



Choosing the Right Instrument

for Cell Counting and Particle Characterization

Applications Across Sizes

Dynamic & Static Light Scattering

Coulter Principle (Electrical Impedance)

	DelsaMax Core DelsaMax Pro					Multisizer 4e													
						Multisizer 3													
Size (nM)	10	50	100	500	1000	Aperture Sizes Available (µM)													
Size (µM)						10	20	30	50	70	100	140	200	280	400	560	1000	2000	
						Effective Aperture Upper and Lower Range (µM)													
						0.2	0.4	0.6	1.0	1.4	2.0	2.8	4.0	5.6	8.0	11.2	20	40	
						8	16	24	40	56	80	112	160	224	320	448	800	1600	
Applications	Nanoparticles																		
	Proteins/Macromolecules																		
	Protein Aggregates																		
	Virus Particles																		
	Liposomes					Liposomes													
	Bacterial Studies																		
	Lg. Protein Aggregates																		
	CMP Slurries																		
	Yeast																		
	Diagnostic Beads																		
	Ink Toners																		
	Mammalian Cells																		
	Fermentation/Bioreactors																		
	Plankton/Cyanobacteria-Biofuels																		
	Abrasives - ASTM Methods																		
Plant and Plankton Cells																			
Blood Clots																			
Stem Cell/Tissue Aggregates																			



CHARACTERIZED
by ingenuity.



Cell Counting and Particle Characterization Decision Tree — Research

