

APPLICATIONS ENABLED BY MULTI-WAVELENGTH ANALYSIS WITH THE OPTIMA AUC



Nanoparticles

- Core size determination
- Drug payload

Quantum Dots

RNA Aptamers

Polysaccharides

Viral Vectors for Gene Therapy

- Empty versus full capsids

Membrane Protein-Lipid Interactions

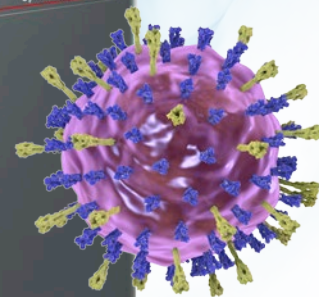
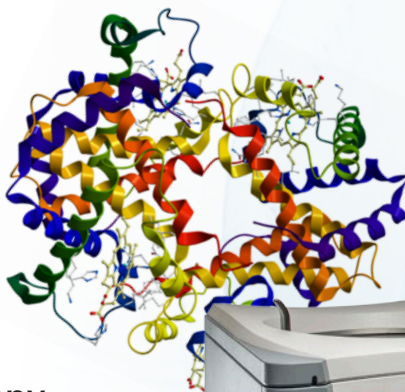
Protein-Protein Interactions

- Fusion proteins (YFP, GFP, SNAP)
- Heme Proteins
- Flavins
- Metalloproteins

Protein-Nucleic Acid Interactions

- Recombinases
- Terminases
- ATPases

...and Many More!



INSTRUMENTS & COMPONENTS

Ordering Information

Component	Part No.	Description
Optima AUC - Absorbance Only	B86438	Hardware only with Absorbance optics
Optima AUC - ABS/INT	B86437	Hardware only with Absorbance + Interference optics
Optima AUC - ABS Bundle + An-60	C00707	Optima AUC (A) + Starter Kit (1 CB, 1 torque stand, 1 Quartz Cell + 1 accessory kit) + An-60
Optima AUC - ABS/INT Bundle + An-60	C00708	Optima AUC (A/I) + Starter Kit (1 CB, 1 torque stand, 1 Quartz Cell + 1 accessory kit) + An-60
Optima AUC - ABS Bundle + An-50	C00709	Optima AUC (A) + Starter Kit (1 CB, 1 torque stand, 1 Quartz Cell + 1 accessory kit) + An-50
Optima AUC - ABS/INT Bundle + An-50	C00710	Optima AUC (A/I) + Starter Kit (1 CB, 1 torque stand, 1 Quartz Cell + 1 accessory kit) + An-50
An-60 rotor	361964	4-hole rotor rated to 60,000 rpm
An-50 rotor	363782	8 hole rotor rated to 50,000 rpm
2-sector charcoal Quartz	392772	SV cell with charcoal flow-through centerpiece and quartz windows
Torque stand Assembly	361318	Required to torque cells to prevent sample leaking
Counterbalance Kit	360219	1 counterbalance with mask + weights
Accessory Kit	392777	6 brass housing plugs
Cell housing kit	334606	2 window assemblies + cell housing



Learn more at info.beckmancoulter.com/OptimaAUC



© 2016 Beckman Coulter, Inc. All rights reserved.
Beckman Coulter, the stylized logo and the Beckman Coulter product and service names mentioned herein are trademarks or registered trademarks of Beckman Coulter, Inc. in the United States and other countries.

For Beckman Coulter's worldwide office locations and phone numbers, please visit "Contact Us" at beckman.com

CENT-2214FLY12.16