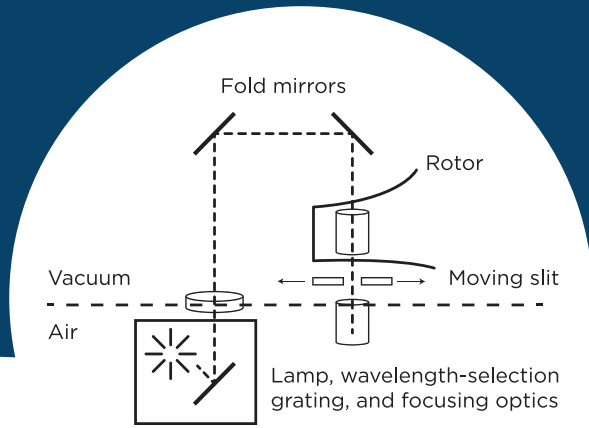




# ANALYTICAL ULTRACENTRIFUGE PRODUCT SPEC COMPARISON



## PROTEOMELAB XL-A / XL-I

## OPTICAL SYSTEMS

## OPTIMA AUC\*\*

ABS: 90sec/cell INT: 5 sec/scan	FASTEST DATA ACQUISITION RATE	ABS: <7 sec/sector* INT: <5 sec/scan
3	MAX # OF WAVELENGTHS	20
+/- 3 nm	WAVELENGTH PRECISION	+/- 0.5 nm
30 µm	LOWEST RADIAL RESOLUTION	10 µm
50 Hz	ABSORBANCE FLASH LAMP FREQUENCY	300 Hz
2048 x 96 pixels	CCD CAMERA SPECIFICATIONS	2048 x 1088 pixels
≥ 4 fringes/cell	INTERFERENCE FRINGES	≥ 10 fringes/cell
ABS: .005 - 1.5 mg/mL INT: .025 - 3-4 mg/ml	USABLE CONCENTRATION RANGES	ABS: .005 - ~1-2 mg/mL Luteinizing Hormone INT: .025 - ~4-5 mg/mL BSA**

\* At Optimal rotor speeds, at 10 µm radial resolution

\*\*Product under development - specifications pending system verification



© 2016 Beckman Coulter, Inc. All rights reserved. Beckman Coulter, the stylized logo, and the Beckman Coulter product and service marks 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](http://beckman.com)

CENT-2023IFG10.16