

Virus Production Workflow – Centrifugation Highlights

70 YEARS OF CENTRIFUGATION

Produce ▶ Harvest ▶ Isolate ▶ Lyse ▶ Clear ▶ Purify ▶ Characterize



Preparation of production cell line



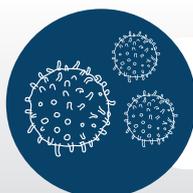
Preparation of plasmid DNA



Maintain and separate live and dead cells in a temperature controlled environment



Transformation
~ 3 days



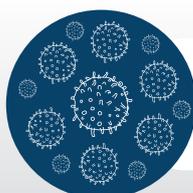
Viral Amplification
~ 2 days



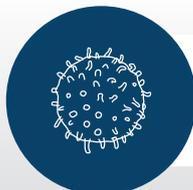
Isolate clean and high number of DNA copies



Test viral expression
~ 4 days



Large Scale Production
~ 5 days



Virus purification
~ 2 days



Avanti JXN-30



Allegra X-15R