


# OPTIMA X SERIES ULTRACENTRIFUGES

PREPARE TO BE MOVED.

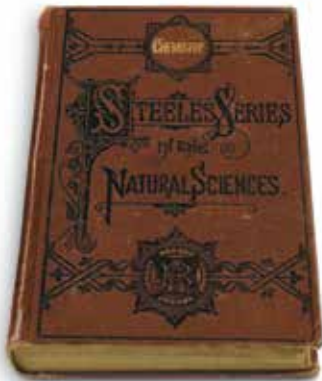


Brilliance  
*at every turn.*





ADVANCED INTELLIGENCE  
OVERSIZED TOUCH-SCREEN  
SO QUIET MULTILINGUAL  
ENERGY EFFICIENT SECURE  
CENTRIFUGATION REDEFINED  
REVOLUTIONARY



**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

## > And it all began with a dusty, dog-eared book.

Founder Arnold O. Beckman's first major discovery? It happened to be an old chemistry textbook, Steele's **Fourteen Weeks in Science**, which he found while rummaging around as a boy up in the family attic. After that, his curiosity was insatiable.

It's in that same spirit Beckman Coulter continues to transform the realm of centrifugation. How can we make it easier to use, run quieter, and offer a more intelligent interface? What is the most advanced, yet intuitive, instrument we can possibly design?

It's the Optima X Series. By challenging convention and raising the bar, we're setting an all-new standard. Again. The most sophisticated centrifuge in the industry, the Optima X Series will revolutionize your world.

At Beckman Coulter, we have a ravenous appetite for the advancement of scientific discovery. How else could fourteen weeks turn into over three-quarters of a century of innovation?



# OPTIMA X SERIES

IT'S BECKMAN BUILT.

Our history is rife 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

## Optima XE and XPN

Welcome to a whole new level.

With all of the Optima X Series' features and enhancements, you'll feel like a kid in a candy shop. Where to begin?

How about starting with the visibility and control of our intelligent user interface that makes the Optima X Series a breeze to operate. Our entry-level model, the XE, delivers all of the features required to quickly set up and complete a run. Our premium model, the XPN, boasts networking capability and a variety of customizable security and

tracking features, ensuring chain of custody and compliance. Both models feature multi-layered BioSafety\* features in order to maintain a safe working environment. And, of course, our promise of total package satisfaction is backed by our exclusive rotors, labware, and unparalleled service.

As you can see, the Optima X Series is Beckman built, with total-system design in mind from the start.



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



with planned durability, so you can trust in reliable performance year after 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.



**1 Large touch-screen display**

Easy to read and navigate. Viewable from across the lab. Intuitive graphical interface with on-screen help. Multilingual software for error-free operation.

**2 Remote monitoring and control (XPN only)**

Saves time. Start, stop, and monitor from computer or smart phone. Receive diagnostic alerts via email.

**3 Energy efficient**

Maintains runs during power fluctuations. Regenerative braking returns energy to local circuit. Highly efficient thermoelectric cooling results in lower power consumption.

**4 Power on-board simulation and calculation tools (XPN only)**

Saves time and samples. Proprietary software optimizes protocols before running to ensure maximum efficiency. Quickly perform common calculations and conversions.

**5 Security and tracking features (XPN only)**

Manage rotor life by serial number. Track detailed usage history. Password protection and electronic signatures help you maintain chain of custody and compliance.

**6 Quiet-drive technology**

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





# Optima X Series

The Beckman of centrifuges.

---

Making your job easier and more fruitful from the start—now that's an unorthodox idea, but we wouldn't have it any other way. Beckman Coulter helps create the world's most productive and efficient laboratories by providing the finest ultracentrifuges available.

The Optima X Series is no exception. It's revolutionary. It has become the all-new standard. And like all of our instruments, each one is a precision-made beauty worthy of the Beckman Coulter name.

Prepare to be moved. Experience perfection. Meet the Optima X Series and let the accolades begin.

“There is no satisfactory substitute for excellence.”

Arnold O. Beckman, PhD





## A global presence. Focused on individuals.

We may not be a household name. 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 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](http://www.beckmancoulter.com)

B2011-12558-10K-LC

© 2011 Beckman Coulter, Inc.

PRINTED IN U.S.A.