip, 0.2ml (Applied Biosystems, Cat.# 4316567)
- Strip Septa Mat, 8-Well (Cat.# CE2308)
- Hi-Di™ formamide (Applied Biosystems, Cat.# 4311320)

For additional information on performing spectral calibration, refer to the *Spectrum CE System Operating Manual #TMD058*.



The quality of formamide is critical. Use Hi-Di™ formamide. Freeze the formamide in aliquots at –20°C. Multiple freeze-thaw cycles or long-term storage at 4°C can cause breakdown of formamide. Poor-quality formamide can contain ions that compete with DNA during injection, which results in lower peak heights.



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.

Notes:

- a. Only use MicroAmp® Optical 8-Tube Strips, 0.2ml (Applied Biosystems, Cat.# 4316567) as a source of 8-well strip tubes. Use of other 8-well strip tubes may affect performance or damage the Spectrum Compact CE System.
- b. Wear gloves when handling consumables and sample cartridges.

4.2 Matrix Sample Preparation

1. At the first use, thaw the 5C Matrix Mix and Matrix Dilution Buffer completely. After the first use, store the reagents at 2–10°C, protected from light.
2. Vortex the 5C Matrix Mix for 10–15 seconds prior to use. Add 10µl of the 5C Matrix Mix to one tube of the Matrix Dilution Buffer. Vortex for 10–15 seconds. Label the tube with the date of dilution. The diluted 5C Matrix Mix can be stored for up to 1 week at 2–10°C.
3. Vortex the diluted 5C Matrix Mix prepared in Step 2 for 10–15 seconds, then add 10µl to 500µl of Hi-Di™ formamide.

- Vortex the diluted 5C Matrix Mix with formamide prepared in Step 3 for 10–15 seconds, then add 15µl to each of the first four wells of an 8-well strip tube. After placing the septa on the 8-well strip tube, briefly centrifuge the plate to remove bubbles. Do not heat denature.

		Wells							
		1	2	3	4	5	6	7	8
Lane	A	Matrix Mix							
	B								
	C								
	D								

4.3 Assembling Sample Cartridge

- Place the 8-well strip tube into the strip base in Lane A with the samples in positions 1–4 (Figure 29).

Note: Lane names A to D and well numbers 1 to 8 are embossed on the strip base. Be sure to check the lane name and well numbers when placing the 8-well strip tube into the base.

- To complete the assembly, place the retainer over the strip in the strip base, aligning the lane names A to D and well numbers 1 to 8 on the retainer to those on the strip base and pressing until the retainer clicks into the strip base (Figure 30).

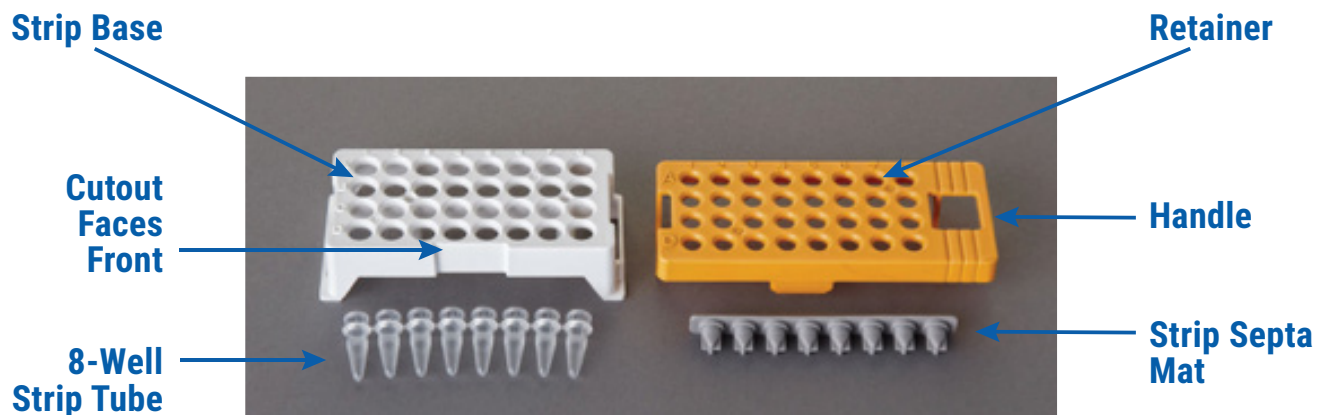


Figure 29. Assembling the Spectrum Compact Strip Base and Retainer.

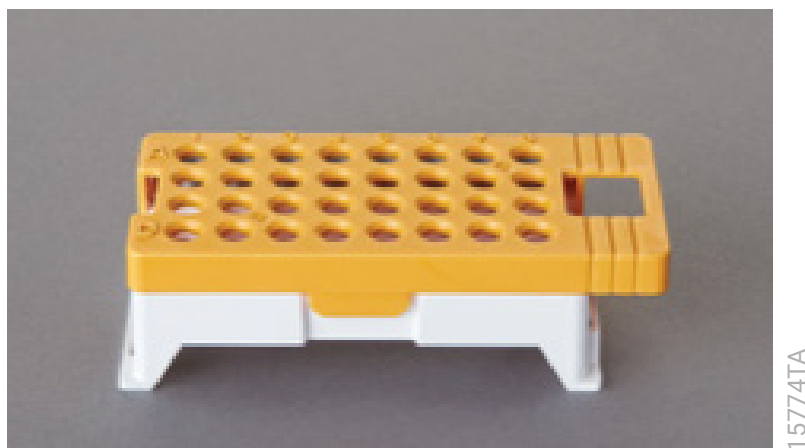


Figure 30. Assembled Spectrum Compact Sample Cartridge.

4.4 Instrument Preparation and Spectral Calibration

These instructions are intended as a guide for running PowerPlex® Matrix Standards on the Spectrum Compact CE System. They are not intended as comprehensive instructions for using the Spectrum Compact CE System. Refer to the *Spectrum Compact CE System Operating Manual #TMD058* for more details on performing spectral calibration.

Notes:

- a. We have found that the use of fresh polymer and new capillary array results in an optimal spectral calibration.
- b. We do not recommend performing spectral calibration with expired reagents. Expired reagents should be replaced before performing a spectral calibration.
- c. Refer to the *Spectrum Compact CE System Operating Manual #TMD058* for more details on installation of consumables, instrument maintenance and spatial calibration.

1. Select the **Consumables** icon in the Header on the 'Main Menu' screen (Figure 31). Ensure that the consumables are not expired and that adequate injections remain for consumables installed.



Figure 31. Spectrum Compact CE System Software 'Main Menu' screen.

2. Select the **Oven Temperature** icon in the Header on the 'Main Menu' screen, as shown in Figure 32, to start preheating the oven temperature to 60°C. The temperature displayed will change as the temperature of the oven increases. When 60°C is reached, a check mark will appear adjacent to the temperature.

Note: We recommend you preheat the oven for at least 30 minutes prior to starting a run. The oven will automatically turn off after 2 hours if a run is not started.

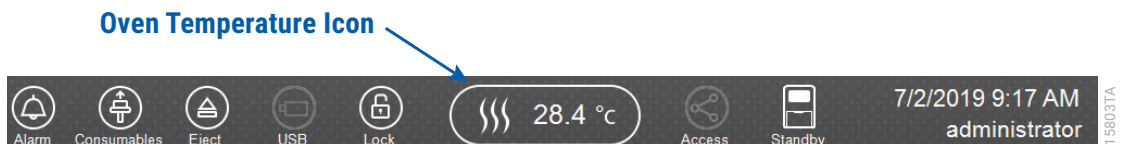


Figure 32. Preheating Oven.

3. Select **Calibration** on the maintenance portion of the 'Main Menu' screen (Figure 31) then select **Spectral Calibration** on the 'Maintenance Calibration' screen (Figure 33).

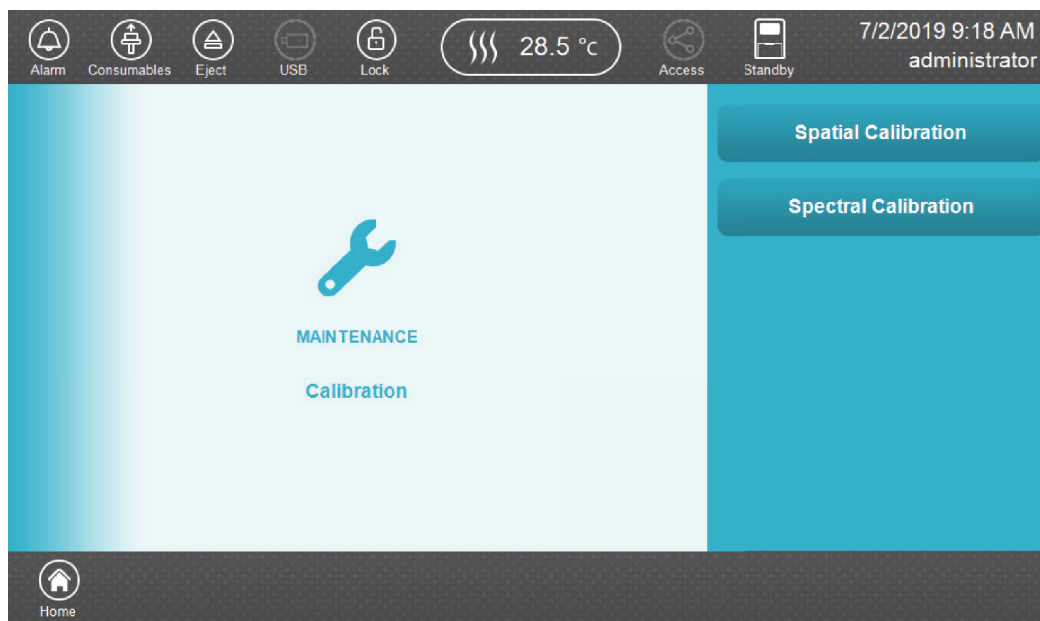


Figure 33. 'Maintenance Calibration' screen.

4. Use the scroll arrows on the right-hand side of the "Dye Set List" screen (Figure 34) to find the correct Dye Set/Application Type/Polymer combination from the displayed list. To perform a spectral calibration using the PowerPlex® 5C Matrix Standard on Polymer4, select **Promega 5-dye** with "Fragment" and "Polymer4" as application and polymer types, respectively, then select **Calibration**. The 'Assemble the Cartridge' screen will open (Figure 35).

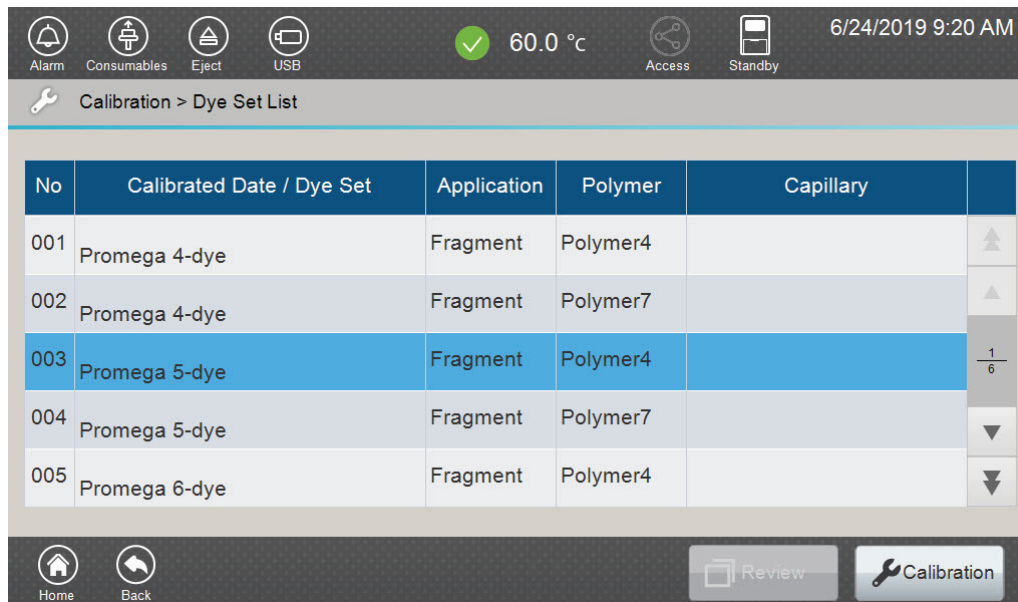


Figure 34. 'Dye Set List' screen.

- Select **Next** on the 'Assemble the Cartridge' screen (Figure 35). A message window will open indicating that the autosampler is moving and telling the user to not open the door. In addition, the status indicator flashes green while the autosampler is moving. After autosampler movement is complete, the message window closes and the status indicator returns to a steady green.

Note: Do not open the front door of the Spectrum Compact CE System while the autosampler is in motion.

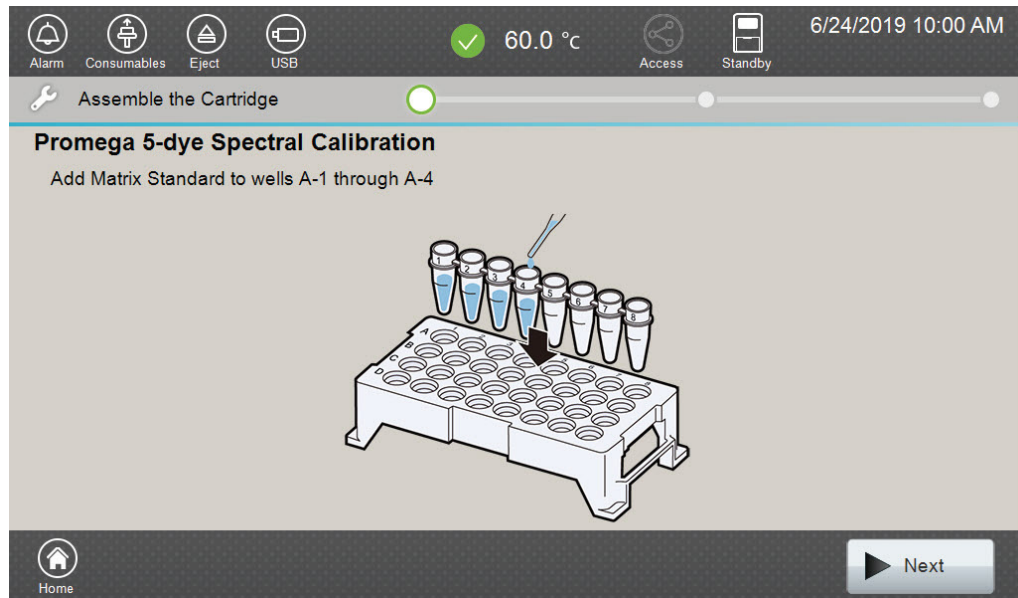


Figure 35. 'Assemble the Cartridge' screen.

- Open the front door of the Spectrum Compact CE System and mount the sample cartridge on the autosampler by following the instructions displayed on the 'Install the Cartridge' screen (Figure 36).

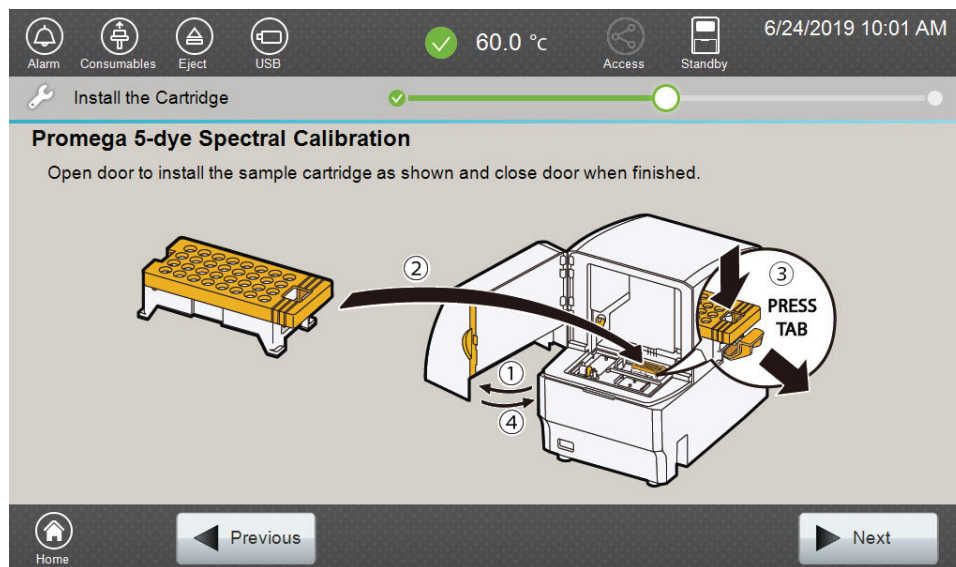


Figure 36. 'Install the Cartridge' screen.

- After mounting the sample cartridge on the autosampler, close the front door of the Spectrum Compact CE System and wait for the status indicator to stop flashing amber and turn steady green.

Note: Do not open the front door of the Spectrum Compact CE System while the autosampler is in motion.

- After the autosampler has returned to its home position, the 'Spectral Calibration' screen will automatically be displayed (Figure 37). Select **Run** to start the spectral calibration.

Note: The 'Raw Data' tab can be used to monitor the run.

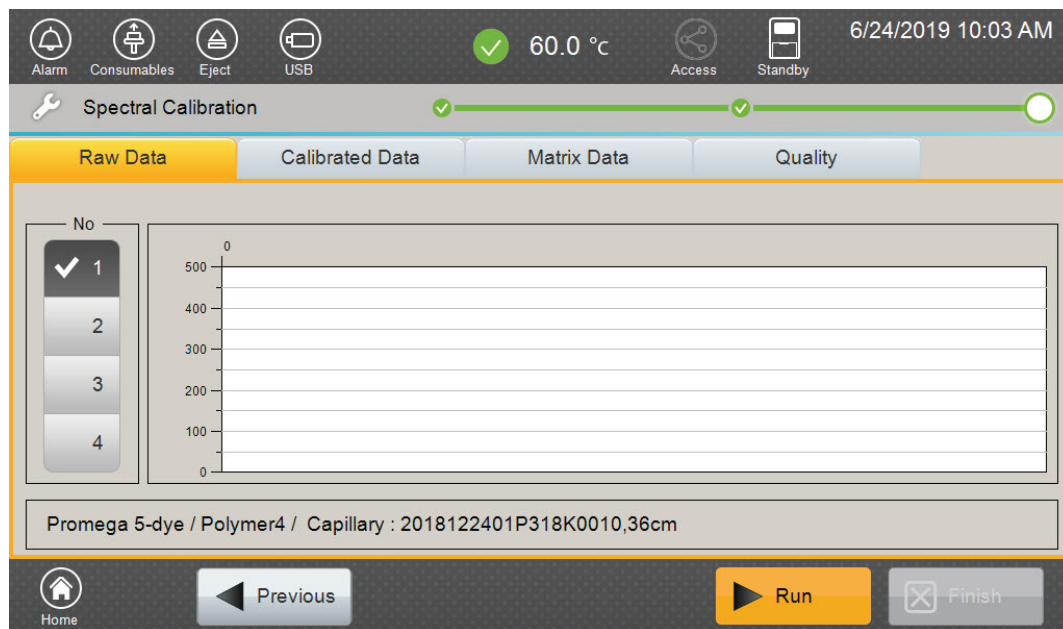


Figure 37. 'Spectral Calibration' screen.

4.5 Results

1. Following the run, the 'Raw Data' tab (Figure 38) will be displayed. The minimum peak height for spectral calibration is 500 relative fluorescent units (RFU) and the maximum peak height is 32,767RFU.

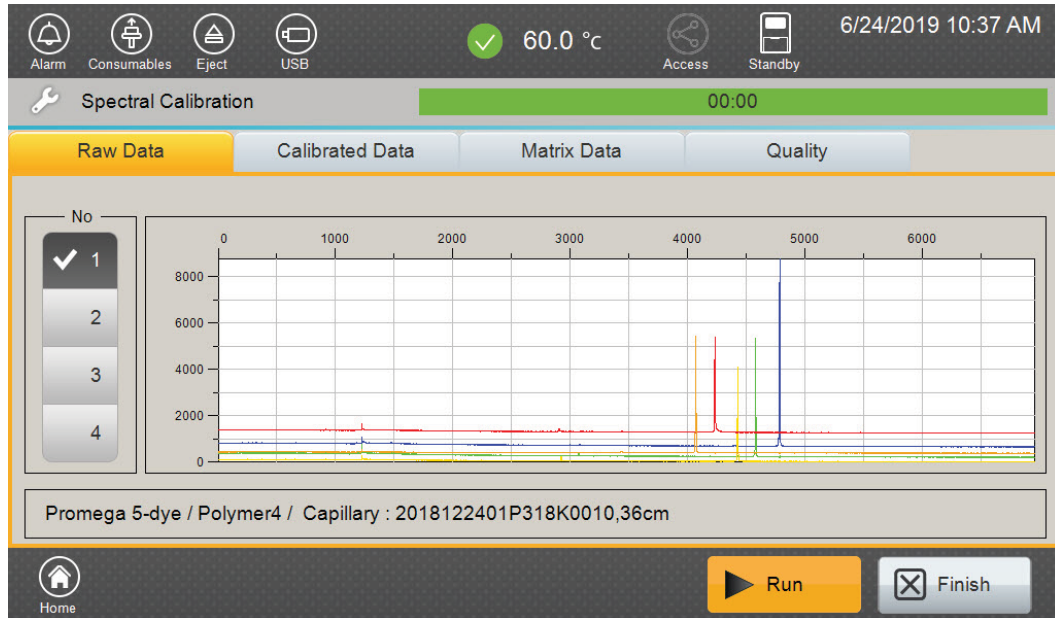


Figure 38. Spectral Calibration 'Raw Data' tab screen.

2. The 'Calibrated Data' tab can be used to view the matrix peaks with both baseline and spectral applied for each capillary (Figure 39).

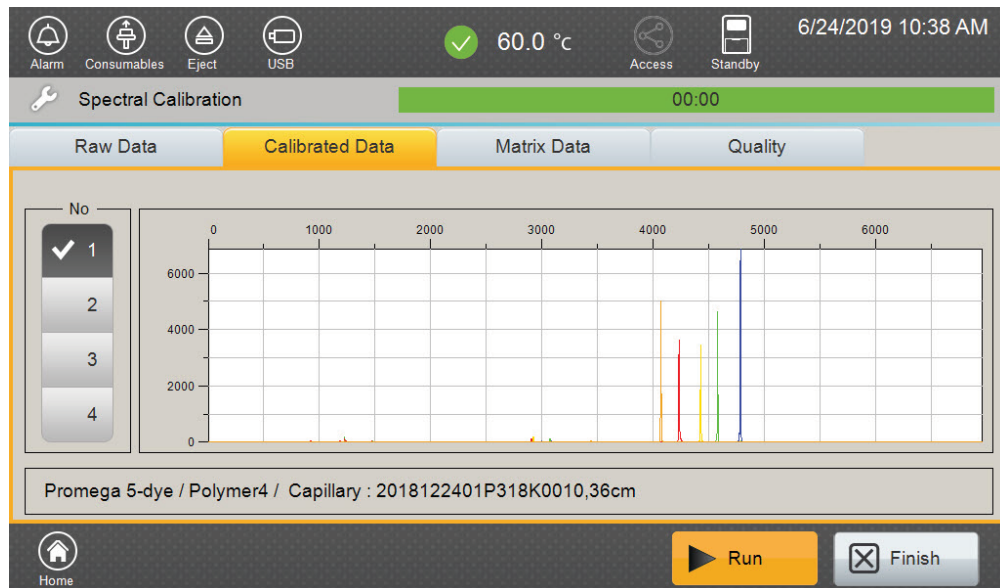


Figure 39. Spectral Calibration 'Calibrated Data' tab screen.

3. The 'Matrix Data' tab can be used to view emission spectra for each capillary (Figure 40).

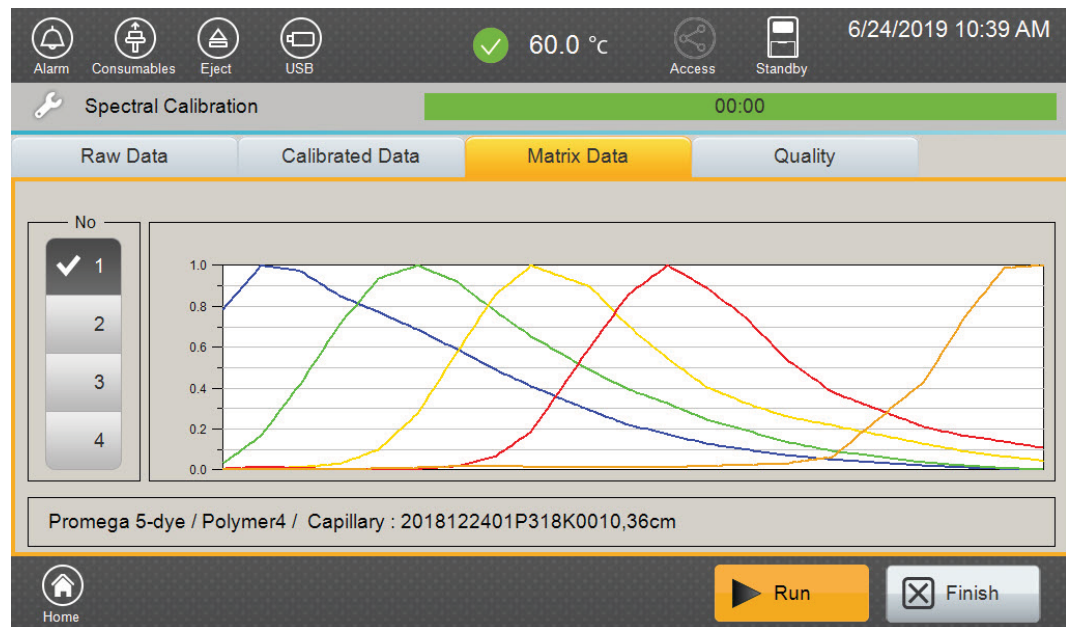


Figure 40. Spectral Calibration 'Matrix Data' tab screen.

4. Review the quality of the spectral calibration by selecting the 'Quality' tab (Figure 41).

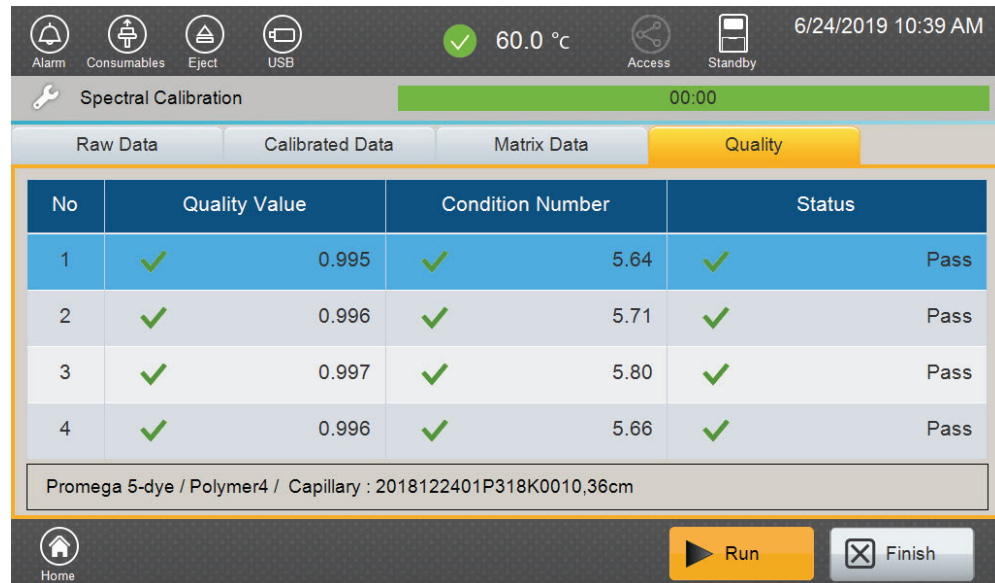


Figure 41. Spectral Calibration 'Quality' tab screen.

5. Each capillary must meet the passing criteria of ≥ 0.95 for the Quality Value and < 13.5 for the Condition Number.

6. If one capillary fails to meet the criteria, it is possible to borrow spectral data from an adjacent capillary. Refer to the *Spectrum Compact CE System Operating Manual #TMD058* for details. If more than one capillary fails, the spectral must be rerun.

Notes:

- a. Selecting **Run** will rerun the spectral calibration.
 - b. Refer to Section 5 for troubleshooting if more than one capillary fails to meet the criteria.
7. After reviewing the results, select **Finish**. This will open a confirmation window. Select **Yes** to apply the spectral calibration results (Figure 42). The spectral calibration results will not be saved unless you select **Yes** on this window.

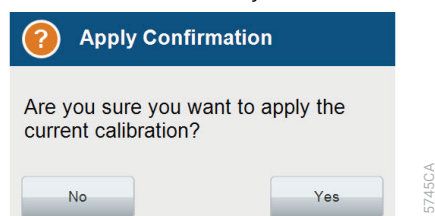


Figure 42. Apply spectral calibration confirmation window.

PowerPlex® 4C Spectral Calibration

5

5.1 Product Components and Storage Conditions

PRODUCT	SIZE	CAT.#
PowerPlex® 4C Matrix Standard	5 preps	DG4800

Not For Medical Diagnostic Use. Includes:

- 150µl 4C Matrix Mix
- 5 × 200µl Matrix Dilution Buffer

Storage Conditions

Upon receipt, store all components at -30°C to -10°C in a nonfrost-free freezer, protected from light. Do not store reagents in the freezer door, where the temperature can fluctuate. After the first use, store the PowerPlex® 4C Matrix Standard components at $2-10^{\circ}\text{C}$, protected from light. We strongly recommend that the PowerPlex® 4C Matrix Standard be stored with the post-amplification reagents. The PowerPlex® 4C Matrix Standard is light-sensitive; dilute the 4C Matrix Mix in the Matrix Dilution Buffer in the provided amber tube. Store the diluted 4C Matrix Mix at $2-10^{\circ}\text{C}$ for up to 1 week.



Do not refreeze the PowerPlex® 4C Matrix Standard components.

Materials to be Supplied by the User

- centrifuge compatible with 8-tube strips
- aerosol-resistant pipette tips
- Spectrum Compact Capillary Cartridge, 4-Capillary 36cm (Cat.# CE2340)
- Spectrum Compact Polymer4 (Cat.# CE2304, CE2404)
- Spectrum Compact Buffer (Cat.# CE2300)
- Spectrum Compact Cathode Septa Mat (Cat.# CE2301)
- Spectrum Compact Cathode Retainer (Cat.# CE2302)
- Spectrum Compact Strip Base & Retainer, 32-Well (Cat.# CE2332)
- MicroAmp® Optical 8-Tube Strip, 0.2ml (Applied Biosystems, Cat.# 4316567)
- Strip Septa Mat, 8-Well (Cat.# CE2308)
- Hi-Di™ formamide (Applied Biosystems, Cat.# 4311320)

For additional information on performing spectral calibration, refer to the *Spectrum CE System Operating Manual #TMD058*.



The quality of formamide is critical. Use Hi-Di™ formamide. Freeze the formamide in aliquots at -20°C. Multiple freeze-thaw cycles or long-term storage at 4°C can cause breakdown of formamide. Poor-quality formamide can contain ions that compete with DNA during injection, which results in lower peak heights.



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.

Notes:

- a. Only use MicroAmp® Optical 8-Tube Strips, 0.2ml (Applied Biosystems, Cat.# 4316567) as a source of 8-well strip tubes. Use of other 8-well strip tubes may affect performance or damage the Spectrum Compact CE System.
- b. Wear gloves when handling consumables and sample cartridges.

5.2 Matrix Sample Preparation

1. At the first use, thaw the 4C Matrix Mix and Matrix Dilution Buffer completely. After the first use, store the reagents at 2–10°C, protected from light.
2. Vortex the 4C Matrix Mix for 10–15 seconds prior to use. Add 10µl of the 4C Matrix Mix to one tube of the Matrix Dilution Buffer. Vortex for 10–15 seconds. Label the tube with the date of dilution. The diluted 4C Matrix Mix can be stored for up to 1 week at 2–10°C.
3. Vortex the diluted 4C Matrix Mix prepared in Step 2 for 10–15 seconds, then add 10µl to 500µl of Hi-Di™ formamide.

- Vortex the diluted 4C Matrix Mix with formamide prepared in Step 3 for 10–15 seconds, then add 15µl to each of the first four wells of an 8-well strip tube. After placing the septa on the 8-well strip tube, briefly centrifuge the plate to remove bubbles. Do not heat denature.

		Wells							
		1	2	3	4	5	6	7	8
Lane	A	Matrix Mix							
	B								
	C								
	D								

5.3 Assembling Sample Cartridge

- Place the 8-well strip tube into the strip base in Lane A with the samples in positions 1–4 (Figure 43).

Note: Lane names A to D and well numbers 1 to 8 are embossed on the strip base. Be sure to check the lane name and well numbers when placing the 8-well strip tube into the base.

- To complete the assembly, place the retainer over the strip in the strip base, aligning the lane names A to D and well numbers 1 to 8 on the retainer to those on the strip base and pressing until the retainer clicks into the strip base (Figure 44).



Figure 43. Assembling the Spectrum Compact Strip Base and Retainer.

15773TA

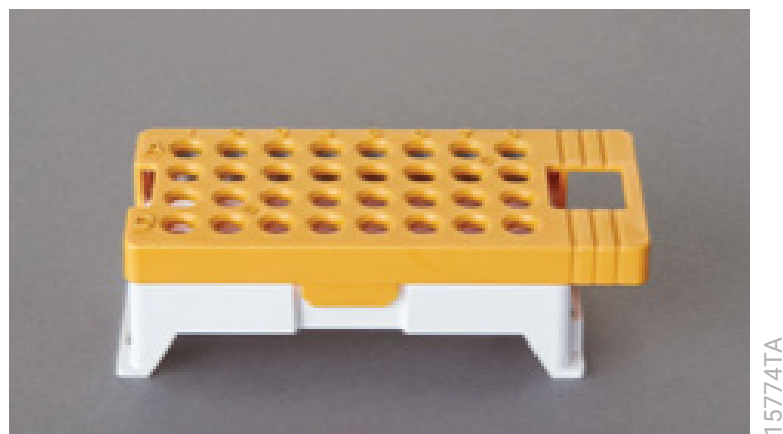


Figure 44. Assembled Spectrum Compact Sample Cartridge.

5.4 Instrument Preparation and Spectral Calibration

These instructions are intended as a guide for running PowerPlex® Matrix Standards on the Spectrum Compact CE System. They are not intended as comprehensive instructions for using the Spectrum Compact CE System. Refer to the *Spectrum Compact CE System Operating Manual #TMD058* for more details on performing spectral calibration.

Notes:

- a. We have found that the use of fresh polymer and new capillary array results in an optimal spectral calibration.
- b. We do not recommend performing spectral calibration with expired reagents. Expired reagents should be replaced before performing a spectral calibration.
- c. Refer to the *Spectrum Compact CE System Operating Manual #TMD058* for more details on installation of consumables, instrument maintenance and spatial calibration.

1. Select the **Consumables** icon in the Header on the 'Main Menu' screen (Figure 45). Ensure that the consumables are not expired and that adequate injections remain for consumables installed.



Figure 45. Spectrum Compact CE System Software 'Main Menu' screen.

2. Select the **Oven Temperature** icon in the Header on the 'Main Menu' screen, as shown in Figure 46, to start preheating the oven temperature to 60°C. The temperature displayed will change as the temperature of the oven increases. When 60°C is reached, a check mark will appear adjacent to the temperature.

Note: We recommend you preheat the oven for at least 30 minutes prior to starting a run. The oven will automatically turn off after 2 hours if a run is not started.

Oven Temperature Icon

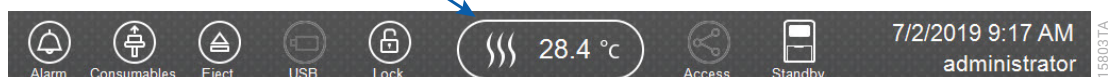


Figure 46. Preheating Oven.

3. Select **Calibration** on the maintenance portion of the 'Main Menu' screen (Figure 45) then select **Spectral Calibration** on the 'Maintenance Calibration' screen (Figure 47).

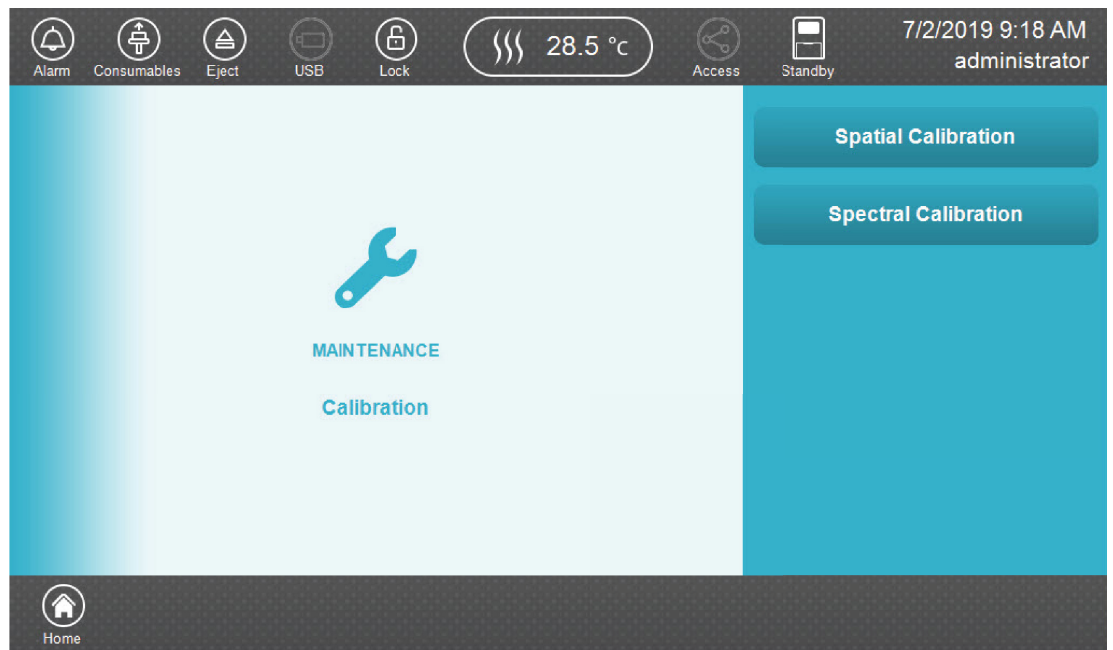


Figure 47. 'Maintenance Calibration' screen.

4. Use the scroll arrows on the right-hand side of the 'Dye Set List' screen (Figure 48) to find the correct Dye Set/Application Type/Polymer combination from the displayed list. To perform a spectral calibration using the PowerPlex® 4C Matrix Standard on Polymer4, select **Promega 4-dye** with "Fragment" and "Polymer4" as application and polymer types, respectively, then select **Calibration**. The 'Assemble the Cartridge' screen will open (Figure 49).

No	Calibrated Date / Dye Set	Application	Polymer	Capillary
001	Promega 4-dye	Fragment	Polymer4	
002	Promega 4-dye	Fragment	Polymer7	
003	Promega 5-dye	Fragment	Polymer4	1/6
004	Promega 5-dye	Fragment	Polymer7	
005	Promega 6-dye	Fragment	Polymer4	

Figure 48. 'Dye Set List' screen.

5. Select **Next** on the 'Assemble the Cartridge' screen (Figure 49). A message window will open indicating that the autosampler is moving and telling the user to not open the door. In addition, the status indicator flashes green while the autosampler is moving. After autosampler movement is complete, the message window closes and the status indicator returns to a steady green.

Note: Do not open the front door of the Spectrum Compact CE System while the autosampler is in motion.

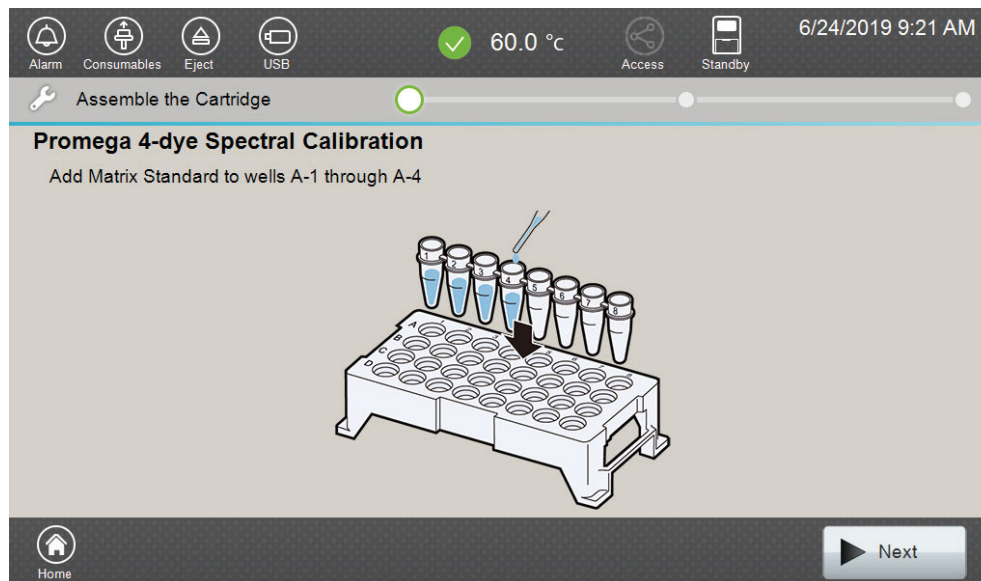


Figure 49. 'Assemble the Cartridge' screen.

6. Open the front door of the Spectrum Compact CE System and mount the sample cartridge on the autosampler by following the instructions displayed on the 'Install the Cartridge' screen (Figure 50).

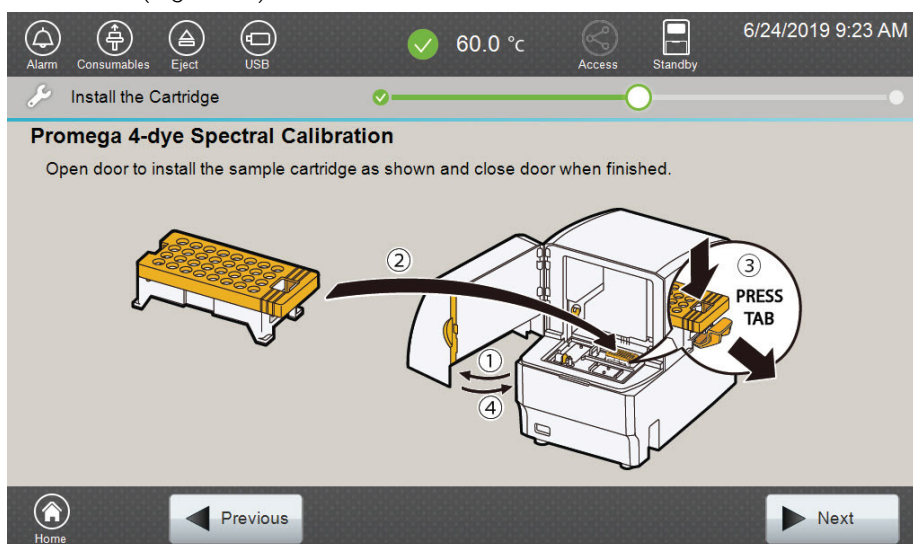


Figure 50. 'Install the Cartridge' screen.

- After mounting the sample cartridge on the autosampler, close the front door of the Spectrum Compact CE System and wait for the status indicator to stop flashing amber and turn steady green.

Note: Do not open the front door of the Spectrum Compact CE System while the autosampler is in motion.

- After the autosampler has returned to its home position, the 'Spectral Calibration' screen will automatically be displayed (Figure 51). Select **Run** to start the spectral calibration.

Note: The 'Raw Data' tab can be used to monitor the run.

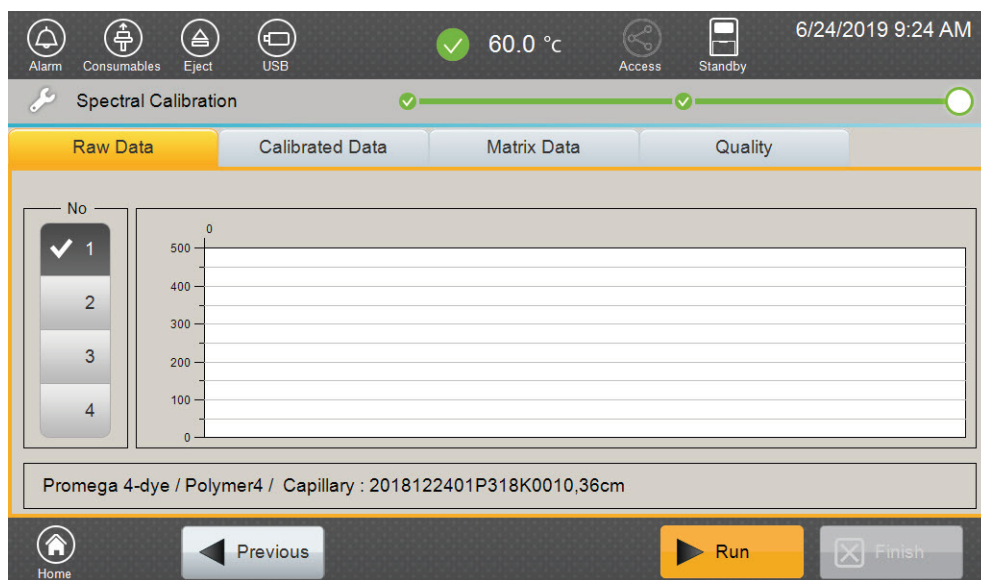


Figure 51. 'Spectral Calibration' screen.

5.5 Results

1. Following the run, the 'Raw Data' tab (Figure 52) will be displayed. The minimum peak height for spectral calibration is 500 relative fluorescent units (RFU) and the maximum peak height is 32,767RFU.

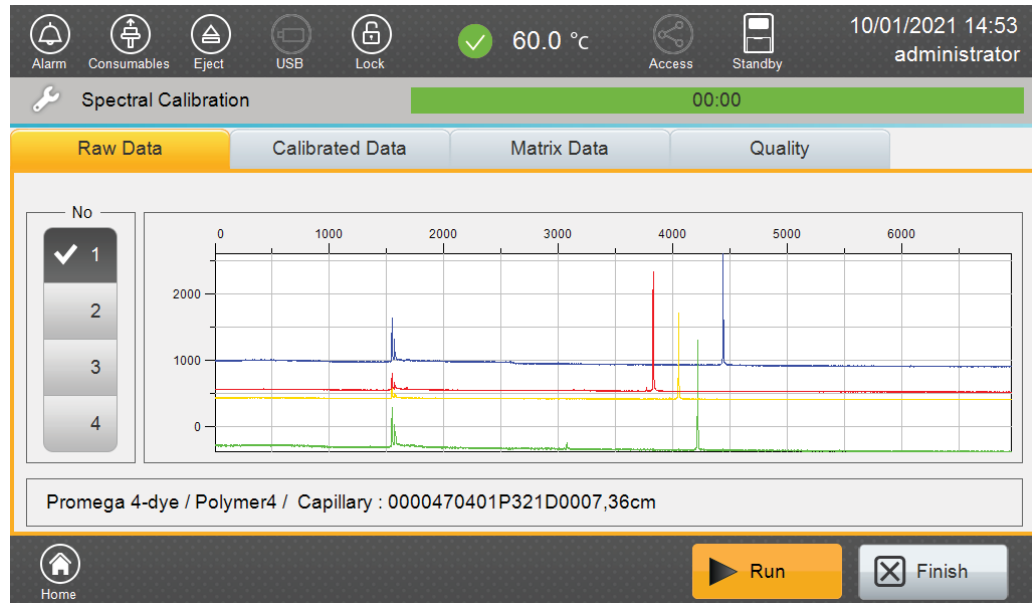


Figure 52. Spectral Calibration 'Raw Data' tab screen.

2. The 'Calibrated Data' tab can be used to view the matrix peaks with both baseline and spectral applied for each capillary (Figure 53).

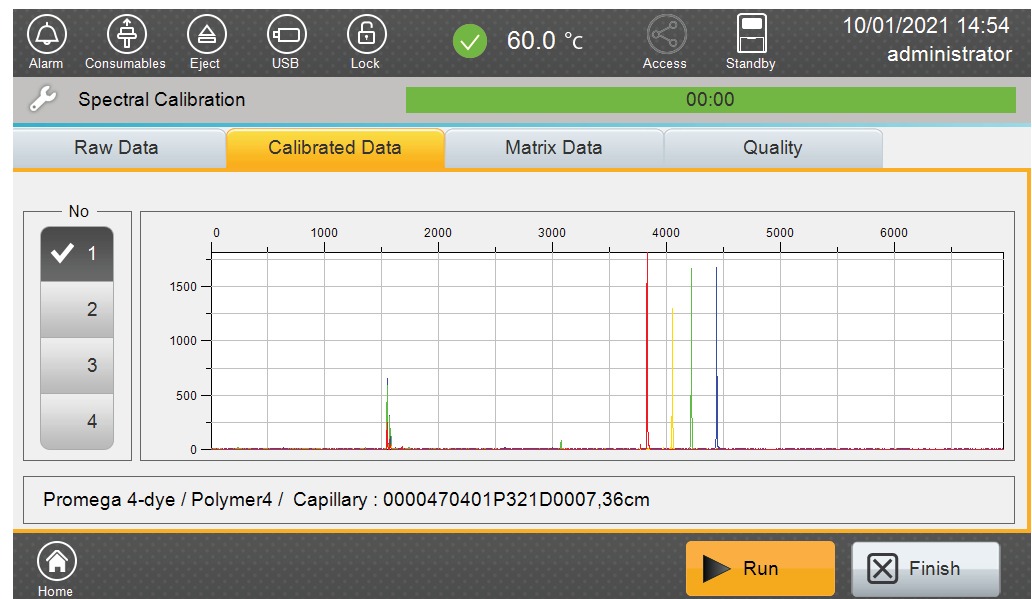


Figure 53. Spectral Calibration 'Calibrated Data' tab screen.

- The 'Matrix Data' tab can be used to view emission spectra for each capillary (Figure 54).

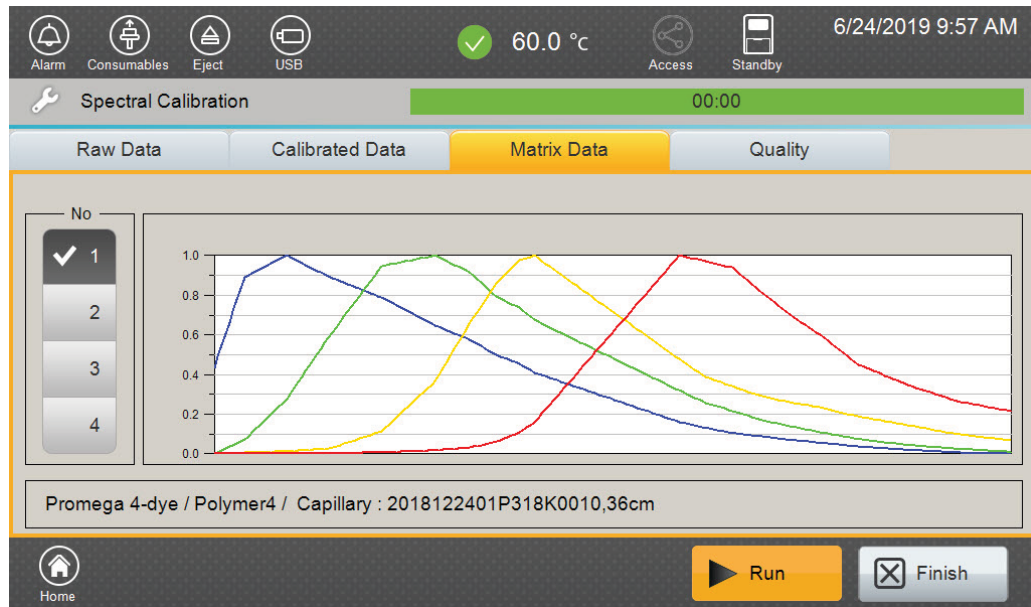


Figure 54. Spectral Calibration 'Matrix Data' tab screen.

- Review the quality of the spectral calibration by selecting the 'Quality' tab (Figure 55).

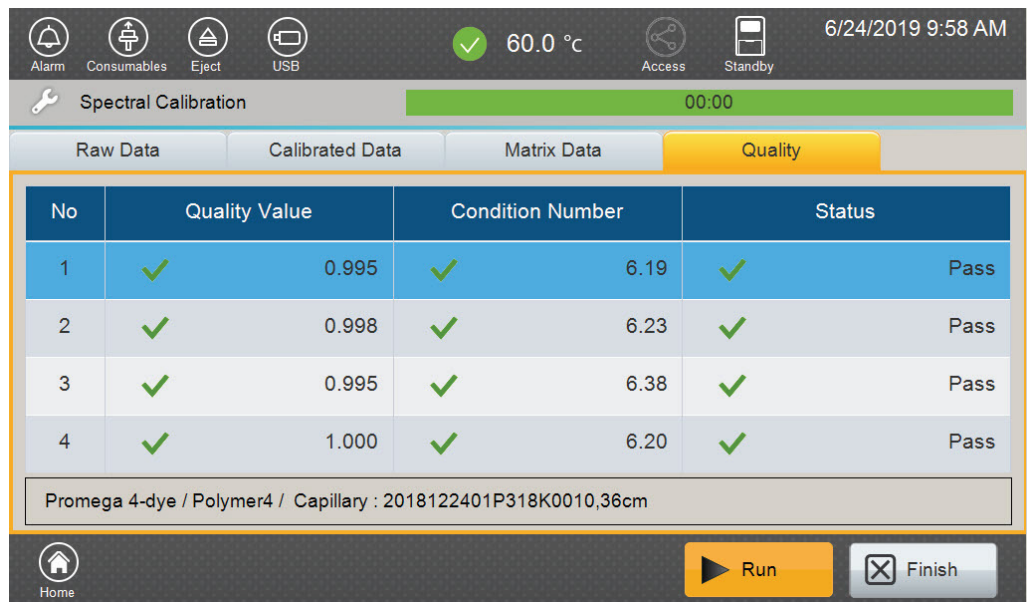


Figure 55. Spectral Calibration 'Quality' tab screen.

- Each capillary must meet the passing criteria of ≥ 0.95 for the Quality Value and < 8.5 for the Condition Number.

6. If one capillary fails to meet the criteria, it is possible to borrow spectral data from an adjacent capillary. Refer to the *Spectrum Compact CE System Operating Manual #TMD058* for details. If more than one capillary fails, the spectral must be rerun.

Notes:

- a. Selecting **Run** will rerun the spectral calibration.
 - b. Refer to Section 5 for troubleshooting if more than one capillary fails to meet the criteria.
7. After reviewing the results, select **Finish**. This will open a confirmation window. Select **Yes** to apply the spectral calibration results (Figure 56). The spectral calibration result will not be saved unless you select **Yes** on this window.

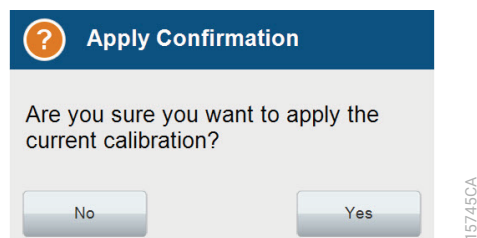


Figure 56. Apply spectral calibration confirmation window.

Troubleshooting

For questions not addressed here, please contact your local Promega Branch Office or Distributor. Contact information available at: www.promega.com E-mail: genetic@promega.com

Symptoms	Causes and Comments
Fewer than three capillaries passed the spectral calibration	Poor-quality formamide was used. The quality of formamide is critical. Freeze formamide in aliquots at -20°C . Multiple freeze-thaw cycles or 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.
	Matrix Mix was too dilute. Matrix Mix that is too dilute will result in low spectral calibration peak heights ($<500\text{RFU}$), which may result in spectral calibration failure. Increase the volume of diluted Matrix Mix added to the formamide during sample preparation.
	Diluted Matrix Mix stored longer than one week at 4°C or at the incorrect temperature. Prepare a fresh dilution.
	Matrix Mix was too concentrated. Matrix Mix that is too concentrated may result in spectral calibration failure due to saturated peaks, bleed-through or over-subtraction in other dye colors. Decrease the volume of diluted Matrix Mix added to the formamide during matrix sample preparation. Ensure you are using matrix standard that has been diluted appropriately.
	Carryover from previous injection detected as matrix peak. Replace cathode buffer cartridge septa. Use fresh reagents.
	For best spectral calibration results, use fresh polymer and fresh buffer.
	Use an array with fewer than 300 injections.
	Poor quality 8-well strip tubes. For best results, use MicroAmp [®] Optical 8-Tube Strip, 0.2ml (Applied Biosystems, Cat.# 4316567).

<p>Elevated spectral bleedthrough in one or more capillaries</p>	<p>If elevated spectral bleedthrough is observed in one or more capillaries after installing a new capillary cartridge, reinstall the capillary cartridge. Completely remove the capillary cartridge from the oven and reinstall as indicated in Section 3.2 of the <i>Spectrum Compact CE System Operating Manual #TMD058</i>. Simply lift the capillary cartridge completely out of the oven by its yellow knob and reinstall immediately back in the oven. Repositioning the detection unit of the capillary cartridge into the detection window of the oven during reinstallation can improve spectral performance.</p> <p>Note: Perform a new spatial and spectral calibration after uninstalling and reinstalling the capillary cartridge.</p>
--	---

Summary of Changes

The following changes were made to the 9/23 revision of this document:

1. Added new Section 2 for Cat.# DG5010 and renumbered subsequent sections and associated figures.
2. Added Spectrum Compact Polymer4 Cat.# CE2404.
3. Edited Section 7.
4. Updated disclaimers and document font.

^(a)U.S. Pat. No. 9,139,868, European Pat. No. 2972229, Japanese Pat. No. 6367307 and other patents pending.

^(b)AQA-8C, TMR-8C, CXR-8C, TOM-8C, WEN-8C and CCO-8C dyes are proprietary.

^(c)TMR-6C, CXR-6C, TOM-6C and WEN dyes are proprietary.

^(d)TMR-ET, CXR-ET and WEN dyes are proprietary.

© 2020–2023 Promega Corporation. All Rights Reserved.

PowerPlex is a registered trademark of Promega Corporation.

Applied Biosystems is a registered trademark of Applied Biosystems, L.L.C. MicroAmp is a registered trademark and Hi-Di is a trademark 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. 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.