

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
04/02**

Application # 21

applicable for: PSS WinGPC Version 6 and higher

More flexibility with the WinGPC Relay control

The PSS WINCHROM and LAN-Interfaces are equipped with three output relays for controlling and saveguiding instruments, e.g. you can:

- turn off the pump after the analysis
- switch a valve after the analysis
- switch transfer valves (for 2D analysis)
- control a fraction collector
- protect the pressure transducers of your viscometer and more.

The relays can be programmed in WinGPC. The user needs to activate the relays, deactivation can be done manually by the user or automatically at the end of the measurement. Use [Definition][timed events] [acquisition stop] in the methods window to define what should be deactivated automatically after measurement stop.



Flexible Control - for every challenge the right solution

The relays control during the measurement can be set for each relay separately and depending on the functionality the relays have to do. The settings are defined in the method window using [Definition] [timed events] [Relay No.].

The selectable parameters in "signal type" depend on the chosen option in "dependent from.."

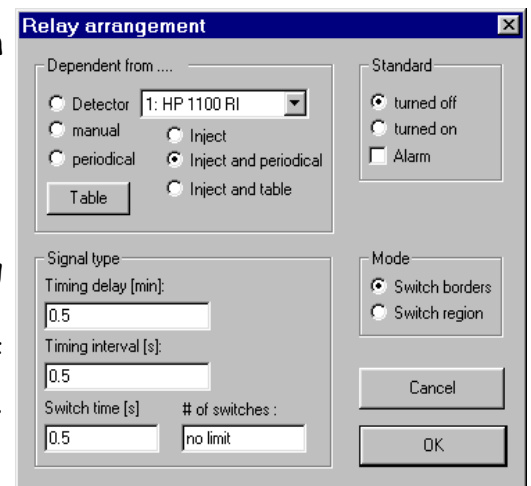
The following selections are available:

1) Detector dependent relay control

e.g. for fraction collectors or for protecting viscometers

Here you can define if within switch borders or if within a switching region the relay state is inverted.

2) Manual relay control



e.g. to turn the laser of a light scattering detector on or off
The relays are activated/deactivated on a mouse click on the relay symbols (O1,O2,O3) in the instrument information box.

3) Periodic relay control

The output relays state is inverted dependent on a given timing interval for a given switching time. The standard state can be selected.

4) Inject dependent and periodic relay control

e.g. for fraction collectors or 2D-transfer valves
The timing interval given in "signal type" is used to invert the relay state.

5) Inject dependent and relay table control

e.g. for fraction collectors or 2D-transfer valves
The WinGPC data editor can be used to edit and save tables for the relay control. Use [File] [Relay table export]) to save the table and the button *Table* in the section "dependent from.." in the window "relay arrangement".

Example relay table:

the first column contains the time in minutes after which the state of the relay is inverted, the second one the switch time in seconds (default value: 1 sec).

Row	Time [min]	Switch Time [sec]
1	4.90000E-1	1.00000E+0
2	7.00000E-1	1.00000E+0
3	1.00000E+0	1.00000E+0
4	1.50000E+0	1.00000E+0
5	2.50000E+0	1.00000E+0
6	3.00000E+0	1.00000E+0
7	3.25000E+0	1.00000E+0