

InfinityLab Pro iQ Series

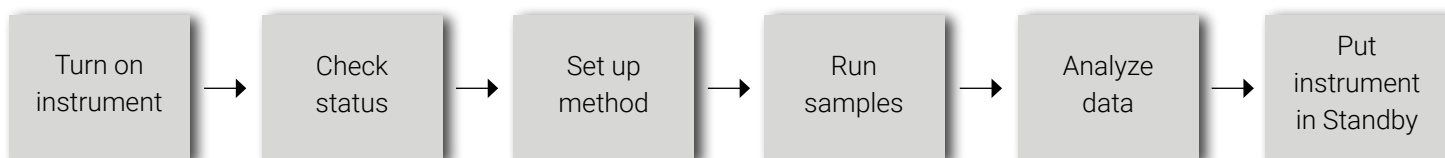
This Quick Reference Guide contains instructions for using the Agilent InfinityLab Pro iQ Series single quadrupole instruments with OpenLab CDS acquisition and data analysis software.

See the SQ Driver online Help for detailed instructions on each workflow step shown below.



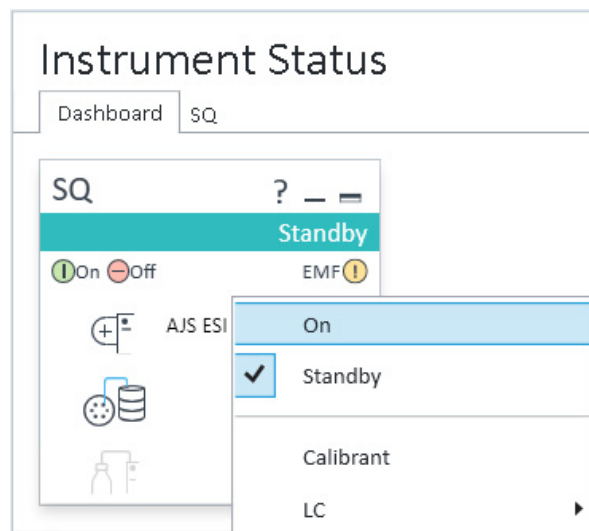
Workflow Overview

The following figure shows the daily LC/MS workflow when using the OpenLab CDS software.



Turning on the LC/MS instrument

In the **Instrument Status** window, right-click the **SQ** device, and select **On**.



Checking instrument status

1. In the **Instrument Status** window, click the **SQ** tab.
2. Scroll through the list of values to monitor the health of the SQ.

Instrument Status	
Dashboard SQ	
Parameter	Actual
Connection State	Online
Error State	No error
Instrument State	Background acquisition
Run State	Idle
Ready State	Ready
Ready Type	
Not Ready Text	
Error Text	
Run Time(min)	0.00
Calibrant	Off
Quantify Value	To MS

3. Click the **Dashboard** tab.
4. Hover over the EMF icon to see any Early Maintenance Feedback (EMF) messages for the SQ.

Instrument Status

Dashboard **SQ**

SQ ? _ =

Idle

On Off EMF

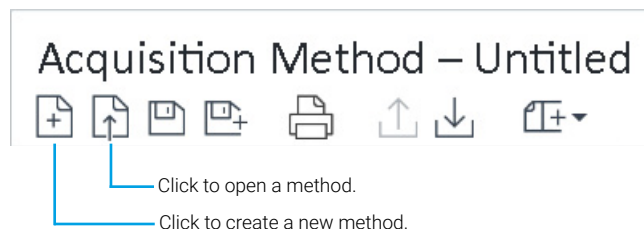
AJS ESI

Checktune expired; Autotune expired.

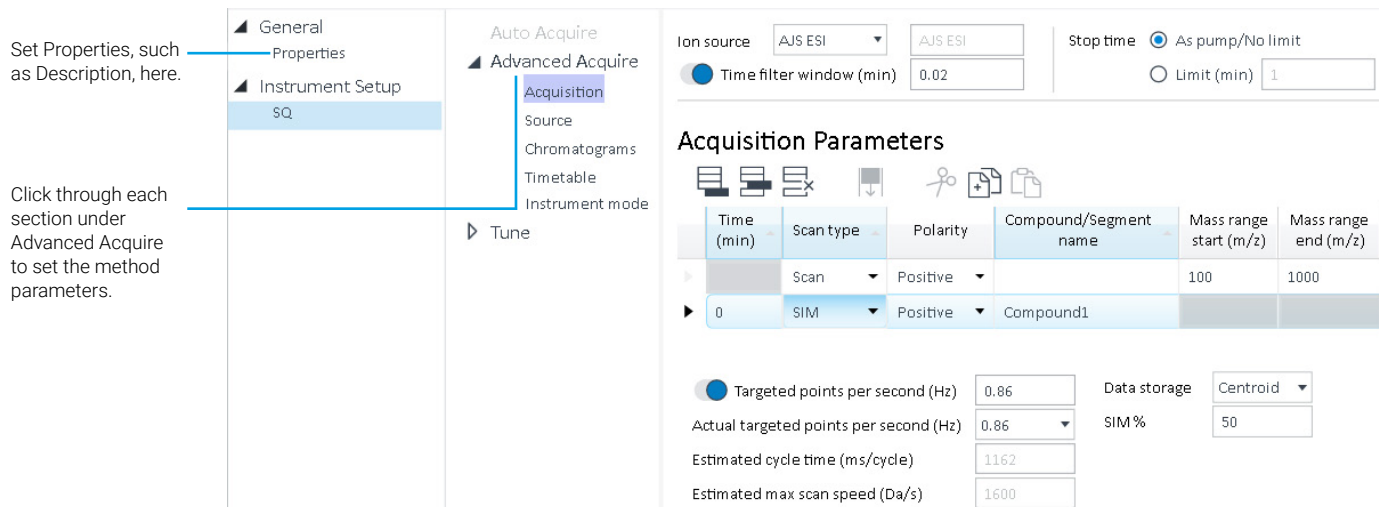
5. Right-click the **SQ** device, and select **Maintenance**. See the LC/MS SQ Driver online Help for detailed instructions on the Maintenance dialog.

Setting up a method

1. On the **Method** layout, create a new method or open an existing method.



2. Edit a method to configure method parameters and instrument configuration settings.



Running samples

Acquire data by running a single sample or running a sequence of samples.

Running a single sample

1. Select the **Single Sample** layout.
2. Complete or edit each section as applicable.
3. If an error is present in one or more parameters, those parameters will be highlighted in red. Hover over a highlighted box to display the error information, or look to the bottom of the Single Sample Analysis window for a description of any errors.
4. Click **Run**.

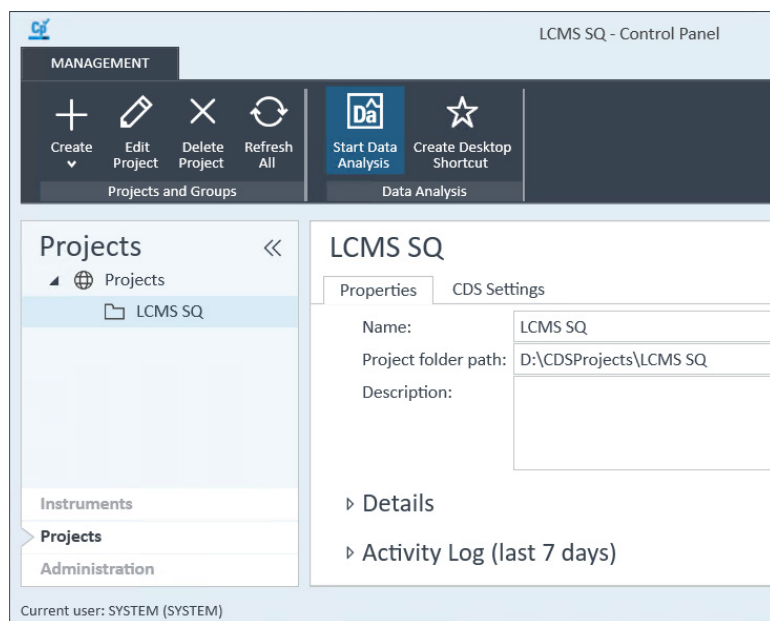
Running a sequence

1. Select the **Sequence** layout.
2. Complete the sequence table. You can:
 - Reuse and modify a saved sequence
 - Create a sequence manually
 - Create a Dual Simultaneous sequence
 - Apply a sequence template
 - Import a CSV or TXT file
3. Select the rows to submit as part of the run. By default, all rows are selected.
4. Click **Validate** to validate the sequence against the specified acquisition method. If invalid entries are found, correct those entries.
5. Click **Run**.

Analyzing data with OpenLab CDS Data Analysis

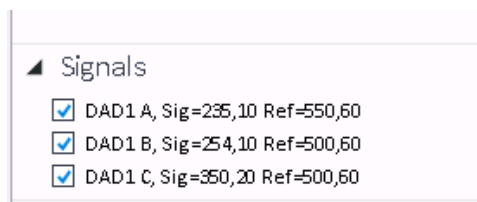
Use Data Analysis to review the sequences and runs completed for your LC/MS project.

1. In OpenLab CDS Control Panel, click **Projects** and select your LC/MS project.
2. Click **Start Data Analysis** in the ribbon. The Data Analysis program opens.



3. In the **Data Selection** view, select your completed sequence result set or single sample and load the associated data.
4. In the navigation pane, select or clear the check boxes for signals in the Signal Selector.

Refer to the online Help for the Data Analysis program for detailed instructions on how to analyze your data.



Putting your instrument in Standby

In the **Instrument Status** window, right-click the **SQ** device, and select **Standby**.