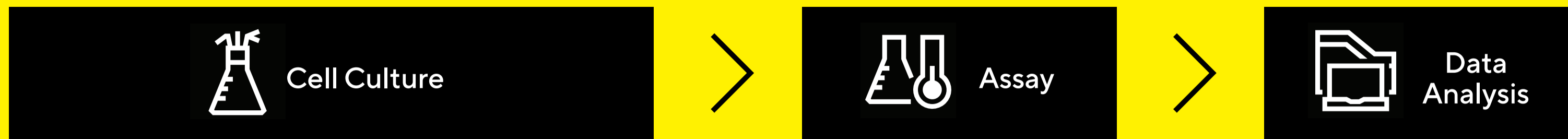
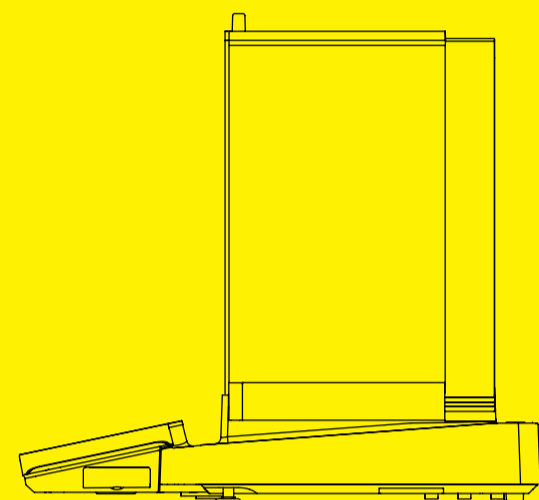


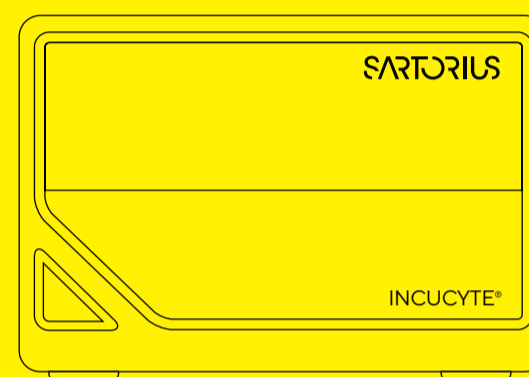
Cell Culture Workflow Solutions



RPMI 1640 Medium



Cubis® II Premium Lab Balances



Incucyte® SX5



Incucyte® Software

Growth and Maintenance (Powder or Liquid)

- RPMI 1640 Medium
- MEM-Eagle, Earle's salts
- DMEM
- F-10 Nutrient Mix
- F-12 Nutrient Mix

Supplementation

- Fetal Bovine Serum (FBS)
- Newborn Calf Serum (NBCS)

Mycoplasma Detection

- Microsart® Mycoplasma qPCR Kit

Media Preparation

- Arium® Water Purification Systems
- Cubis® II Premium Lab Balances
- Minisart® Syringe Filters
- Sartolab® RF | BT Vacuum Filters

Liquid Media Handling

- Picus® and Picus® Nxt Electronic Pipettes
- Tacta® Mechanical Pipettes
- Safetyspace® Filter Tips

Cryopreservation

- NutriFreez® D10 Cryopreservation Medium

Cell Health and Viability, Apoptosis, Cytotoxicity

- Incucyte® Live-Cell Analysis Instruments

Protein Purification and Concentration

- Sartoclear Dynamics® Lab
- Vivaspin® Turbo 15 RC
- Vivaflow Crossflow

Image Acquisition, Viewing and Analysis

- Incucyte® Base Analysis Software