


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.



Working More Efficiently with the PSS WINGPC Button Bar

PSS WINGPC Application Note #13:

applicable: WINGPC versions 6

WINGPC 6 features a new button bar, which shall improve the accessibility of WINGPC functions and commands by minimizing the number of mouse clicks. The icons in the button bar follow the Windows programming rules and are available in all WINGPC sub-windows. In contrast to the WINGPC menus, which change with the context (which sub-window is active), the button bar is static and always identical. This allows direct point-and-shoot access to the most often used functions. The use of an icon can dramatically reduce the number of mouse clicks (e.g. opening a project file using the "file open" icon allows immediate file selection, whereas the menu commands require the selection of [Raw Data] [Load...] [File] [Load...]). Additionally all sub-windows can be toggled between foreground and background just with a mouse click.



<-----Function Range-----> <-----Window Range----->

How is the button bar used?

The button bar is divided in two segments. The segment on the left organizes often used WINGPC commands (e.g. file, print, overlay), the right range summarizes sub-window access.

A mouse click on an icon in the function range forces the software to issue the function without any delay or user selection (e.g. print report from active sub-window, save file). The icons in the window range toggle the sub-windows between foreground and background. However, a window can be hidden by other overlapping sub-windows. In such cases the icon might not know about it and a second mouse click on the button is necessary to bring the sub-window back to the foreground (the first click would bring it to the background, while the second one will bring it back to the foreground).

The overlay icon also works as an independent toggle for chromatogram overlays and calibration overlays. The toggle-state "overlay on/off" is shown as a pressed/released

Polymer Standards Service GmbH

POB 3368 D-55023 Mainz





Phone: +49-6131-9 62 39-0 Fax: +49-6131-9 62 39 11

email: info@polymer.de <http://www.polymer.de>



icon. This means that the chromatogram overlay can be active ("on") if a chromatogram window is active, while the overlay icon is "off" if the calibration window is active and no calibration curves are overlaid (and vice versa).

The following icons in the function range work context sensitive:

-  opens a project or calibration file (if calibration window is active)
-  closes the active raw data or calibration file (if calibration window is active)
-  prints report of currently active window (e.g. elugram, raw data, calibration)
-  creates page preview and clipboard copy of currently active window

