

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
04/01

Application # 17

for: PSS WINGPC Version 6.2 and higher

1 for All - Multi-area Data Processing saves Effort

This new WINGPC feature allows to subdivide a chromatogram into several ranges, which will be processed simultaneously. This is extremely useful, if a sample consists of several peaks or peaks with shoulders, which shall be evaluated individually. Up to now, this had to be done interactively by setting the integration limits for each peak. Now, this can be done by setting just one WINGPC option.

Performing a multi-area analysis:

1) Definition of peak range limits and selection of results

Up to 8 chromatogram ranges can be defined for which GPC results will be calculated simultaneously; all result parameters are available. The definition of peak range limits and results for the report is done best with a representative sample which has already been run. A dialog box allows to set all parameters in the raw data window menu [Options] [Analysis...] [Multi area settings] (Cf. Fig. 1).

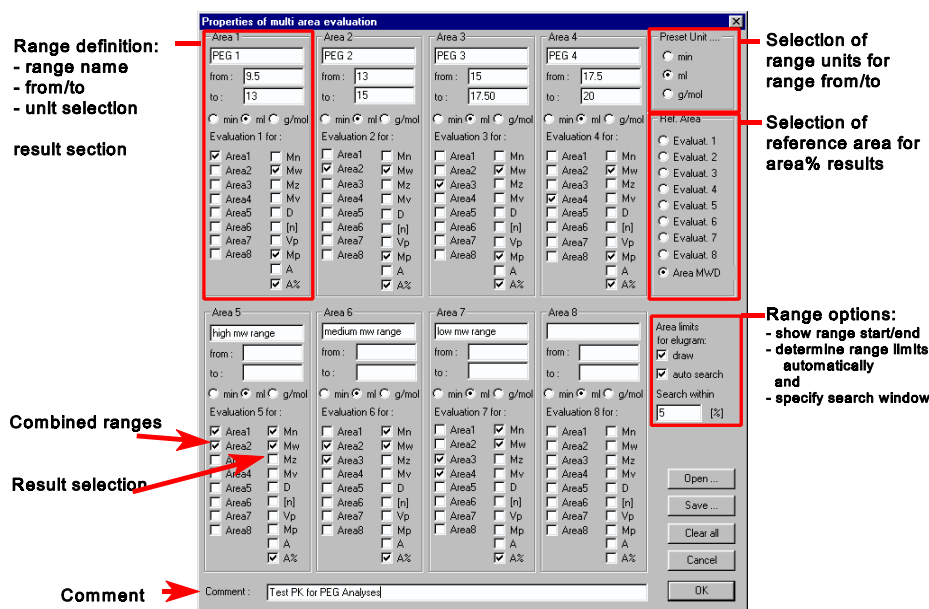


Fig. 1: Definition of multi-area ranges and result parameters

Individual peak ranges can be combined and result parameters for reporting can be selected to further increase flexibility.

