

Optima MAX-XP

The premier tabletop ultracentrifuge.



Brilliance
at every turn.



Optima MAX-XP

Unprecedented. Uncomplicated.

How does the Optima MAX-XP compare to other tabletop ultracentrifuges? It doesn't. The MAX-XP features all the functionality and efficiency of our entry-level model, plus unprecedented speed and technological advancements you simply will not find elsewhere. The premier model in the

Remote capabilities

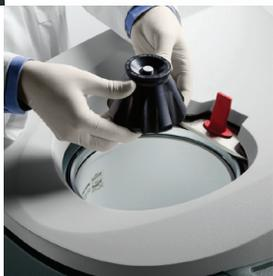
Advanced networking capabilities allow you to operate and monitor instruments from outside the lab.

Security/tracking features

Password-controlled access, user and run data tracking, audit trail documentation, rotor logs and more. A USB port simplifies export of run histories and reports.

User-friendly, multilingual interface

Full-color LCD touch-screen with nine languages—the most available in the industry—for error-free operation. Highly flexible options, enabling users to set multiple programs of up to five steps each or delay start times.



Optima MAX Series, the MAX-XP certainly lives up to its top-of-the-line status.

The MAX-XP boasts the most sophisticated software platform available in a tabletop, and yet it's stunningly simple to use and customize. Every detail is designed to protect and streamline your process—from security and tracking features to remote

capabilities that allow you to have full confidence in your laboratory's operations, even when you're away.

This robust, comprehensive platform is a fantastic addition or upgrade to your lab. And the ingenuity built into the MAX-XP speaks volumes about the total-system design that has come to define Beckman Coulter solutions.

Complete compatibility

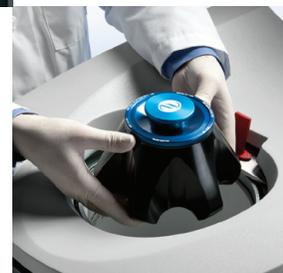
Fully compatible with our extensive existing library of tabletop rotors and labware, including fixed-angle, swinging bucket and unique near-vertical tube (NVT) rotor models.

Multi-level approach to BioSafety

HEPA filtration, BioSafe* tubes, fluid containment annulus rotors, fits in standard BioSafety* hood.

Widely trusted

Relied upon by scientists worldwide; 95% of the published protocols citing tabletop or micro-ultracentrifuges reference a Beckman Coulter instrument or rotor:**



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

**Based on 2009 Google Scholar searches.

Optima MAX-XP

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

	MAX-TL	MAX-XP
User-defined programs	Virtually Limitless	
Available rotors	9	16
Intuitive, multilingual interface	■	■
Numerical vacuum display	■	■
Context-sensitive, on-screen help	■	■
RPM or RCF user selectable		■
Password protection/User ID		■
User customizable		■
Electronic record-keeping support		■
Delay run		■
Run hold		■
Pulse function		■
Detailed run histories		■
Rotor tracking by serial number		■
Export data via USB port		■
On-screen rotor library		■
Remote monitoring & control option		■

SPECIFICATIONS

	MAX-XP
Maximum Speed (rpm)	150,000
Maximum RCF (x g)	1,019,000
Speed Control	±50 rpm of set speed
Set Temperature	0° C to 40° C in 1° C increments
Temperature Control	±2° C of set temperature
Ambient Operating Range	15° C to 35° C
Accel/Decel Profiles	10/11
Operating Languages	English, French, Italian, German, Spanish, Russian, Japanese, Portuguese, and Chinese (Simplified)
Heat Output	2400 BTU/hr (0.7 kW)
Refrigeration System	Thermoelectric—no CFCs, ODCs
Sound Level	<47 dBA
HEPA Filter	Available (Sold Separately)
Remote Control	Available (Sold Separately)
Power Requirement	100-240V, 50/60 Hz
Dimensions	73.9W x 61.7D x 34.5H cm (29.1W x 24.3D x 13.6H in)
Weight	105 kg (230 lb)
Order Number	393315



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

B2012-12945-7K-LC

© 2012 Beckman Coulter, Inc.

PRINTED IN U.S.A.