

Protein Purification Workflow – Centrifugation Highlights



Centrifuge to harvest cells containing protein



• 200 xg 15min

Harvest cells producing proteinfrom media

Cell lysis and treatment

Speed and Volume

Centrifuge to clear cell debris

- 3000 xg 15min
- keep supernatant

Separate insoluble (junk) from soluble protein

0.58*

Protein Isolation



Optima XPN 100

Density gradients to separate native protein, protein-ligand, oligomers or aggregates

- Time
- Speed
- Gradient



Hermatic sealing minimize air-liquid interface

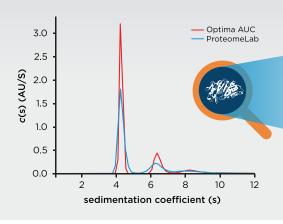
Ultracentrifugation is efficient as it can most effectively separate different protein products

Optima AUC

Protein characterization

• Biophysical Experiments with AUC

Analytical Ultracentrifugation can uniquely provide biophysical parameter and characterize protein-based products











© 2017 Beckman Coulter, Inc. All rights reserved. Beckman Coulter, the stylized logo, and the Beckman Coulter product and service marks mentioned herein are trademarks or registered trademarks of Beckman Coulter, Inc. in the United States and other countries.