

Errata Notice

This document contains references to PSS or Polymer Standards Service. Please note that PSS is now Agilent. This document will be republished as an Agilent document in the future.



**WINGPC
Applications**

Application # 16

POB 3368
Phone: +49-6131-96239-0
D-55023 Mainz
www.polymer.de

**POLYMER
STANDARDS
SERVICE GmbH**



**WINGPC
Newsletter
WINGPC6.2**

applicable for: PSS WINGPC Version 6

Automatic Recalibration with PSS WINGPC

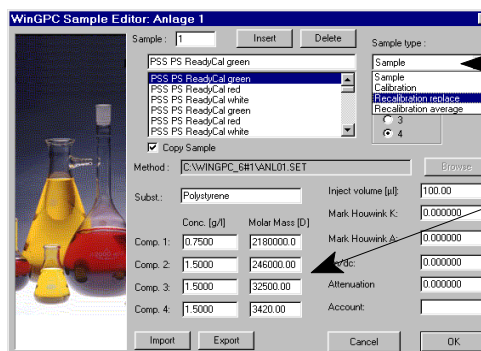
WINGPC 6 can automatically process samples, establish a new calibration curve or recalibrate an existing calibration file.

The sample type selection (*Raw data window*: [Editor] [Samples]) defines how WINGPC processes the actual sample:

Sample type	WINGPC
<i>Sample</i>	Default Automatic processing of the sample (see Appl. #14).
<i>Calibration</i>	WINGPC searches for the peak maximum and adds the calibration point to the calibration table of the calibration file selected in the WINGPC Automation Settings (see Appl. #15).
<i>Recalibration replace</i>	WINGPC searches for the peak maximum and compares the sample name with the names in the calibration table of the calibration file selected in the WINGPC Automation Settings. The entry of the sample with identical name is replaced by the new one.
<i>Rekalibration average</i>	WINGPC searches for the peak maximum and compares the sample name with the names in the calibration table of the calibration file chosen in the automation settings. WINGPC calculates the arithmetic average of the elution volumes of samples with identical sample name.

Automatic Recalibration

1. Enter the sample names and molecular weights in the sample editor (*Raw data window*: [Editor] [Samples]). If you are using more than one component per sample, the molecular weights have to be entered in descending order: component 1 highest, component 4 lowest molecular weight.



Sample type

Molecular weights of all components (decreasing order)

Note: The name of the new sample and that of the sample in the already existing calibration table have to be the same. The best way to make sure that the sample names are identical is to work with the import/export option in the sample editor.

select sample type "Recalibration replace" or "Recalibration average" for all recalibration samples.

2. Open the dialog *WINGPC Automation Settings* in the *method window* ([Definition] [Automation properties]).

Load the calibration file for recalibration using the *Browse* button in *Calibration*. If an internal standard correction is required, activate **internal standard correction** and specify the detector for the correction. For a negative signal the option **Negative Peak** has to be activated. Set the baseline limits and the integration limits and close the dialog.

Note: Please make sure that no "system peaks" or solvent peaks are within the integration limits. Otherwise the molar mass assignment might be wrong.

4. Activate automatic processing with [Definition] [Automation activated] and start the measurement.

Hints:

The default value of the allowed deviation for recalibration is set to 5%. If the deviation is larger this sample will not be recalibrated; the original peak positions will be kept. You can change this value in the dialog *WINGPC Automation Settings* (Max. Deviation %).

Note: The maximum deviation is also valid for the internal standard correction.

When you are working with "Recalibration average" you can check the number of values used for the average value in the calibration window. A right mouse click on a column header in the calibration table opens a menu. Select *Columns* and activate **Average** in the dialog.