

# CellCelector Automated Cell -**Imaging and Picking System**

- High-throughput single plasma B cell and hybridoma clone screening for antibody discovery
- Massively parallel cocultivation of different cell types for functional studies
- Rare single cell detection and isolation

## iQue<sup>®</sup> Advanced Flow **Cytometry Platform**

- High content, multifactorial results for greater biological insight
- Unparalleled speed allows library screening in days, not months

# Incucyte<sup>®</sup> Live-Cell Analysis System

- Remote image acquisition and analysis while cells remain safely within their tissue culture environment
- More biologically and physiologically relevant insights

#### **Octet® Label-Free Biomolecular Interaction Analysis**

- Parallel processing of up to 96 samples; non-destructive and in real-time with BLI
- Quantitation or binding kinetics measured directly in unpurified samples with BLI
- High-quality hit characterization based on SPR

# Sartoclear Dynamics<sup>®</sup> Lab V Kits

Harvesting of proteins in one step

### Vivaspin<sup>®</sup> | Vivaflow<sup>®</sup> Ultrafilters and Vivapure<sup>®</sup> Membrane Adsorbers

Concentration, diafiltration and purification of antibodies and antigens

# **Biologics Molecule Development** Workflow—Sartorius Solutions for Every Step of the Process



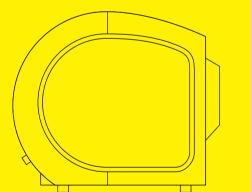


### **Target Identification**

Test target hypothesis to establish the proof of concept by modulating or silencing a protein of interest under different conditions

**Library Screening** 

library to identify positive hits



SVIPCTRV3



Simplify and Accelerate Your Development Process From Ideation to the Candidate Molecule





Screen whole-cell binding hybridoma, single B cells or phage



**Hit Confirmation** 

Confirm hits in kinetic binding assays



Lead Evaluation and Optimization Assess and optimize properties of lead candidates: titer, affinity to target, epitope binning, cross-reactivity



Candidate Characterization Characterize lead candidates for receptor binding and other functional properties

