

Agilent 7100 Capillary Electrophoresis System in Empower

Technical Note

This technical guide describes the configuration and implementation of the G7100A Capillary Electrophoresis in Empower 3.



Agilent Technologies

Contents

Introduction	3
Set up and Configuration of CE in Empower 3	5
CE Instrument set up	5
The CE Driver	5
Various Configuration Steps for CE in Empower	5
Using the CE in Empower	12
The CE Status Dashboard in Empower – Direct Control	12
Creating an Instrument Method	15
Replenishment and Conditioning	25
Indirect detection	28
Application of external pressure	28
Empower features for CE Instrument Control	33
Known Limitations/Not tested configurations:	33
Extended CE driver features available in Agilent OpenLAB CDS ChemStation Edition	34
Additional Information	36
Waters	36
APPENDIX - Examples	37
Example: Single test run	37
Example: Multiple injections using the IQ Kit for CE	39
Example: Indirect detection using the Plating Bath Buffer	41

Introduction

Waters Instrument Control Software (ICS) is Waters Corporation's adoption of the Agilent Instrument Control Framework (ICF) for their data systems. This guide describes how to configure and use the 7100 Capillary Electrophoresis (CE) in Waters Empower environment.

Empower 3 offers all key functions to control the G7100A Agilent Capillary Electrophoresis system. All established techniques are feasible, such as the use of direct and indirect detection modes, use of special electrolyte systems (containing micelles or gel matrices) and also capillary isoelectric focusing (cIEF) or capillary electrochromatography (CEC). The application of external high pressure is supported.

For data analysis, Empower offers processing tools for peak integration and calculation of concentrations in CE specific mode.

Table 1 Supported and unsupported configurations using CE in Empower

Waters Instrument Control Packages	Agilent ICF/ Agilent LC Driver	CE support in Waters Empower 3
ICF Support v2.2 #667005584	A.02.04 A.02.14	<i>Supported</i> G7100A Capillary Electrophoresis, with the internal components: <ul style="list-style-type: none"> • G7150A CE Mainframe • G7151A DAD Min. Firmware B.06.73 (older firmware does not support the use of RFID Tags for the DAD UV-Lamp)
ICF Support v2.2 #667005449	A.02.03 DU2 A.02.13	<i>Not supported</i> G1600 Capillary Electrophoresis
And any lower integration		

Introduction

NOTE

Ensure that the Agilent LC/CE modules in the system *meet or exceed* the minimum firmware requirements specified by the 3rd-party CDS software vendor and Agilent's firmware set/firmware interoperability requirements. Agilent recommends using the latest available firmware set.

<http://www.agilent.com/en-us/firmwareDownload?whid=69761>

Set up and Configuration of CE in Empower 3

CE Instrument set up

- 1 Close Empower.
- 2 Set up the CE system.
- 3 The CE uses LAN communication, e.g. via the multiport ethernet card. Connect the CE to the LAC/E box. If another instrument is connected to the LAC/E, connect via switch or hub.
- 4 Switch the instrument on.

The CE Driver

The CE driver is *part of* the LC Driver and not listed as a separate component. Therefore, Agilent ICF consists of two components; the Waters integration adds one additional component. All three components must be present in **Control Panel > Programs and Features**:

 Agilent Instrument Control Framework - LC Drivers A.02.14	Agilent Technologies	12/21/2016	111 MB	2.14.115
 Agilent Instrument Control Framework A.02.04	Agilent Technologies	12/21/2016	48.2 MB	2.4.124
 Agilent LC	Waters Corporation	12/20/2016	16.8 MB	2.2.0.0

Figure 1 Components in Programs and Feature

Various Configuration Steps for CE in Empower

To run CE in Empower the following items are required:

- A CE specific project for data collection
- The configuration of the CE instrument

Creating a CE Project

Configure a specific CE project to review the acquired data later.

- 1 Select **File > New > Project** to open the **New Project Wizard**.
- 2 Follow the instructions in the wizard screen and proceed with default settings if no other values are given via Standard Operating Procedure (SOP).

- 3 In the **New Project Wizard > Options** window make sure to enable **CE/CIA**.

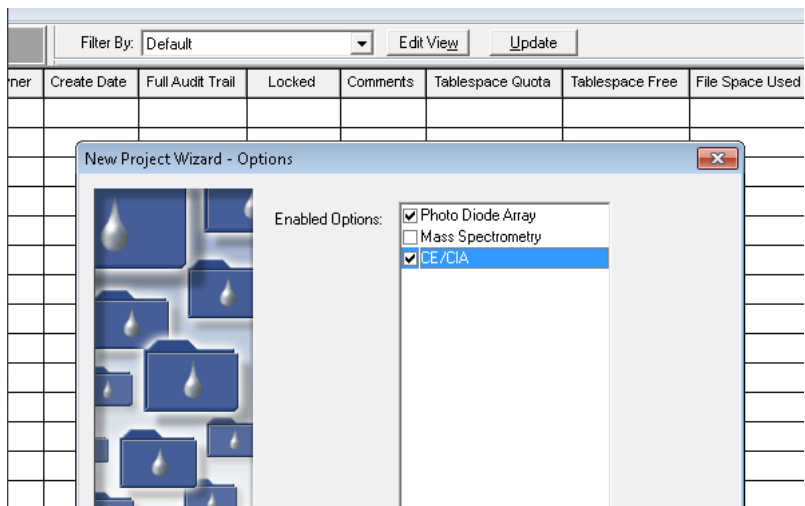


Figure 2 Enabled Options for the project

PreConfiguration of CE in Empower

For the following configuration options the PreConfiguration Utility is required to access the CE configuration in the CE configuration editor:

- Defining the **Pressure Unit** (see also chapter Known Limitations on page 33)
- Defining the **Temperature Control** mode (**Temperature Control** switched on by default)
- Defining the **Analog In Signal** (see also chapter Known Limitations on page 33)

NOTE

Selecting PSI as pressure unit, the CE status dashboard shows the pressure in PSI while the online plot and the resulting Aux Traces on the Chromatogram offer the values in mbar/bar.

Software required ICF Support v2.2 (See chapter Additional Information on page 36 for Waters support documentation TECN134936402).

- 1 In the Empower Configuration Manager, select **Tools > Agilent PreConfiguration**.

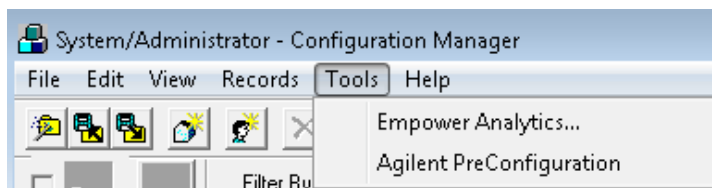


Figure 3 Configuration Manager

- 2 Enter the IP address or host name of the LAC/E box that your instrument is connected to into the pop-up screen **Configuration Directory** and click **Connect**.

Note

Do not enter the IP address of the instrument here.

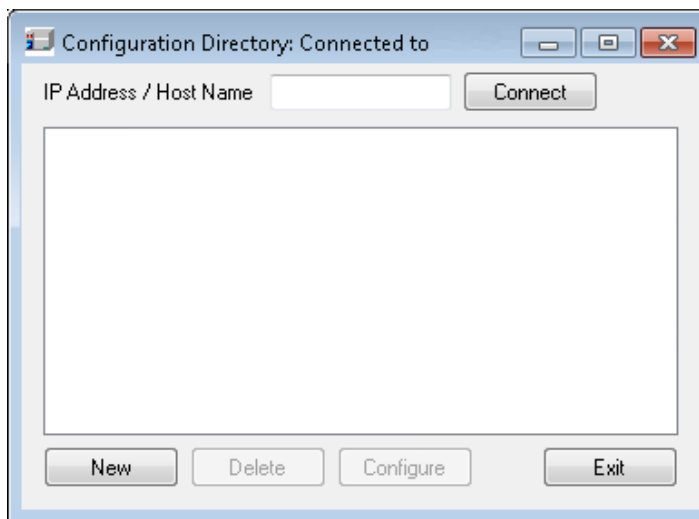


Figure 4 IP address to connect

Set up and Configuration of CE in Empower 3

- 3 Once the IP address is connected, click **New** to open the PreConfiguration Utility.

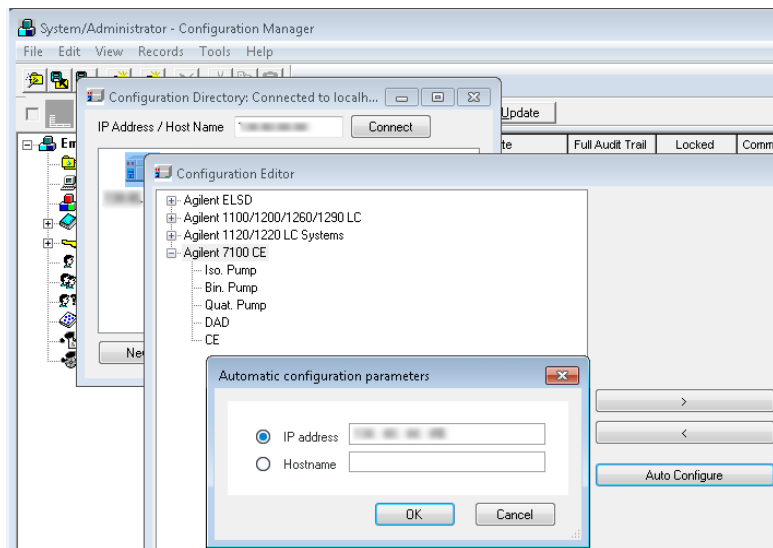


Figure 5 Configuration of the CE

- 4 In the Configuration Editor select the Agilent 7100 CE and click **Auto Configure**.
- 5 Enter the IP address of the instrument and click **OK**. The instrument is detected and the CE and DAD appear on the right-hand side of the configuration window.

Hint

The default IP address is 192.168.254.11. Please refer to the *G7100A User Manual* if an IP address change is required.

- 6 Double-click CE G7150A on the right side of the window or select it and click **Configure** at the bottom of the screen. The CE configuration window opens.

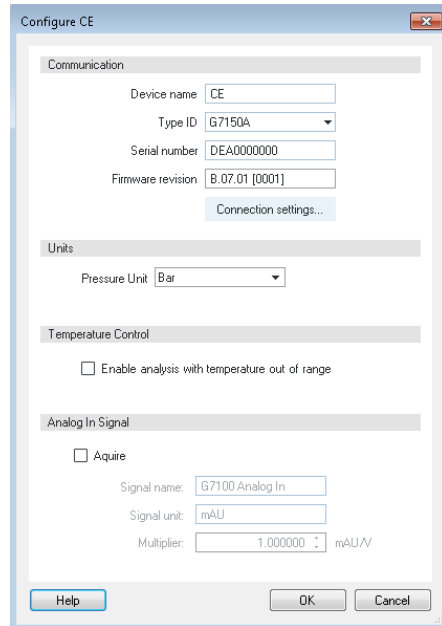


Figure 6 Possible configuration options for CE

- a Select the preferred **Pressure Unit** in the CE status dashboard for use in the data and the pressure auxiliary channel
- b If **Temperature Control** is not required, select **Enable analysis with temperature out of range**.

NOTE

Temperature control is enabled by default. The run only starts once the given temperature is reached. You can change this behavior by deselecting the box. A reboot (and reconfiguration via the PreConfiguration Utility) is required.

- c Enable **Acquire** for the **Analog In Signal** to collect data.

- 7 Click **OK** to close the screen.
- 8 Click **OK** to close the configuration editor.
- 9 Click **OK** to close the PreConfiguration Utility.

Set up and Configuration of CE in Empower 3

Configuration of CE in Empower

Refer to the Waters Empower documentation for installation and configuration of the LC/CE system in Empower.

- 1 Open the **Waters DHCP Server Configuration** window by entering the following command in the windows command line:

C:\Empower\Instruments\Waters DHCP Server Configuration.exe.

OR

In the Empower Configuration Manager window, select **Node** in the tree on the left-hand side, then right-click on the node you want to add the instrument to and select **Properties**. Select the **Configure DHCP** tab and click **Configure DHCP**.

- 2 Add the **IP Address** and **MAC Address** manually for the CE mainframe.
- 3 Select **Instrument Type AgilentLC** and click **OK** to leave the screen.

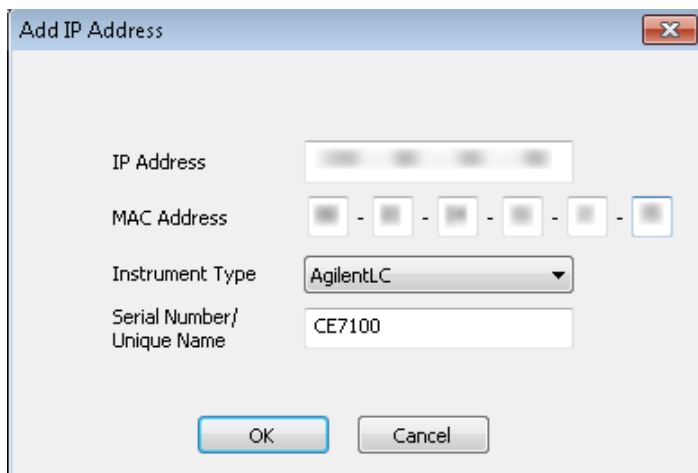


Figure 7 Entering communication details for CE

- 4 Access the **Nodes Properties** in the Empower Configuration Manager and verify that the Instrument is shown as OK.

- 5 Generate a new chromatographic system with the newly configured instrument using **File > New > Chromatographic System**. Follow the instructions on the screen.

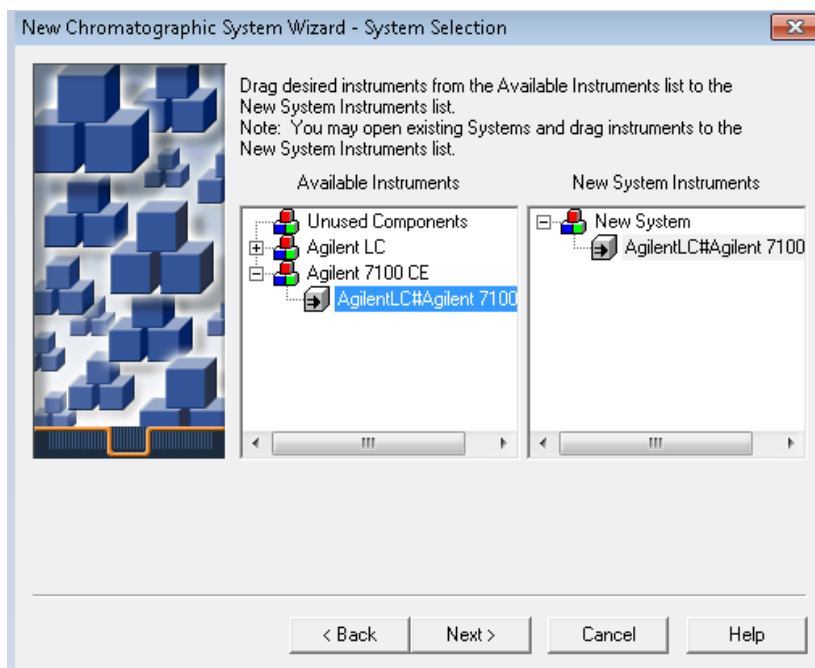


Figure 8 Generating a new chromatographic system

Using the CE in Empower

The CE Status Dashboard in Empower – Direct Control

- 1 Start Empower and open the **Run Samples** Screen.
- 2 Select the CE/CIA specific project for data collection together with the CE system.
- 3 The CE Status window displays all available modules with their status information. Details can be seen by hovering over the status bar with the mouse.

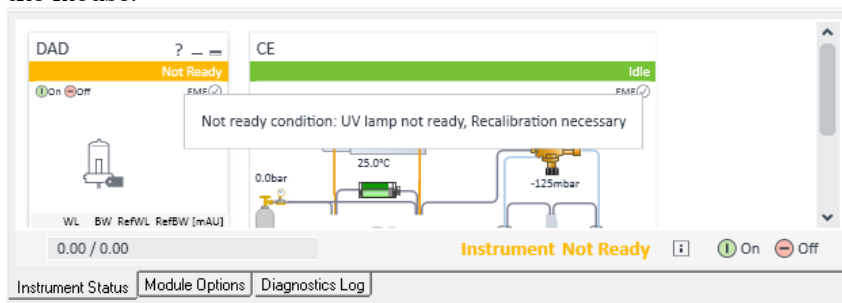


Figure 9 CE dashboard status information

- 4 The given space does not allow enlarging the CE status window to full size, therefore a vertical scrollbar is present to adjust the required section of the status dashboard.

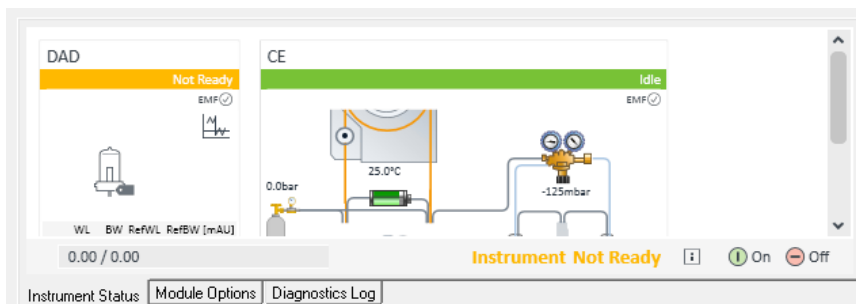


Figure 10 CE Status dashboard – upper part

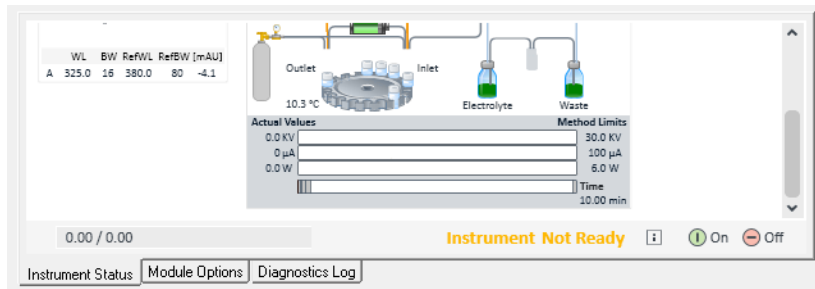


Figure 11 CE Status dashboard – lower part

Each component of the instrument is represented by an icon on the CE status dashboard. Right-click an icon to access direct control. A context menu appears and an action can be performed. Some contextual actions require further user interaction. Click **OK** to trigger the action.

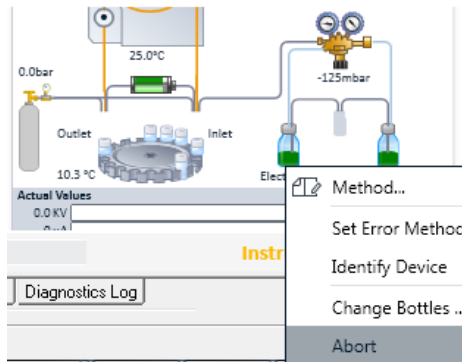


Figure 12 Example of direct control - actions available for bottle icon

Following major actions are accessible via the CE Status dashboard only (list does not claim to be exhaustive):

- Change Cassette
- Switch on lamp/Switch off lamp
- Set Inlet Vial
- Set Replenish Vial
- Unload Replenish Lifter
- Unload Inlet Lifter
- Set Outlet Vial
- Unload Outlet Lifter
- Flushing Capillary by Flush
- Injection by Apply Pressure
- Apply Voltage
- Replenish Vials
- Change Bottles
- Get Vial

All actions are performed immediately and the changes are reflected in the graphical user interface (GUI). The system is in **not ready** state (yellow), while performing these actions. After the action finishes, the instrument resumes an **idle** state (green). Any direct change of a parameter in the direct control menu (control and/or method) does not change the current instrument method.

NOTE

It is not possible to see the detector signal while flushing the capillary. A feasible workaround is to check the numerical value of the detection wavelength, which is displayed in the DAD part of the GUI (left side).

Creating an Instrument Method

The instrument method contains all the parameters necessary to perform the sample acquisition.

- 1 Open the instrument method editor, for example via **Edit > Instrument Method**.
- 2 Click through the various tabs and enter the instrument method parameters.
- 3 Save the instrument method under a dedicated name. This method can now be used as part of a method set.

The instrument method editor offers various tabs:

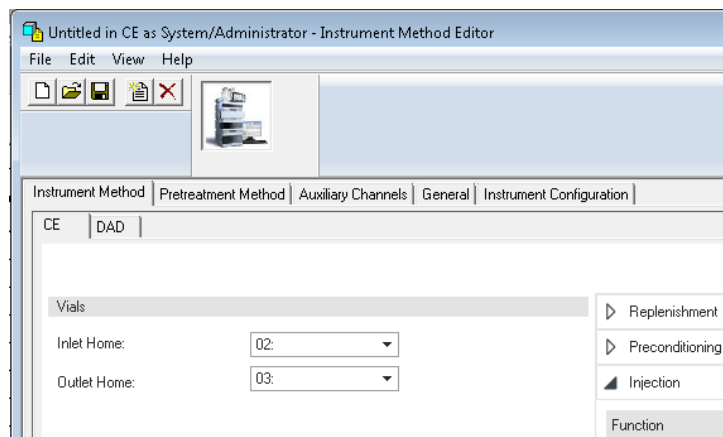


Figure 13 Instrument Method Editor

- Instrument Method tab:** The instrument method tab provides access to all method parameters of the CE system, one tab per module, **CE** and **DAD**. Select the appropriate tab to enter the method parameters. For details on the parameters refer to the online help and the CE or DAD user manual.
- Pretreatment Method tab:** There are no pretreatment steps available for CE.
- Auxiliary Channels tab:** Offers selection of auxiliary traces during a run, e.g. the CE leak current.

- d General** tab: Offers option to mark this method as shutdown method, see page 20. Allows selective shutdown of the CE and DAD.
- e Instrument Configuration** tab: The CE system configuration is available in the **Instrument Configuration** tab. All settings in this dialog must be defined using the PreConfiguration Utility. Changes in this section of the method are not applied to the system; they are saved with the instrument method only. Any changes, such as enabling temperature control, need to be done in the PreConfiguration Utility to apply the new configuration to the physical instrument. In addition, the LAC/E box has to be rebooted and a new autoconfiguration is required in order to adjust the new configuration with the LC Driver and Empower.

Instrument Method tab

The CE method Screen

The CE method screen contains all parameters used to adjust the separation, meaning the home vials can be defined as well as **Cassette Temperature**, **Voltage**, and **Stoptime**. Press F1 on the keyboard to access the online help, offering explanations for each method parameter. The left-hand side of the method screen includes the initial starting parameters.

CE (G7100A)

Vials

Inlet Home: 01

Outlet Home: 02

Cassette Temperature

Not controlled

25.0 °C

High Voltage System

Enable High Voltage

Voltage: 0.0 KV positive Polarity

Current: 300 µA

Power: 6.0 W

Low Current alarm limit: Off

100 µA

Stoptime

As Injector/No Limit

4.00 min

Off

1.00 min

Replenishment

Preconditioning

Injection

Timetable (1/100 events)

Time [min]	Function	Parameter
0.30	Change Voltage	30 KV

Add Remove Clear all

Cut Copy Paste

Postconditioning

Figure 14 CE method screen

NOTE

Enter a **Stoptime** for the CE and/or the DAD. For run execution a Run Time is specified in Empower later, ensure that both values are synchronized. For more details, see section Executing Runs on page 21.

Using the CE in Empower

Hint

Depending on the used SOP, the following best practices are used.

- a Ensure that Empower **Run Time** and Agilent Module **Stoptime** (for one module) are synchronized.
- b Set an Agilent Module **Stoptime** of 1 min, and set the real sample time in the sample set screen.

If an entered parameter is out of range, a warning sign is shown. By hovering with the mouse pointer over the parameter, the message displays the possible range, e.g. **Voltage is out of range [-30kV to 30 kV]**.

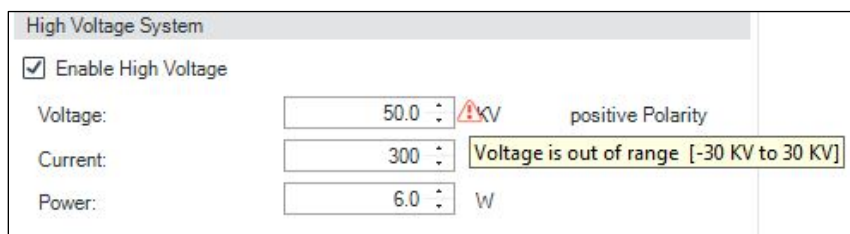


Figure 15 Out of range handling

The right-hand side of the method screen offers additional features. The online help explains each of the features:

- Replenishment
- Preconditioning
- Injection
- Timetable
- Postconditioning

The DAD method screen

On the left side, the DAD method offers the main method parameters to acquire up to eight signals. Refer to the online help to learn more about each parameter. Remember to set the stop time either to CE or to the same time as the Empower run time.

The right side offers additional method parameters, e.g. for spectra collection and a timetable, in order to change parameters during a run.

Unlike with other Agilent DAD, there is no VIS-lamp available.

The screenshot shows the 'DAD (G7151A)' method configuration screen. The 'Signals' section contains the following table:

Signal	Acquire	Wave length	Band width	Reference Wavelength	Reference Bandwidth
Signal A	<input checked="" type="checkbox"/>	200.0	5.0	360.0	100.0 nm
Signal B	<input checked="" type="checkbox"/>	210.0	4.0	360.0	100.0 nm
Signal C	<input checked="" type="checkbox"/>	214.0	4.0	360.0	100.0 nm
Signal D	<input checked="" type="checkbox"/>	230.0	4.0	360.0	100.0 nm
Signal E	<input checked="" type="checkbox"/>	260.0	4.0	360.0	100.0 nm
Signal F	<input type="checkbox"/>	273.0	4.0	360.0	100.0 nm
Signal G	<input type="checkbox"/>	280.0	4.0	360.0	100.0 nm
Signal H	<input type="checkbox"/>	250.0	100.0	360.0	100.0 nm

The 'Peakwidth' section is set to '> 0.025 min (0.5 s response time) (10 Hz)'. The 'Stoptime' section has 'As CE' selected. The 'Posttime' section has 'Off' selected. The 'Advanced' section includes 'Spectrum' (Store: All, Range from: 190.0 to 400.0 nm, Step: 2.0 nm), 'Analog Output' (Zero Offset: 5%, Attenuation: 1000 mAU), 'Indirect UV - Margin for negative Absorbance' (100 mAU), and 'Autobalance' (Prerun and UV Lamp checked, Postrun unchecked).

Figure 16 DAD method screen

General Method Tab

Empower offers the option of selecting a **System Shutdown Method**, which is executed at the end of the sequence.

- 1 To generate a shutdown method, define the desired shutdown conditions, such as vials and times in the CE/DAD method screen first and select **Shutdown after Run**. It is possible to perform a selective shutdown, for example only the DAD.
- 2 To execute the shutdown method, either select the shutdown method in the Run Sample set window or load the shutdown method as the last line of the sequence.

The screenshot displays the 'Instrument Configuration' tab with the following settings:




- Fraction Collector Options:** Start at Location: 1
- System Shutdown Method:**
 - Shutdown after Run  Do not use this Instrument Method in a Sample Set for your regular analysis. Instead, specify this method as the Shutdown method on the Run Samples Defaults screen found under the Customize/Defaults menu selection of Run Samples or QuickStart.
 -  Default behavior: the system will attain initial conditions and run until the Stoptime. At the Stoptime, pump flow will go to zero, column temperature will be turned off, detector lamps will be turned off, and autosampler temperature will remain unchanged.
- Selective Shutdown:**
 -  Selecting any of the modules below overrides the default shutdown behavior and ignores Stoptime. The selected modules will turn off after initial conditions are met. Unselected modules will remain on at the initial conditions specified in this Method.
 - CE (CE0) - Off
 - DAD (DAD0) - Off

Figure 17 Shutdown Method screen

Executing Runs

NOTE

The Empower **Run Time** is shown in the **Single Run** tab and in the sample sets. The Agilent **Stop Time** is an instrument method parameter.

- If the **Run Time** is *longer* than the instrument method's **Stop Time**, the **Run Time** is used for the measurements, regardless of what is stated in the instrument method. The online plot and the resulting electropherogram end with the **Run Time**.
- If the **Run Time** is *shorter* than the instrument method's **Stop Time**, the online-plot stops at the **Run Time** but the run continues until the end of the method's **Stop Time**. The resulting electropherogram ends at the method's **Stop Time**. The longer time is used for recording data.

The method's **Post Time** is independently added after the **Stop Time**, and no data is recorded during the post time. The next sequence run starts after the **Post Time** of the previous run is finished.

Single Runs

- 1 Enter or select the required values and parameters in the **Single** tab for a single run execution. Click the **Inject** icon to start the single run.
- 2 During the run execution, you can stop the run at any time by clicking the red Stop button.

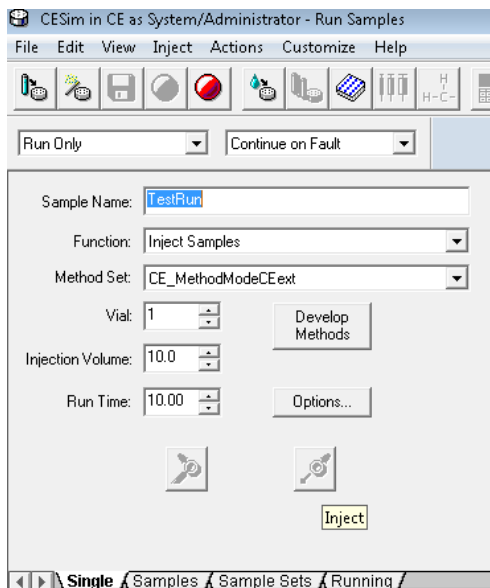


Figure 18 Single Run tab

Sequence Execution/Sample sets

Empower offers multiple injections from the same vial as well as consecutive injections from different injection vials. The details have to be added to the sample/sample set table (sequence), depending on the required workflow.

- 1 Click the **Sample** tab to open the table.
- 2 Click into the first line to fill in the appropriate values. For the following values the entries are mandatory:
 - Vial position
 - Injection volume
 - Number of injections
 - Sample name
 - Method set
 - Run time (use the same time as in the instrument method)

NOTE

The parameter **Inj. Vol (µl)** requires a value. This value is *not* used for the injection, but it is reported. The instrument method parameters (for example: pressure and time) are used for the injection.

(Instead of volume the quantifiable parameters are pressure × time for hydrodynamic injection, or voltage × time for electrokinetic injection).

Entering an injection volume of 0 µl does not result in a blank run. The value is ignored for the run execution, but it is reported.

-
- 3 Start the sequence. Reminder: During run execution, the red Stop button allows you to stop the run at any time.

NOTE

If a vial is missing, the current run is stopped and the whole sequence is aborted. The option **Next sample on Fault** does not apply.

- 4 It is possible to shut down the CE/DAD at the end of the sample set. Load the shutdown method (for details see section General Method Tab on page 20) as the last line of the sequence.

Replenishment and Conditioning

Replenishment allows to automatically change the buffer of a vial either in serial or parallel mode.

Preconditioning and **Postconditioning** offers conditioning of the capillary.

These settings are method settings and therefore expected each time the method is running. If this is not wanted, make sure to generate methods with and without replenishment and/or pre/postconditioning.

During multiple analyses the buffer must be refreshed after a certain number of runs. Typically, the exchange of the buffer is done every 3 - 10 runs. Depending on the stability of the buffer, it is sometimes necessary to refresh it before each run.

NOTE

Using the replenishment as a part of the method, the replenishment is carried out for each run where this is part of the method.

If you want the refresh to be done less, ensure to create at least two methods.

Example:

Run 1 method A with replenishment
Run 2 method
Run 3 method
Run 4 method with replenishment
Run 5 method
Run 6 method
Run 7 method with replenishment
Run 8 method
Run 9 method

The replenishment system provides a quick way to change the buffer automatically. The system removes the used buffer from the vials and transfers it into the waste bottle. Then the vials are filled with fresh buffer from the electrolyte reservoir.

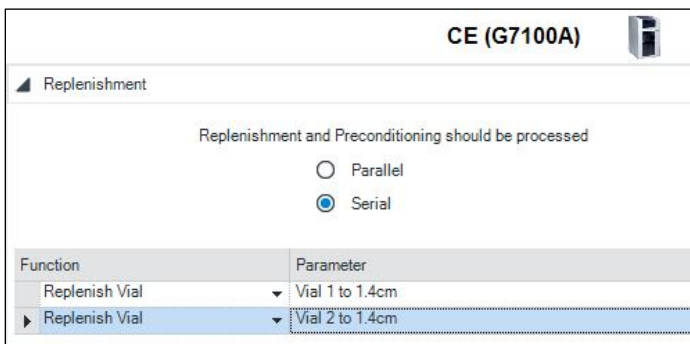
NOTE

Refer to the G7100A User manual on the requirements for the bottles, as only specific bottles can be used. For example, they need to be pressure-stable.

The waste bottle should be 500 ml, whereas the electrolyte bottle can also be a 100 ml bottle. Both bottles are included in the accessory kit.

Following replenishment, tasks can be set up in the CE replenishment table in the Instrument method:

- **Empty vial:** empties a vial into the waste bottle
- **Fill vial:** fills a vial to a user-selectable level from the electrolyte reservoir
- **Clean tubes:** flushes the replenishment system to clean the tubes
- **Replenish vial:** empties a vial into the waste and fills it up from the electrolyte reservoir
- **Wait:** waits for a specified time



CE (G7100A)

Replenishment

Replenishment and Preconditioning should be processed

Parallel

Serial

Function	Parameter
Replenish Vial	Vial 1 to 1.4cm
Replenish Vial	Vial 2 to 1.4cm

Figure 19 Replenishment Table

It is possible to execute replenishment in parallel, which decreases the overall analysis time as replenishment and pre- or postconditioning are executed simultaneously.

NOTE

For parallel replenishment, the vials for the first run must be filled with the correct solutions, because the replenishment starts at the end of the run.

NOTE

To efficiently clean the replenishment system after use, the special functions **Flush Tubes** and **Clean Level Sensor** should be performed. This advanced cleaning is only available in the Agilent LabAdvisor software.

Indirect detection

The indirect detection mode allows the application of indirect photometric detection of non-UV absorbing analytes like inorganic cations, inorganic anions and small carbohydrates.

To attain indirect photometric detection, an ionic compound with a high UV absorbance intensity is used as the background electrolyte (BGE). During the separation any non-absorbing analyte causes a reduction of the high background signal, resulting in negative peaks.

Empower can record and calculate positive and negative signals. If most of the analytes are non-absorbing, the detector signals for signal and reference wavelength can be inverted, so that most of the peaks are positive.

Application of external pressure

For some special applications using high viscosity buffer solution (e.g. a gel matrix in the CGE and cIEF) the capillary can be conditioned using external high pressure to ensure that the capillary is filled in adequate time.

Two additional modes are available for configuration:

- **CEC:** Capillary Electrochromatography allows high pressure to be applied on both electrolyte vials in the run during the high-voltage application to prevent outgassing and bubble formation.

A connected external pressure supply is required for this mode. When this mode is selected, additional pressure options are available in the instrument method:

- High pressure flush in **Preconditioning, Injection** and **Postconditioning**.
- External pressure as method setpoint and timetable entry.
- **CE/p:** additional external pressure connected

Configure CEC or CE/p mode

If an external pressure source is available (pressure 2 - 12 bar) and connected to the CE device, these modes can be configured in the **Module Options** tab, which can be found below the CE status dashboard. Select the required CE Mode.

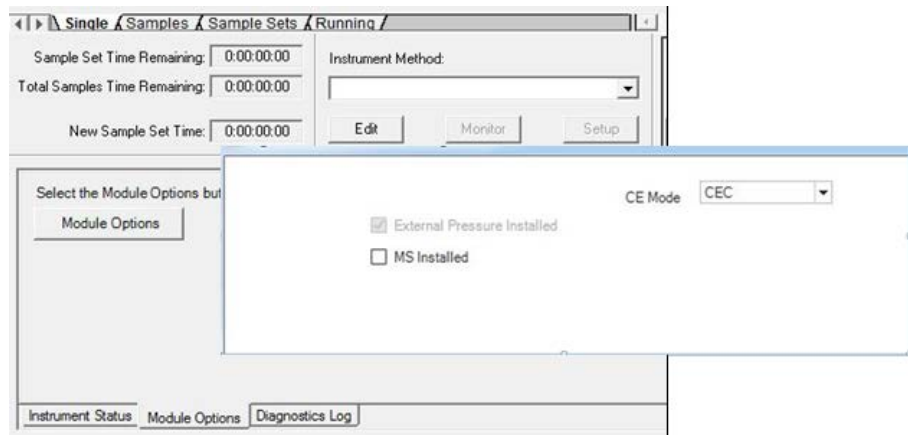


Figure 20 Additional CE mode

When the **External Pressure Installed** check box is selected, the gas cylinder is shown in the CE status dashboard accessible via the **Instrument Status** tab. High pressure can now be applied by right-clicking on the gas cylinder.

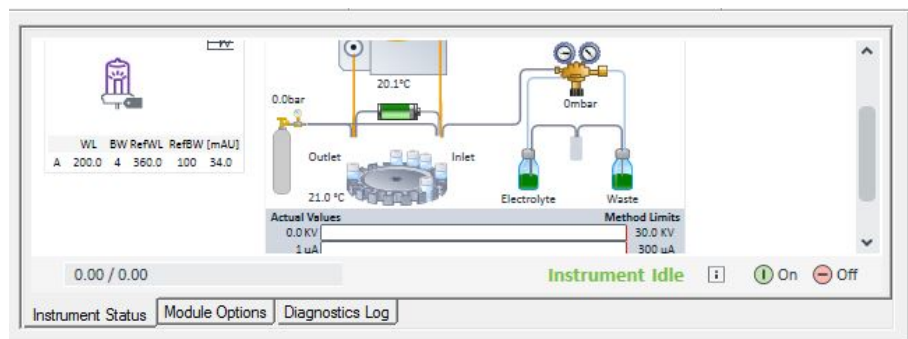


Figure 21 Access to external pressure

Using the CE in Empower

NOTE

Make sure to reboot the LAC/E box after any mode change in the Module Option tab. Without a reboot (*OR* close Empower and end the processes **InstrumentServer.exe** and **AgilentPlugInServer.exe** and reopen Empower) the configuration changes are not correctly applied and can result in various issues.

- If the PreConfiguration Utility was used, it is essential to create a new autoconfiguration.
 - Verify that other instruments attached to the Workstation or LAC/E are not collecting data. Rebooting the LAC/E will interrupt data collection on all instruments attached and the lost data will not be able to be recovered.
-

Method impact for CEC or CE/p mode**NOTE**

If the new pressure section does not appear in the method section, the configuration mode has not been applied correctly. Confirm that the correct mode is defined in the **Module Option** tab and then reboot the LAC/E. If a PreConfiguration exists, run a new autoconfiguration in addition.

- If **CE/p** is selected, the instrument method offers high pressure settings as additional method parameters on the right-hand CE method screen:
 - Preconditioning (right-hand CE method screen)
 - Injection (right-hand CE method screen)
 - Timetable (right-hand CE method screen)
 - Postconditioning (right-hand CE method screen)
- If **CEC** is selected, the instrument method offers high pressure settings as additional method parameters on the right-hand CE method screen:
 - Preconditioning (right-hand CE method screen)
 - Injection (right-hand CE method screen)
 - Timetable (right-hand CE method screen)
 - Postconditioning (right-hand CE method screen)

Only in CEC mode, the instrument offers an additional initial value **Pressure** on the left-hand CE method screen. Selecting **External**, the pressure is applied to both home vials during run execution (see Figure 23).

Using the CE in Empower

Instrument Method | Pretreatment Method | Auxiliary Channels | General | Instrument Configur

CE | DAD

Vials

Inlet Home: 01

Outlet Home: 02

Cassette Temperature

Not controlled

25.0 °C

High Voltage System

Enable High Voltage

Voltage: 0.0 KV positive Polarity

Current: 300 µA

Power: 6.0 W

Low Current alarm limit: Off 100 µA

Pressure

Off

Internal 0 mbar

Raw External Pressure

External 2.0 bar

Stop time | Post time

As Injector/No Limit Off

5.00 min 1.00 min

Figure 22 Applying external pressure – new menu

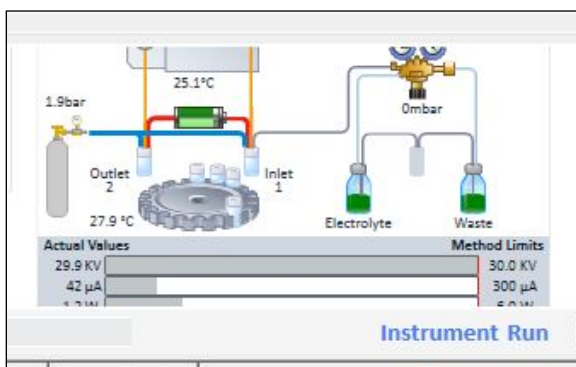


Figure 23 Applying external pressure during a run

Empower features for CE Instrument Control

Empower 3 offers all key functions to control the G7100A Agilent Capillary Electrophoresis system. All established techniques are feasible, such as the use of direct and indirect detection modes, use of special electrolyte systems (containing micelles or gel matrices) and also capillary isoelectric focusing (cIEF) or capillary electrochromatography (CEC). The application of external high pressure is supported.

For data analysis, Empower offers processing tools for peak integration and calculation of concentrations in CE specific mode.

Known Limitations/Not tested configurations:

- **Pressure Unit** configuration
Selecting PSI as pressure unit, the CE status dashboard shows the pressure in PSI, while the online plot and the resulting Aux Traces on the Chromatogram offer the values in mbar/bar. The correct pressure is applied.
- **Analog In**
The feature **Analog In** has not been explicitly tested, but the functionality is present.
- **MS installed**
The feature MS installed has not been tested.



Figure 24 MS installed

Extended CE driver features available in Agilent OpenLAB CDS ChemStation Edition

The following items are not CE driver features, hence they are not available in Empower environment.

Control and Action Menus

In Non-Agilent CDS all actions and direct controls can only be accessed via the CE Status dashboard. There are no menu items for instrument control in the Empower menu bar.

Fraction Collection

The system offers **Set up for Peak Fraction Collection** whether a fraction collector is present or not. This Macro is only available in OpenLAB CDS ChemStation Edition.

User vials

The CE driver offers to select user vials (1-10) as part of the method. User vial parameters are defined in the sequence table as the hosting CDS interlinks method and sequence. User vials are not available in Non-Agilent CDS environment.

Capillary Catalog, Capillary Handling

OpenLAB CDS ChemStation Edition offers a capillary catalog database. This database is not available in Non-Agilent CDS environment.
Workaround: Add the capillary information to the **Method Comments**.

Sample Diagram

OpenLAB CDS ChemStation Edition offers a sample diagram informing about each vial's purpose. This table is not available in Non-Agilent CDS environment.

Vial Table

OpenLAB CDS ChemStation Edition offers a vial table, giving an overview of vial position, name and comment.

Calibration Curve options

The following Calibration Curves are unavailable in Empower:

- Mobility correction
- Calibration Type cIEF
- Calibration for determination of isoelectric points or molecular weights

Additional Information

Waters

Additional information on the PreConfiguration Utility is provided on the Waters Support Webpage for registered users:

<http://www.waters.com/waters/support.htm?lid=134936402&cid=511442&type=TECN>

Document reference: TECN134936402

Title: Using the Agilent PreConfiguration Utility with Agilent ICF Support Version 2.2

APPENDIX - Examples

Example: Single test run

For testing the functions of the 7100 CE and Empower 3 a test method can be created. The equipment and method parameters can be summarized as follows:

Capillary:	Fused silica capillary 50 µm ID, 48.5 cm total (40 cm effective), ext. Light path (G1600-60232)
Electrolyte:	20 mM Borate (from IQ-Kit: 5063-6514)
Home vials:	1 and 2 (filled with electrolyte solution)
Sample:	4-Hydroxyacetophenone solution 1 mmol/l (from IQ-Kit: 5063-6514)

NOTE

The sample has to be diluted 1:10 before the run.

Voltage:	+30 kV
Temperature:	25°C
Stop Time:	4 min
Preconditioning:	300 s (from electrolyte vial 3 into waste vial)
Injection:	50 mbar for 10 s (from injection vial to outlet vial)
Post-injection:	50 mbar for 5 s (from inlet vial to outlet vial)
DAD:	Signal: 200 nm (BW 5 nm) Reference: 360 nm (BW 100 nm) Peakwidth: >0.025 min (10 Hz)
DAD Timetable:	0.5 min Balance

The run can be executed as single run (in this case save the instrument method as well as the method set). If defined in the method, the online plot shows the selected signals and the final result electropherogram can be obtained in the **Browse** menu of Empower or via the projects.

NOTE

Ensure that the Run Time is the same as defined in the CE/DAD method.

APPENDIX - Examples

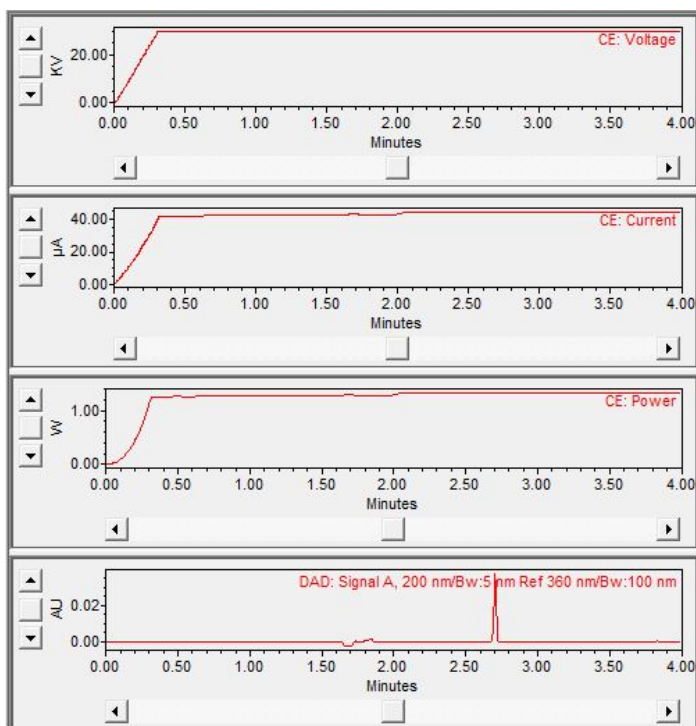


Figure 25 Online plot during the run

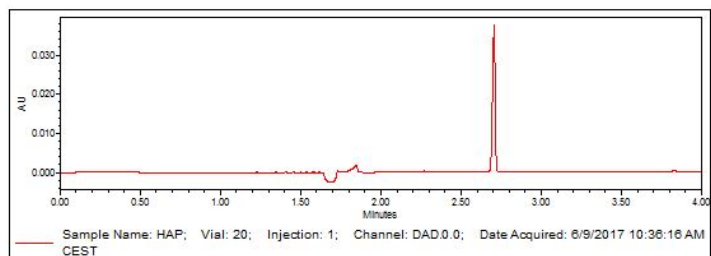


Figure 26 Resulting electropherogram

Example: Multiple injections using the IQ Kit for CE

For testing multiple injections in one sequence, the instrument method and the experimental setup described in section Example: Single test run can be used. See the sequence table (Figure 27) and the overlay of the resulting electropherograms (Figure 28).

A set of six injections was used to calculate system precision. The peak areas and relative standard deviations (Excel: STDEV) for the areas, migration times and the corrected areas are summarized in the following table.

Table 2 Peak areas and relative standard deviations

	Area	t_m [min]	t_m [s]	corr. Area (area/ t_m [s])
Injection 1	50049	2.730	163.80	3.05549E+02
Injection 2	50547	2.723	163.38	3.09383E+02
Injection 3	50189	2.717	163.02	3.07870E+02
Injection 4	49767	2.711	162.66	3.05957E+02
Injection 5	49187	2.706	162.36	3.02950E+02
Injection 6	49579	2.702	162.12	3.05817E+02
Average	49886	2.715	162.89	3.06254E+02
Deviation STDEV:	0.96 %	0.39 %	0.39 %	0.72 %

Vial	Inj Vol (uL)	# of Injs	Label	SampleName	Level	Sample Matrix	Function	Method Set / Report Method	Label Reference	Processing	Run Time (Minutes)	Data Start (Minutes)	Next In Delay (Minute)
11	20	10.0	1	Blank			Inject Samples	20170313		Normal	5.00	0.00	0.1
2	21	10.0	6	4-HAP, 1:10			Inject Samples	20170313		Normal	5.00	0.00	0.1
3	20	10.0	1	Blank			Inject Samples	20170313b		Normal	5.00	0.00	0.1

Sixfold sample injection

Figure 27 Sequence table for multiple injections

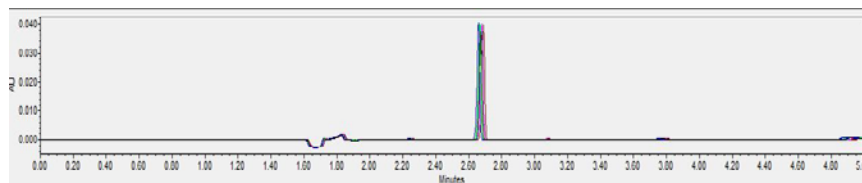


Figure 28 **Overlay of six consecutive injections**

Using Empower 3 in combination with the Agilent 7100 CE, a good relative standard deviation around 1 % for the peak areas and below 0.5 % for the migration times were obtained. These satisfactory values are achieved by a special self-regulating injection procedure, which significantly improves the injection precision. For more information please refer to the *Agilent CE Manual*.

Example: Indirect detection using the Plating Bath Buffer

A test method can be created using the *Plating Bath Buffer* for testing the indirect detection functions. The following reagents and parameters can be used:

Capillary:	Fused silica capillary 50 μm ID, 80.5 cm total (72 cm effective) (G1600-62211)
Electrolyte:	Plating Bath Buffer (Agilent: 5064-8236)
Home vials:	1 and 2 (filled with electrolyte solution)
Sample:	Suitable <i>Plating Bath Mixture</i> (laboratory made)
Voltage:	-30 kV
Temperature:	25°C
Stop time:	15 min
Preconditioning:	240 s (from electrolyte vial 3 into waste vial)
Injection:	1 Injection vial to outlet vial: 50 mbar for 10 s 2 Post-injection: 50 mbar for 5 s (from inlet vial to outlet vial)
DAD:	Signal A: 350 nm (BW 20 nm), reference: 275 nm (BW 10 nm) Signal B: 275 nm (BW 10 nm), reference: 350 nm (BW 20 nm) Peak width: >0.025 min (10 Hz) Timetable: 1.0 min Balance

Comparing the reversed wavelength the peaks can be shown as positive or negative peaks.

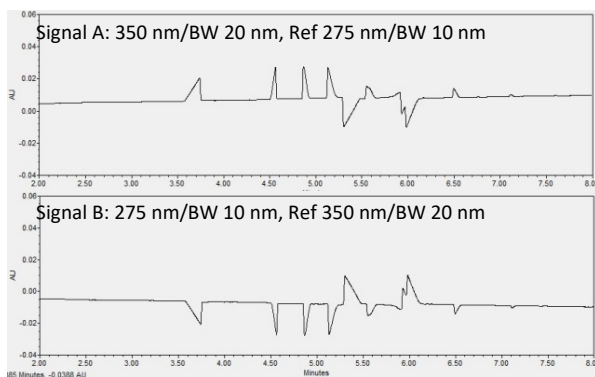


Figure 29 Indirect detection

It is evident from the electropherogram in Figure 30 that the selected wavelength values result in positive peaks for the non-UV absorbing analytes: sulfate, malate, hypophosphite, phosphite and lactate. The nickel-complex has its own UV intensity and results in a negative peak. The figure also shows that the integration and evaluation of the peak areas are possible for positive and negative peaks in the same signal.

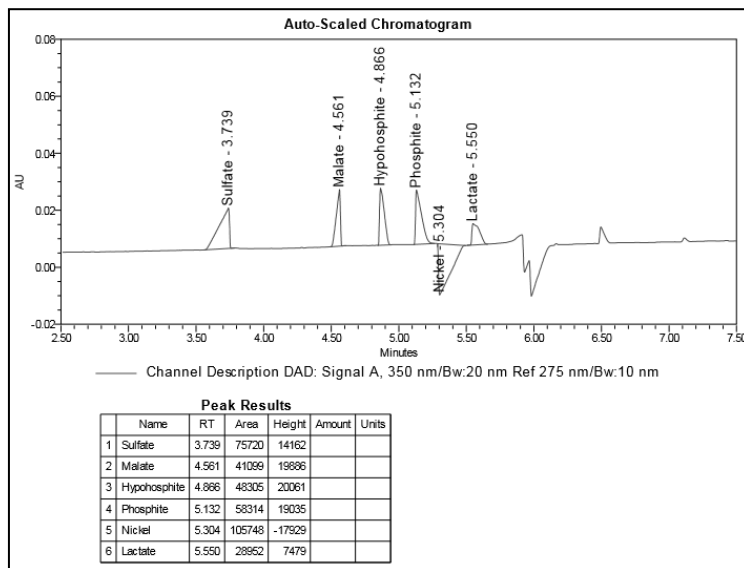


Figure 30 Integration and reporting of positive and negative peaks



G7100-90200

Edition: 01/2018

Edited in Germany

© Agilent Technologies, Inc 2018

Agilent Technologies, Inc
 Hewlett-Packard-Strasse 8
 76337 Waldbronn
 Germany