

VISCOTEK OmniSEC 5.1 Software



MOLECULAR WEIGHT



MOLECULAR SIZE



MOLECULAR STRUCTURE

OmniSEC software is a complete, integrated solution for the advanced analysis of proteins and polymers by GPC/SEC. OmniSEC is used to control instruments and acquire data from all products in the Viscotek range including the TDA, SEC-MALS 20, 270 and GPCmax, as well as the Zetasizer μ V and others.



Key features include:

- Modern look and feel, with optimized protein or polymer user interfaces,
- Calculation of molecular weight and molecular weight distribution from light scattering, or column calibration data.
- Calculation of intrinsic viscosity and the derivation of Mark-Houwink parameters,
- Analysis of multiple peaks,
- Copolymer / conjugate analysis by combining UV and RI detectors,
- Advanced protein analysis using OmniSEC BIO option,
- Overlay multiple chromatograms and distributions,
- Data export,
- Automatic determination of dn/dc and dA/dc .
- Automatically set baselines and limits to go from data to results in just two clicks,
- ...and many more.

OmniSEC can perform multiple types of calculations from conventional calibration, universal calibration and light scattering and can calculate all parameters related to GPC/SEC results.

	Specification
System control	up to 4 systems
Run length	no limit
Data collection rate	5 Hz
Max data channels	6
Peak detection	Manual or automatic
Baseline detection	Manual or automatic
Baseline fitting	linear up to 5th order curve
Primary calculated results	Molecular weight and distribution
	Polydispersity
	Intrinsic viscosity (η)
	Hydrodynamic radius (R_H)
	Radius of gyration (R_G)
	Peak weight fraction
	Composition
	Mark-Houwink a & $\log K$
	Branching number and frequency
	dn/dc and dA/dc
Other features	Overlay view
	Trend view
	Second virial coefficient calculator
	3D UV PDA data viewer
	DLS correlation function view
	Report designer
	OmniSEC BIO for advanced protein analysis
Integration with Waters Empower®	
Minimum System requirements	Windows XP Pro and Windows 7
	Intel dual core or hyper threaded processor
	2 Gb RAM
	150 Mb HDD space
	1024 x 768, 16 bit colour screen
	CD-ROM drive
1 free USB	



Malvern Instruments Limited
Groewood Road, Malvern,
Worcestershire, UK, WR14 1XZ

Tel +44 1684 892456
Fax +44 1684 892789

www.malvern.com

Malvern Instruments Worldwide
Sales and service centers in over
65 countries; for details visit:
www.malvern.com/contact

© Malvern Instruments Ltd 2014

Malvern Instruments is part of Spectris plc, the Precision
Instrumentation and Controls Company.

Spectris and the Spectris logo are Trade Marks of Spectris plc.



All information supplied within is correct at time of publication.

Malvern Instruments pursues a policy of continual improvement due
to technical development. We therefore reserve the right to deviate
from information, descriptions, and specifications in this publication
without notice. Malvern Instruments shall not be liable for errors
contained herein or for incidental or consequential damages in
connection with the furnishing, performance or use of this material.

Malvern and the 'hills' logo are International Trade Marks owned by
Malvern Instruments Ltd.