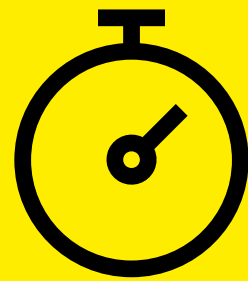


# Five Ways to Accelerate Your Cell Line Development Process

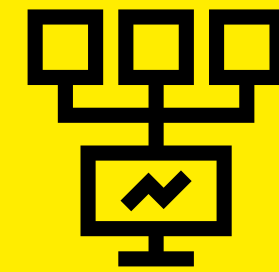
## 01. Perform Early Characterization

Implement comprehensive testing programs as soon as possible in your process development.



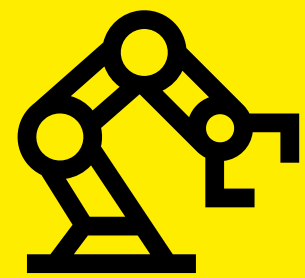
## 02. Make Use of Integrated Platforms

Consolidated systems simplify the development process.



## 03. Access External Services

Cell line development and characterization packages give you access to a suite of assays designed to simplify and fast-track your path to regulatory approval.



## 04. Adopt Automation

Multi-parallel technologies facilitate high-throughput cell culture from initial research phase to preparing for commercial production.



## 05. Take Advantage of Analytic Software

Rapid access to high-quality data is essential for measuring success, reducing risk, and predicting future outcomes.

**SARTORIUS**