



**Vector Construction**

Equipped with a therapeutic antibody's genetic sequence, scientists insert that sequence into a molecular vehicle called a plasmid vector

**Stable Transfection and Selection**

Plasmid vectors deliver the desirable DNA construct to thousands of host cells, spurring antibody expression

*Cell Line Engineering and Selection*

**Single Cell Depositing and Monoclonality Verification**

*Cell Line Characterization*

**Expansion and Characterization**

**Scale-Up**

Rigorous checks and balances are required to ensure cell line efficiency and viability during scale up

*Process Development*

**Master Cell Banking**

Selected cell lines are stored in a master cell bank, where they're cryopreserved until future use

# Cell Line Development:

## Accelerating Answers with Automation

From generating vaccines and checkpoint inhibitors to developing revolutionary gene therapies, cell line development possesses the power to accelerate innovation in the life sciences. It's a critical cog in the biologics production wheel. Despite advances in technology, manual bottlenecks still plague this workflow. At Beckman Coulter Life Sciences, our goal is to empower and outfit your cell line development workflow with leading-edge automation technology that will inform, unify and accelerate your cell line development workflow, allowing you to produce and select efficient cell lines. But we're more than a technological solution for scientific excellence. Our expert architects in cell line development are here to answer your questions. From antibody engineering to Master Bank Selection, we're here for you at all touchpoints. Explore our array of instrumentation, software, and reagent solutions that'll transform your cell line development workflow into a nimble, efficient and results-driven enterprise.