



PSS WINGPC Application Note #2:

Safeguarding Viscometric Detectors Using PSS WINGPC

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.



Pressure transducers used in viscometric GPC detectors are sensitive and expensive. Exceeding the pressure limits for the transducers may result in non-reversible deformations of the membrane. Exchanging the pressure transducer is the consequence. Using PSS WINGPC's output relay it is possible to prevent the transducers to exceed their pressure limits and time and money can be saved without any further investment. PSS WINGPC software allows for signal dependant relay closure. Thus simply use WINGPC's feature to shut off the pump, if the pressure value of your transducer increases above a save value. It is even possible to have an alarm issued in case of an unattended over-pressure condition.

The relay control windows are accessible from the *method window* selecting [Definition][Timed events][Relais No.].

The example shown below closes the output relays to shut of the pump, if the pressure signal exceeds 1000 Pa.

