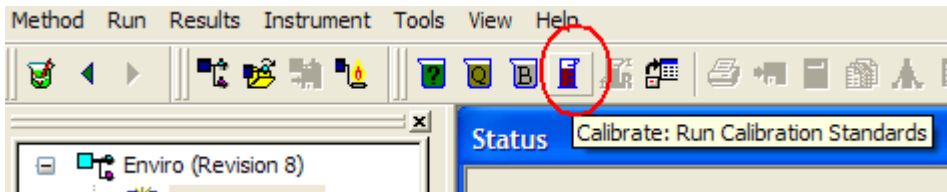


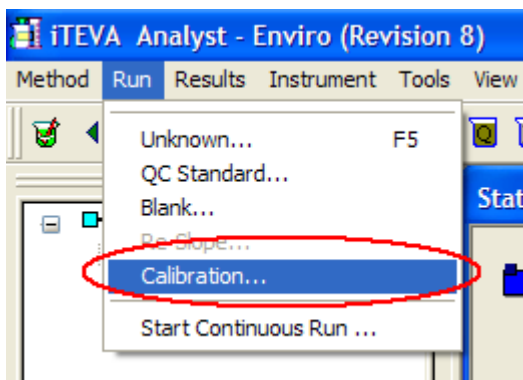
Quick Guides for the iCAP 6000 Series

How to run a manual calibration and unknowns, QCs in iTEVA

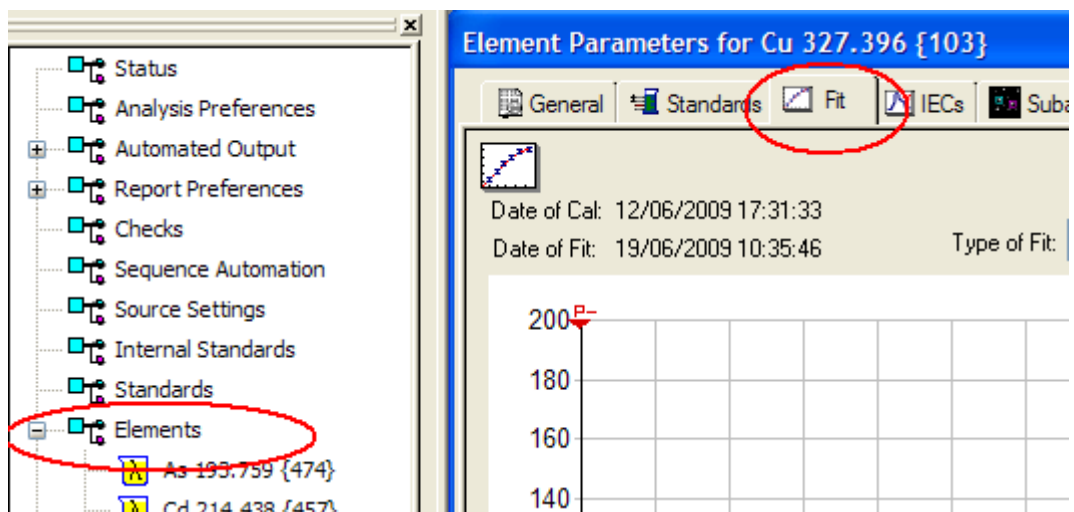
- 1 Open iTEVA and log on
- 2 Single click on *Analyst*, then open the method you require from the list by pointing the cursor on the method name, and then clicking *OK*
- 3 To calibrate the method, either use the *Calibrate* icon on the toolbar



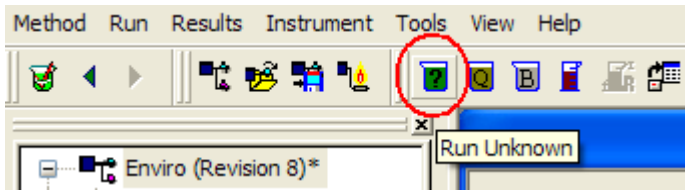
Or the menu option *Run/Calibration*



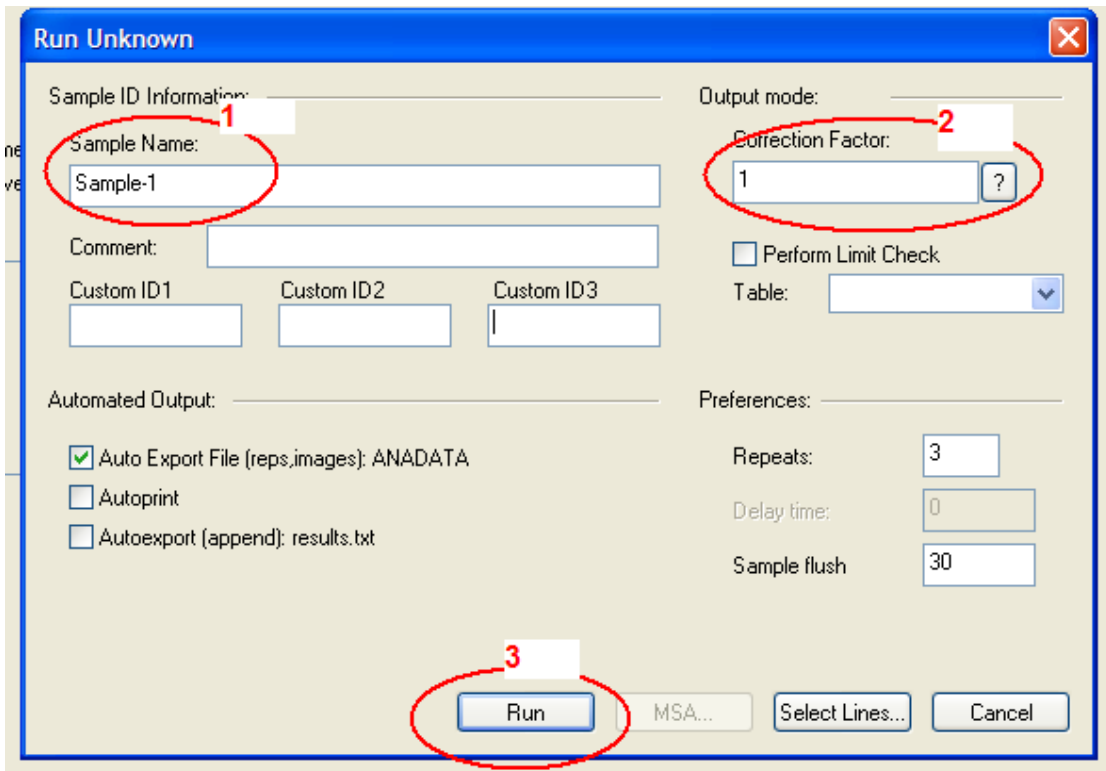
- 4 After the calibration is complete, the method will be updated with the new calibration curve. This can be viewed from *Method/Elements/Fit*



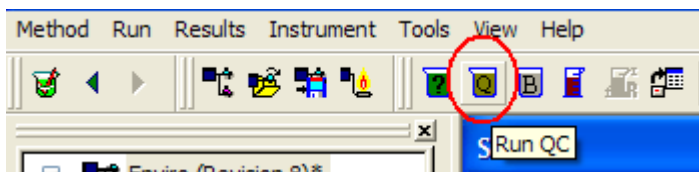
- 5 To run a sample, use the *Run Unknown* icon on the toolbar, or the menu option *Run/Unknown*.



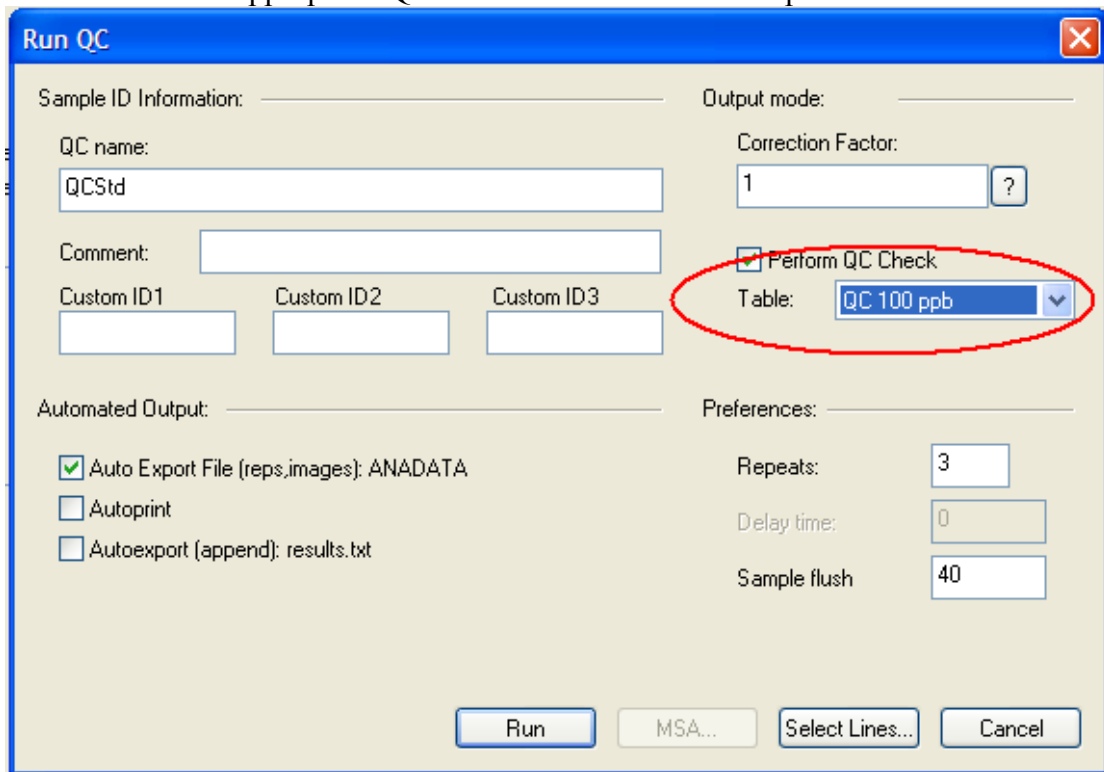
- 6 Enter sample name and any dilution factors, and click *Run*/press enter.



- 7 To run a QC, use the *Run QC* icon on the toolbar, or the menu option *Run/QC*.



8 Select the appropriate QC Check table and click *Run*/press enter.



9 If you need to print the results, use the *Print Sample* icon, to print the selected sample



or the menu option *Results/Print* to print multiple samples.

