

VISCOTEK GPC RI DETECTOR (GPC2502)



MOLECULAR WEIGHT



MOLECULAR SIZE



MOLECULAR STRUCTURE

The Viscotek GPC RI detector is a modular differential refractive index detector for the measurement of sample concentration during GPC/SEC experiments. The detector is heated from ambient temperature to 55°C for improved baseline stability.

The reference cell can be purged manually on the instrument or automatically via software control and has an integrator output for connection to other hardware.

The system includes analogue connections for the measured signal and to initiate data collection.



Viscotek GPC RI Detector (GPC2502)

Parameter	Specification
Light source wavelength	Tungsten lamp
Refractive index range (RIU)	1 to 1.75
Flow rate range (ml/min)	0 to 3
Max pressure (MPa)	0.6
Cell volume (μ l)	9
Linearity ($\Delta\mu$ RIU)	0-1000
Noise (nRIU)	5
Drift (μ RIU/hr)	<1
Integrator out	± 1 V
PC Connection	RS232
Data rate (Hz)	5
Temperature control	Ambient, 35 - 55°C in 1°C steps (normal operation), 10 - 30°C in 1°C steps (cold room operation)
Voltage (V)	100-120/220-240, 50/60 Hz
Power consumption (W)	max 50
Size (w, d, h) (mm)	220 x 350 x 160
Weight (kg)	12

Malvern Instruments Ltd

Enigma Business Park • Grovewood Road
Malvern • Worcestershire • UK • WR14 1XZ

Tel: +44 (0)1684 892456

Fax: +44 (0)1684 892789

Malvern Instruments Worldwide

Sales and service centers in over 50 countries
for details visit www.malvern.com/contact

© Malvern Instruments Ltd 2013

more information at www.malvern.com