



Choosing the Right Instrument

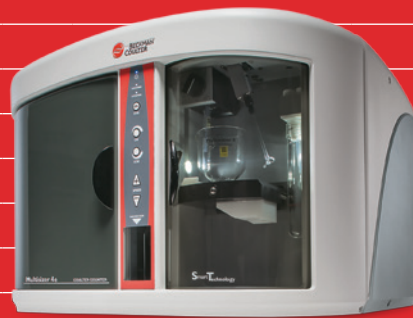
for Cell Counting and Particle Characterization

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

