

THE NEW DELSAMAX™ SERIES

UNPRECEDENTED SIZING AND ZETA POTENTIAL ANALYSIS.



Characterized
by ingenuity



A POWERFUL STEP FORWARD IN YOUR PROCESS.

As an original pioneer in particle characterization, Beckman Coulter is proud to announce the innovative new DelsaMax Series, an unprecedented advancement in the speed, accuracy and outcomes of nanoparticle research. From the parallel technology of the PRO model to the sample handling capabilities of the ASSIST, DelsaMax instruments allow you to delve deeper into discovery than ever before.

DelsaMax PRO light scattering analyzer

Allows simultaneous analysis of both particle size and zeta potential of samples as small as 45 μL in as fast as one second.

DelsaMax CORE light scattering analyzer

Provides dynamic light scattering measurement from 0.4 nm to 10,000 nm, at sample volumes as low as 1 μL —the lowest sample volume available today.

DelsaMax ASSIST cell pressurization system

Optimizes particle preparation by pressurizing the sample cell to reduce bubbling, which can interfere with measurements.

DelsaMax Analysis software

Features an intuitive interface with fully customizable autocorrection analysis and the ability to create overlays, custom user variables, reports and more.

KEY INNOVATIONS

Incredible Speed

Two simultaneous detection systems allow rapid measurements (both size and zeta potential in as little as one second).

Provides 32 simultaneous measurements reducing run time and increasing accuracy.

Precision & Accuracy

Proprietary “normalization” algorithm provides robust distributions.

Edit data points or “flyers” without re-running samples.

Optimized autocorrelation function selects the particles of interest.

Unrivaled Reliability

Get up to 10 protein measurements per sample without sample degradation with both dynamic and “real” static light scattering.

Parallel, independent measuring systems in both systems offer instant cross-checking of results.

Easily overlay results to check for consistency and accuracy.

DELSAMAX PRO

Size Range	0.4 to 10,000 nm, hydrodynamic diameter (limited by particle sedimentation)
Molar Mass Range	5×10^7 g/mol (Da) (dependent on molecular shape model)
Temperature Range	4° C to 70° C
Minimum Sample Volume	45 μL
Minimum Measurement Time	1 second
Zeta Potential Measurement	
Minimum Sample Volume	170 μL , excluding tubing
Ionic Strength Range	0 to 50 mS/cm (4 times the conductivity of physiological saline)
Mobility Range	No practical limit
Mobility Size Range	2 nm to 15 μm diameter
Mobility Sensitivity	1 mg/ μL Lysozyme
Minimum Measurement Time	1 second

DELSAMAX CORE

Size Measurement	
Dynamic Light Scattering Range (Diameter-nm)	0.4 to 10,000 nm
Static Scattering Molecular Weight Range	300 to 10^6 Da (concentration dependent)
Minimum Sensitivity	0.1 mg/ μL Lysozyme
Scattering Angle	90°
Minimum Sample Volume	125 μL standard cuvette, 4 μL disposable cuvette
Correlator	512 channels (100 nsec sampling time in a multi-tau layout)
Data Acquisition Time	1 to 3,600 seconds
Minimum Measurement Time	1 second
Temperature Range	-15° C to 150° C (Quartz Cuvette) -15° C to 80° C (Disposable)

Learn more about the DelsaMax Series, visit www.delsamax.com



DelsaMax, DelsaMax PRO, DelsaMax CORE, DelsaMax ASSIST, and DelsaMax Analyzer are pending trademarks. Beckman Coulter and the stylized logo are trademarks of Beckman Coulter, Inc. and are registered with the USPTO.

For Beckman Coulter's worldwide office locations and phone numbers, please visit "Contact Us" at www.beckmancoulter.com
FL-17889A © 2013 Beckman Coulter, Inc. PRINTED IN U.S.A.