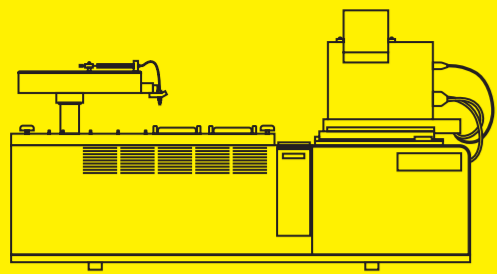


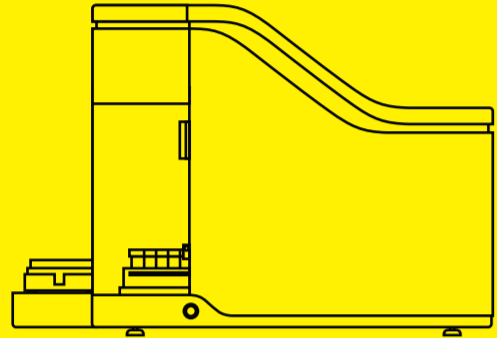
# Innovations for Stable, Scalable and Productive Cell Lines

Simplify Your Development Process from Clone Selection to Working Cell Bank



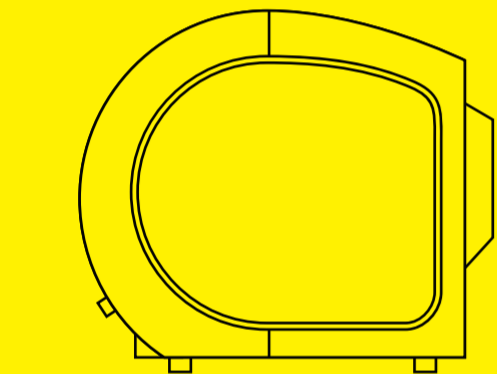
## CellCelector Automated Picking Platform

- Single cell cloning and picking
- Integrated monoclonality, viability and productivity proof
- Characterize tens of thousands of clones in parallel



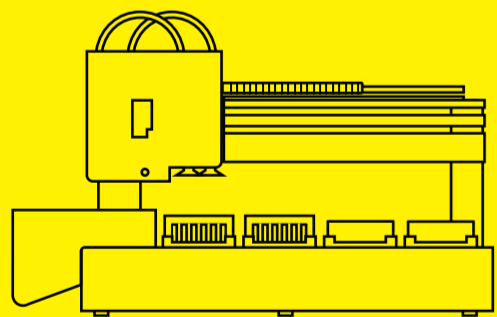
## iQue® Advanced Flow Cytometry Platform

- Viable cell concentration (VCC) determination
- Viability determination
- Productivity screening



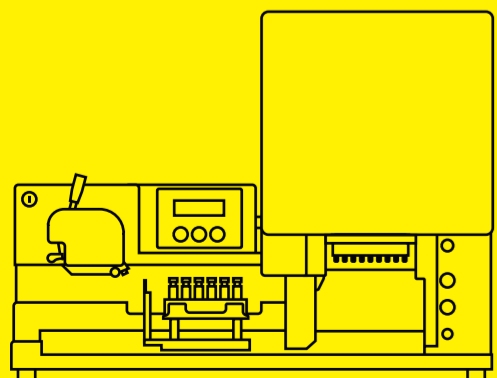
## Octet® Label-Free Biomolecular Interaction Analysis

- Titer determination
- Evaluate critical quality attributes (CQAs)
- Quality and stability monitoring



## Ambr® 15 Cell Culture

- Multiparallel cultivation and clone ranking
- Media screening and process optimization



## Fill-It Automated Cell Banking System

- Cryovial cell banking



### Gene Cloning and Initial Clone Selection

Rapid screening of minipools for product titer, cell health and growth



### Clone Selection and Confirmatory Analytics

Early identification of high productivity clones with optimal critical quality attributes and target specificities



### Cultivation and Media Optimization

Screen culture conditions and media formulations for optimal growth and protein production



### Cell Line Evaluation and Characterization

Characterize top clone for stability, productivity and quality attributes of the protein



### Cell Banking

Research Cell Bank (RCB), Master Cell Bank (MCB), and Working Cell Bank (WCB) preparation

