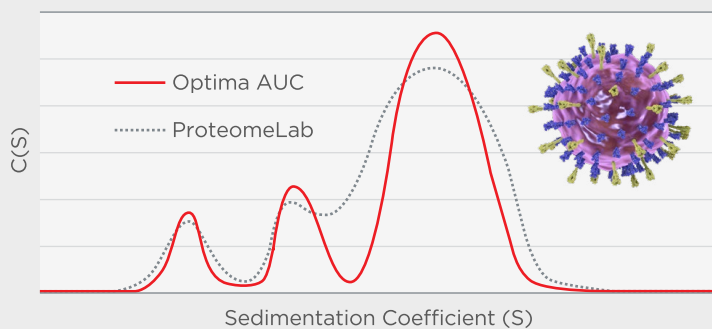


Optima AUC Advantages

Convenient Operation & Control

The Optima AUC includes new features to improve the user experience from method setup to data recovery. Quick-turn absorbance and interference optics, paired with remote instrument setup and viewing, offer streamlined method setup and live data analysis. Importantly, the Optima AUC maintains compatibility with existing Beckman AUC rotors and accessories.



Increased Data Quality & Resolution

Resolving similar components, such as viral particle capsids, can be challenging. However, the faster scan rate and improved radial resolution in the Optima AUC enable enhanced discrimination of such materials. Along with greater wavelength precision, these enhancements yield more reliable and consistent results.

Time-saving Features Improve Efficiency

Using the Optima AUC, time savings is achieved in several ways. For example, multi-wavelength capabilities enable up to 20 absorbance analyses and simultaneous interference detection in one experiment. Further, improved scan rates allow for shorter run times with greater resolution. These time-saving features of the Optima AUC aid in improving sample throughput and workflow efficiency.



Summary

The Optima AUC offers many advancements to provide convenient analytical ultracentrifugation in a modern package. Click [here](#) to learn more about the new features in the Optima AUC or to get started with AUC. Also, visit beckman.com/spINSIGHTS to explore other topics.



© 2019 Beckman Coulter, Inc. All rights reserved. Beckman Coulter, the stylized logo and the Beckman Coulter product and service names mentioned herein are trademarks or registered trademarks of Beckman Coulter, Inc. in the United States and other countries.

For Beckman Coulter's worldwide office locations and phone numbers, please visit Contact Us at beckman.com

CENT-5861FLY08.19