

**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**



**WINGPC**  
Newsletter  
02/01

for: PSS WINGPC Version 6

Application # 14

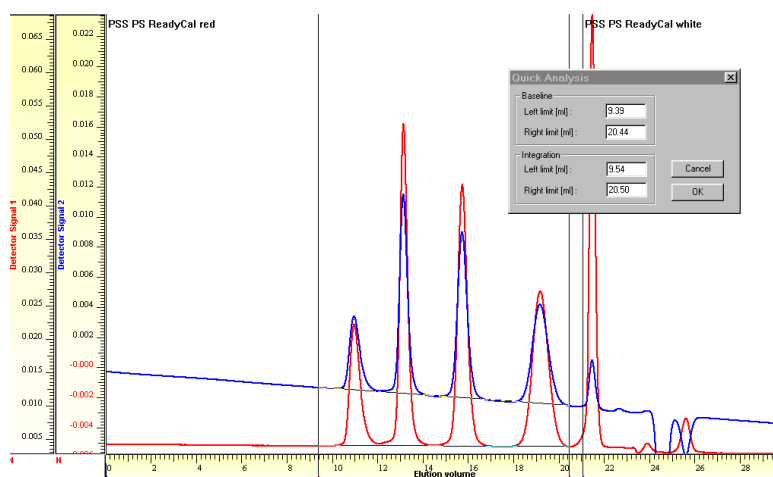
## Automatic Data Evaluation with PSS WINGPC

The five major tasks for any GPC evaluation are the internal standard correction, the assigning of the calibration curve, the setting of the baseline and integration limits and the printout of the results. These tasks are done in the raw data, the elugram and the MWD window. In WINGPC Versions < 6 only the baselines and the integration limits could be set automatically for all samples. The internal standard correction or the printout had to be done one by one. The menu item (Quick analysis) is still available in WINGPC Version 6, but the automatic data evaluation has been improved with the new automation feature.

The following section describes the differences between the quick analysis and the fully automated analysis.

### *Quick analysis*

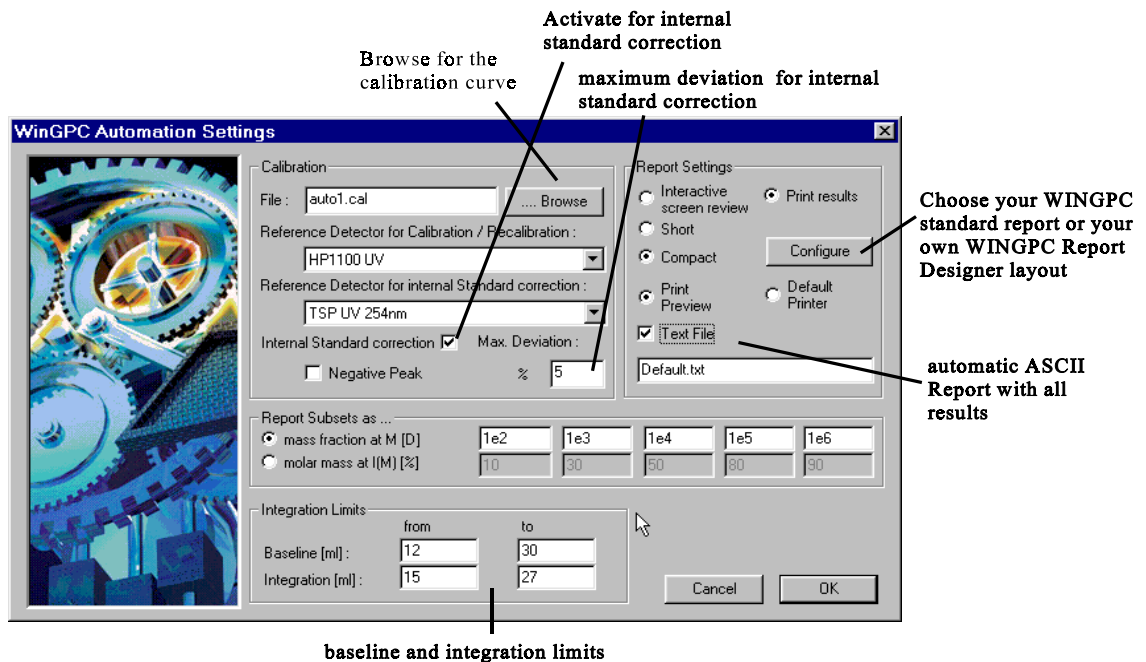
This menu item allows the setting of the baseline and the integration limits for all samples in the whole sample sequence beginning at the currently selected sample. Activate the Raw Data Window; [Options] [Analysis] [Quick analysis] opens a window where the limits for the baseline and the integration limits can be set. If the actual sample is already evaluated you will find the limits of this sample as default values! Leaving the window with OK sets these limits for all following samples. An interactive correction of baselines and/or integration limits is still possible.



### *WINGPC Automation Setup*

For a fully automated data evaluation you can use the new WINGPC Automation Setup while the data acquisition is running. All major tasks for data evaluation

are carried out by the automation feature. The automation settings are defined in your method window. The dialog WINGPC Automation Settings can be found under [Definition] [Automation properties].



Besides loading the calibration curve and setting the baseline and the integration limits you can choose if an **internal standard correction** should be carried out. You can also define the maximum deviation (here: 5%) for which the correction is allowed and which detector has to be used (here: UV detector). A correction with a negative internal standard is also possible if the option "Negative Peak" is activated.

In **Report Settings** you can decide if you want a printout or just an interactive screen review. **Configure** allows to choose the print layout. Here you have two options: a printout with a individual **WINGPC Report Designer** layout or a standard printout where you define what is on **top**, in the **middle** or on the **bottom** of the report (Graphic Elugram, Graphic MWD, Method or Report). Activation of the box **Text file** enables the automatic generation of an ASCII text file with all results and sample information (identical with the ASCII MWD Report). You can enter a descriptive file name; WINGPC adds a sample identifier to that file name automatically. After completing all entries activate the automation using [Definition] [Automation activated], enter the sample names in the sample editor of the raw data window and start data acquisition.