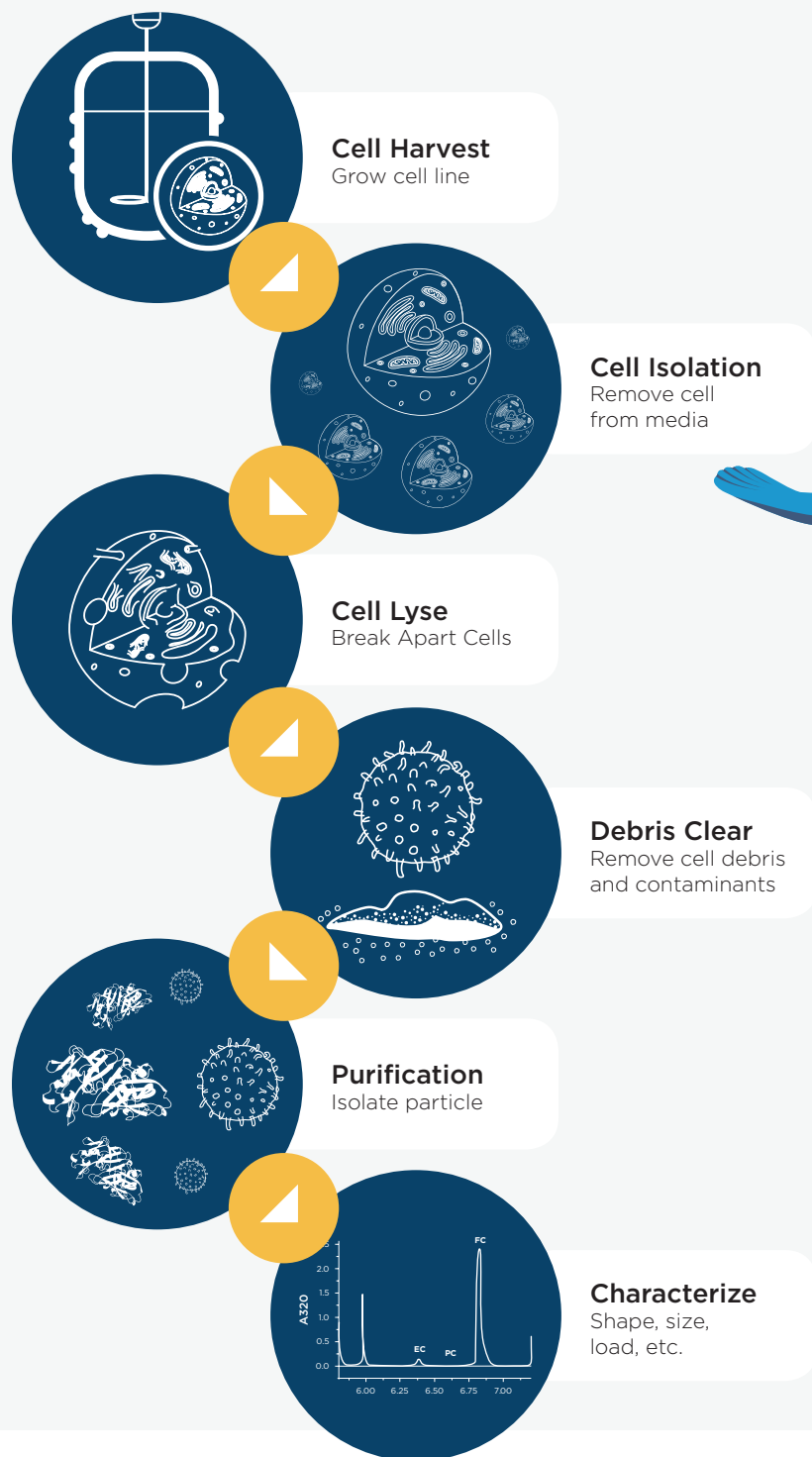


Biologics Preparation Workflow



**2 Main Applications
of Focus: Virus and
Protein Purification**