

**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**

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

**POLYMER  
STANDARDS  
SERVICE GmbH**

**PSS**

**WINGPC  
Newsletter  
WINGPC6.2**

Application # 15

applicable for: PSS WINGPC Version 6

## Automatic Calibration with PSS WINGPC Documentation in the log-book

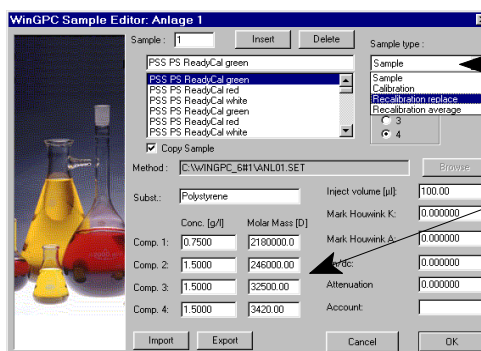
1. Create a new calibration curve in the *calibration window* with [File] [New] or using the icon in the button bar.

*Note:* Select the fit function you want to use for this column set and enter the parameters ([Calibration] [Parameters]). If you want to work with an internal standard correction, enter the value for the internal standard here.

Save the calibration curve and close it using [File] [close] or the corresponding icon in the button bar.

2. Enter the sample names and molecular weights in the sample editor (*Raw data window: [Editor] [Samples]*).

*Note:* 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)

For all calibration samples select sample type "Calibration".

3. Open the dialog *WINGPC Automation Settings* in the *method window* with [Definition] [Automation properties].

Load the new calibration curve using the *Browse* button in **Calibration**. If you want an internal standard correction, activate **internal standard correction** and specify the detector. The option **Negative Peak** has to be activated for a negative signal. Set the baseline limits and the integration limits and close the dialog.

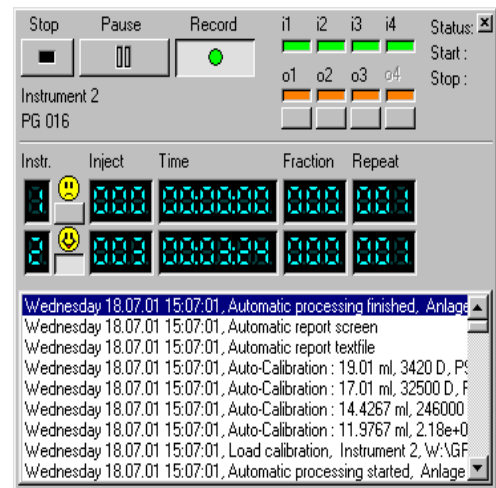
*Note:* Please make sure that no "system peaks" or solvent peaks are within the integration limits. Otherwise the molecular weight assignment might be wrong.

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

### *Documentation*

All steps of the automatic processing are documented in a log-book. The log-book is located at the bottom of the instrument information box where you start and stop measurements.

Errors, like internal standard not found, are also logged here. Entries in the log-book can not be modified by the user.



Documentation of an automatic calibration  
with 4 components in one sample