

Optima XE

The essential ultracentrifuge.



Brilliance
at every turn.



Optima XE

It's Beckman built.

If you're looking for practicality, but desire the functionality and performance of an ultracentrifuge, you've found the perfect instrument: the Optima XE. This ultracentrifuge provides all of the features you'd expect, along with some very beneficial enhancements you won't find anywhere else.

Large touch-screen display

Displays more data. Easy to read and navigate. Viewable from across the lab.

Multilingual software

Includes nine languages for error-free operation.

Zonal and continuous flow operation screens

Intuitive, graphical step-by-step interface simplifies workflow.

Context-sensitive on-screen help

Find answers easily. Reduce training time and user error.



With total-system design in mind from the start, Beckman Coulter created the Optima XE to simplify use, increase productivity, and reduce error. You'll applaud its intelligent user interface, multi-layered BioSafety features, and energy efficiency. And, of course, our promise of total package satisfaction

is backed by our exclusive rotors, labware, and unparalleled service.

By challenging convention and raising the bar, Beckman Coulter is setting an all-new standard. Again. With advanced intelligence and intuitive design, the Optima XE will revolutionize your world.



Energy efficient

Regenerative braking returns energy to local circuit reducing utility costs. Highly efficient thermoelectric cooling results in lower power consumption. Uses the same amount of power as a 60-watt light bulb while idling.

Improved power tolerance

Maintains runs during power fluctuations. Automatically switches transformer taps to handle a wider tolerated range of 180 VAC to 264 VAC. Tracks power interruption data to convey if parameters remained within an acceptable range.

Quiet-drive technology

Designed for lowest drive noise—this is the quietest floor model on the market. Operates at less than 51 dBA.

BioSafe and BioSafety are terms intended to describe the enhanced biocontainment features of our products.



Optima XE

The Beckman of centrifuges.

Beckman Coulter is the global leader in ultracentrifuges. Our quality stands the test of time—every piece of our equipment is engineered with planned durability, so you can trust in reliable performance year after year, after year. And when you invest in Beckman Coulter instruments, you're backed by an incomparable worldwide support organization. All to help you maintain your investment and ensure top performance from your lab.

FEATURES

	XE	XPN
Giant LCD touch-screen display	■	■
Intuitive, multilingual interface	■	■
Improved power tolerance	■	■
Compatible with existing Beckman Coulter rotors	■	■
Step-by-step zonal/CF operation screens	■	■
Rotor tracking by serial number		■
Remote monitoring and control		■
Password protection and three user access levels		■
Exportable run logs and electronic signatures		■
eXpert simulation software		■
Speed/temperature vs. time plot		■
User-defined programs		■

SPECIFICATIONS

	XE-100	XE-90
Maximum Speed (rpm)	100,000	90,000
Maximum RCF (x g)	802,000	694,000
Speed Control	±2 rpm of set speed (above 1,000 rpm)	
Set Temperature	0 to 40° C in 1° C increments	
Temperature Control	±0.5° C of set temperature*	
Ambient Operating Range	10° to 35° C	
Accel/Decel Profiles	10/11	
Braking	Regenerative, power reclamation	
Operating Languages	English, French, Italian, German, Spanish, Japanese, Russian, Korean, and Chinese (Simplified)	
Heat Output	3400 BTU/hr (1.0 kW)	
Drive Cooling	Air-cooled	
Refrigeration System	Thermoelectric—no CFCs, ODCs	
Sound Level	<51 dBa	
HEPA Filter	Available	
Power Requirement	200-240V, 30 A, 50/60 Hz	
Dimensions	94W x 68.1D x 125.7H cm (37W x 26.8D x 49.5H in)	
Weight	485 kg (1068 lb)	
Order Number	A94516	A94471

*±0.8° C when using SW 32 swinging bucket rotor.



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

B2011-12560-2.5K-LC

© 2011 Beckman Coulter, Inc.

PRINTED IN U.S.A.