

Optima MAX-TL

The essential tabletop ultracentrifuge.



Brilliance
at every turn.



Optima MAX-TL

Simple sophistication.

The Optima MAX Series of tabletop ultracentrifuges is the absolute top of the line. And while the MAX-TL tabletop is our entry-level model, it's so much more than just another tabletop. The MAX-TL delivers optimum functionality and efficiency within a compact, quiet package. You can reach the speeds

Multilingual software

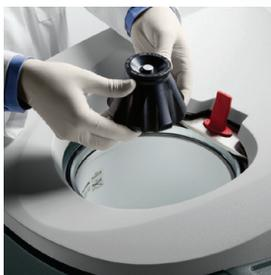
Includes nine languages—the most available in the industry—for error-free operation across the globe.

Full-color LCD touch-screen

Easy to navigate and allows you to quickly set up and complete a run.

Multi-level approach to BioSafety

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



you need to match your separation demands, while your individual experience remains uncomplicated and uncompromised.

The interface makes navigation and operation a breeze and the software includes nine languages, so labs across the globe can reduce errors. And for added

versatility, the MAX-TL is compatible with the existing Beckman Coulter TL-Series rotors and labware.

Performance. Convenience. Usability. All components of the total-system design that has come to define Beckman Coulter solutions. And all components, you'll find, are abundantly supplied in the MAX-TL.

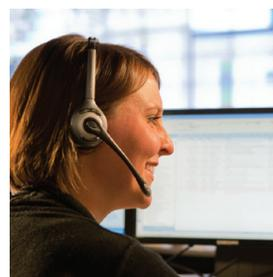


Complete compatibility

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

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-TL

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

	Virtually Limitless	
User-defined programs		
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-TL
Maximum Speed (rpm)	120,000
Maximum RCF ($\times g$)	657,000
Speed Control	± 50 rpm of set speed
Set Temperature	0° C to 40° C in 1° C increments
Temperature Control	$\pm 2^\circ$ 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)
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	A95761



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

B2012-12946-7K-LC

© 2012 Beckman Coulter, Inc.

PRINTED IN U.S.A.