

# OPTIMA MAX SERIES ULTRACENTRIFUGES

TOP-OF-THE-LINE TABLETOPS.



**Brilliance**  
*at every turn.*



ULTRA-POWERFUL INTUITIVE  
DESIGNED TO FIT ANY LAB  
GLOBAL COMPATIBILITY REMOTE  
MONITORING BIOSAFETY SO  
SOPHISTICATED YET SO SIMPLE



**1947** Model E ultracentrifuge

**1963** First titanium ultracentrifuge rotor

**1998** First centrifuge to surpass one million g-force

**2010** Centrifuge Center of Excellence established in Indianapolis, Indiana

### > The Henry Ford of life sciences.

Not long after he invented the now-famous acidimeter, founder Arnold O. Beckman, Ph.D., released the DU Spectrophotometer<sup>†</sup> in 1940. Dubbed “the Model T of life sciences,” Beckman’s spectrophotometer streamlined analytical chemistry, offering more rapid and accurate results than ever before.

More than 70 years later, the tradition of creating solutions that accelerate scientific discovery lives on at Beckman Coulter. And at the top of the line is our Optima MAX Series tabletop ultracentrifuges.

As an inspired addition to the Optima line, Optima MAX instruments are the most sophisticated, comprehensive tabletops in the industry. And like Ford’s Model T, precise engineering is packaged in a stunningly simple, intuitive platform, designed for scientists and laboratories the world over.

It’s this kind of ingenuity that has come to be synonymous with the Beckman Coulter name. And as our founder established, we continue to move the industry forward at the speed of innovation.



<sup>†</sup>Photo courtesy of National Optical Astronomy Observatory/Association of Universities for Research in Astronomy/National Science Foundation.

# OPTIMA MAX SERIES

WORTHY OF THE BECKMAN NAME.

Our history is rich with innovations that have benefited the greater good, having found thousands of applications in science, industry, and medicine. It is this heritage that proves 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.

Our equipment isn't the only thing that's exceptional. When you invest in Beckman Coulter instruments, you're backed by an incomparable worldwide

support organization. You can rely on easy Internet or phone access, field service teams, customizable service plans, original Beckman Coulter parts, and more. All to help you maintain your investment and ensure top performance from your lab.

## Optima MAX-XP and TL

Expanding tabletop possibilities.

Elevating your laboratory's potential starts with the optimal tabletop ultracentrifuge. And the Optima MAX Series is the absolute top of the line. Reaching up to 150,000 rpm and over 1,000,000 x g, Optima MAX instruments set the standard in tabletop speed.

Our entry-level model, the TL, is designed for optimum functionality and efficiency within a compact, quiet package. Our premium model, the XP, features advanced software with password-controlled security features, remote control capabilities, and a variety of tracking and archiving features.

Both models feature multi-layered BioSafety\* measures to ensure an uncompromised working environment, and both work perfectly with our existing library of rotors and labware.

The Optima MAX Series takes sophistication and total system design to a whole new level. And there has never been a tabletop series so worthy of the Beckman name.

### 1 Total-system design

Instruments, rotors and labware are designed to work together perfectly.

### 2 Remote capabilities (XP only)

Operate and monitor instruments from outside the lab.

### 3 Multi-level approach to BioSafety\*

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

### 4 Security and tracking features (XP only)

Password-controlled access, user and run data tracking, audit trail documentation, and rotor logs.

### 5 User-friendly, multilingual interface

Full-color LCD touch-screen with nine languages—the most available in the industry.

### 6 Versatility

The MAX-XP is compatible with our complete existing library of tabletop rotors and labware. The MAX-TL is compatible with our existing TL Series rotors and labware.

### 7 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 Series

Designed for any space. And any demand.

The Optima MAX Series of tabletop ultracentrifuges delivers the power and performance you need in a compact platform that fits seamlessly into any laboratory.

But Optima MAX instruments are far from the average tabletops. They bring

more sophistication than ever, without sacrificing usability. All in an effort to accelerate the rate at which you make new and important scientific discoveries, something to which we at Beckman Coulter have been whole-heartedly dedicated since our inception.

“There is no satisfactory substitute for excellence.”

*Arnold O. Beckman, PhD*





## A global presence. Focused on individuals.

Our innovations happen behind the scenes. Yet we touch people the world over by developing, manufacturing and marketing discovery systems and products that improve human health.

For over 75 years, we've been building a global reputation in hospitals, labs, and universities, where our life science research instruments are relied upon to perform vital roles day in and day out. Whether helping to solve complex biological problems, investigate the causes of disease, or find potential new cures, Beckman Coulter is focused on innovations that ultimately lead to healthier lives.

[www.beckmanbuilt.com](http://www.beckmanbuilt.com)



Beckman Coulter, the stylized logo, and DU 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](http://www.beckmancoulter.com)

B2012-12944-5K-LC

© 2012 Beckman Coulter, Inc.

PRINTED IN U.S.A.