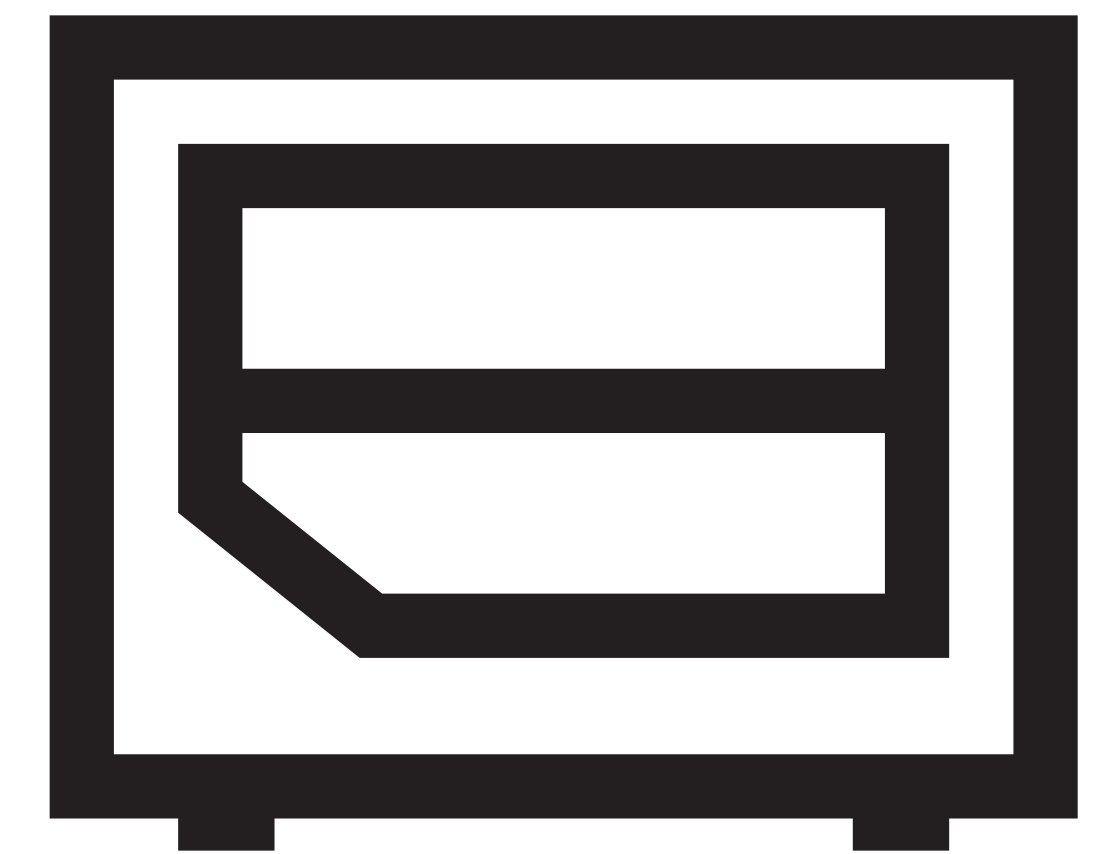
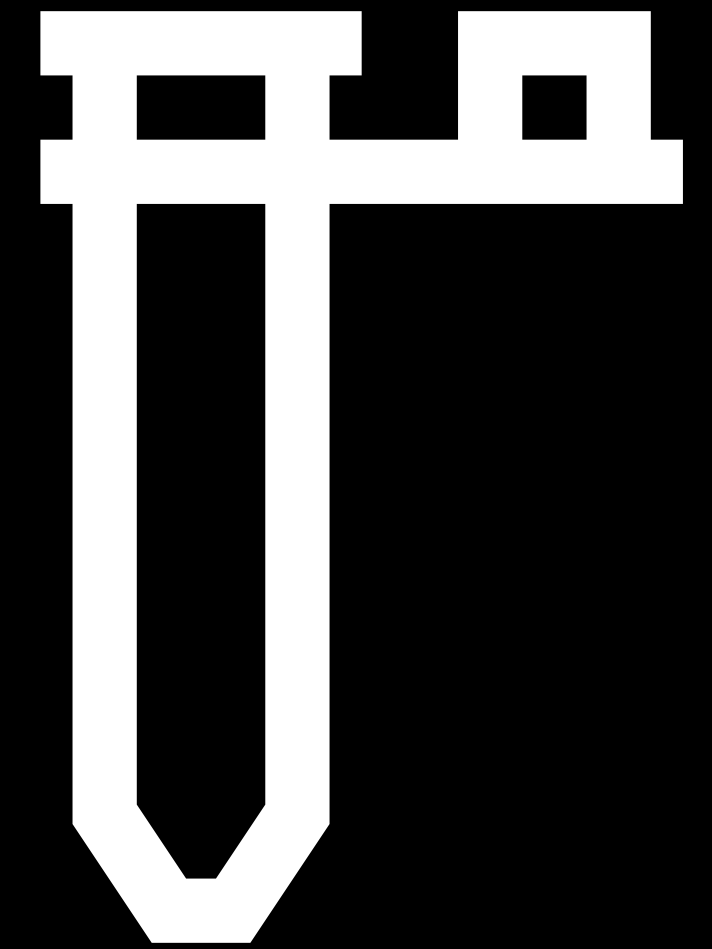
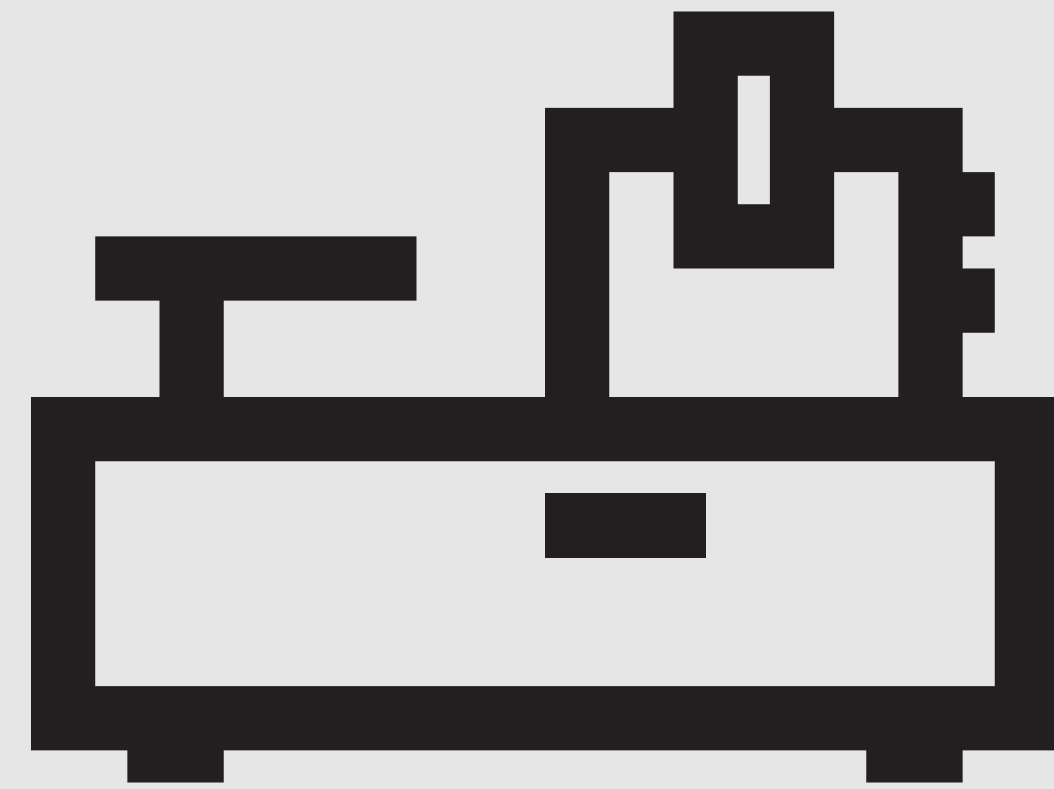


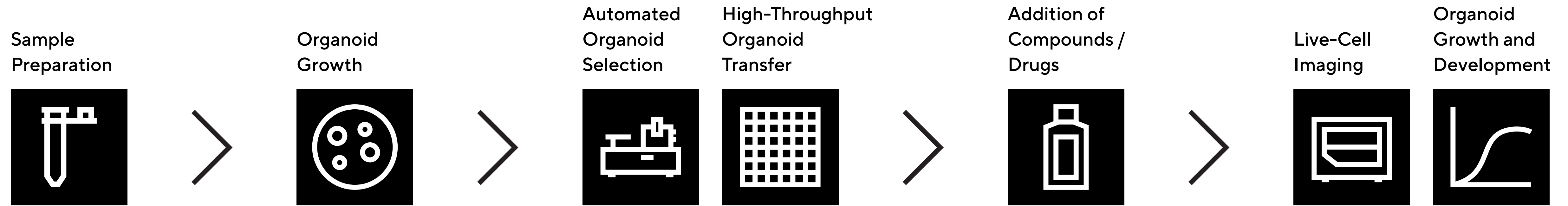
SARTORIUS

Simplifying Progress

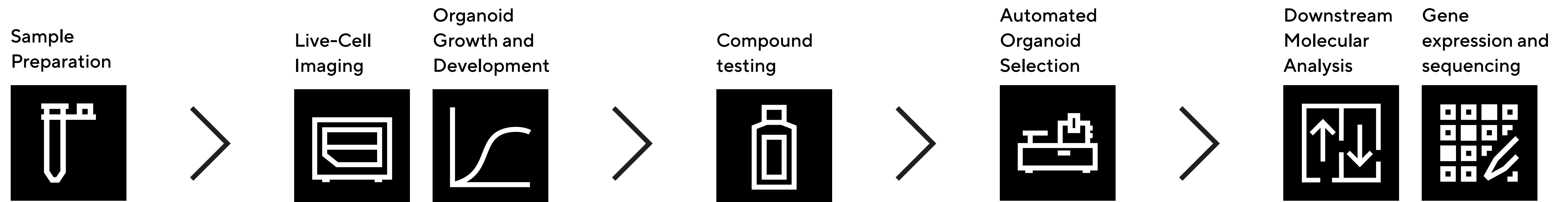
Advanced
Cell Model
Workflow



Live-Cell Imaging as the End-Point Assay



Molecular Analysis as the End-Point Assay



Germany

Sartorius Lab Instruments GmbH & Co. KG
Otto-Brenner-Strasse 20
37079 Goettingen
Phone +49 551 308 0

USA

Sartorius Corporation
565 Johnson Avenue
Bohemia, NY 11716
Phone +1 631 254 4249
Toll-free +1 800 635 2906



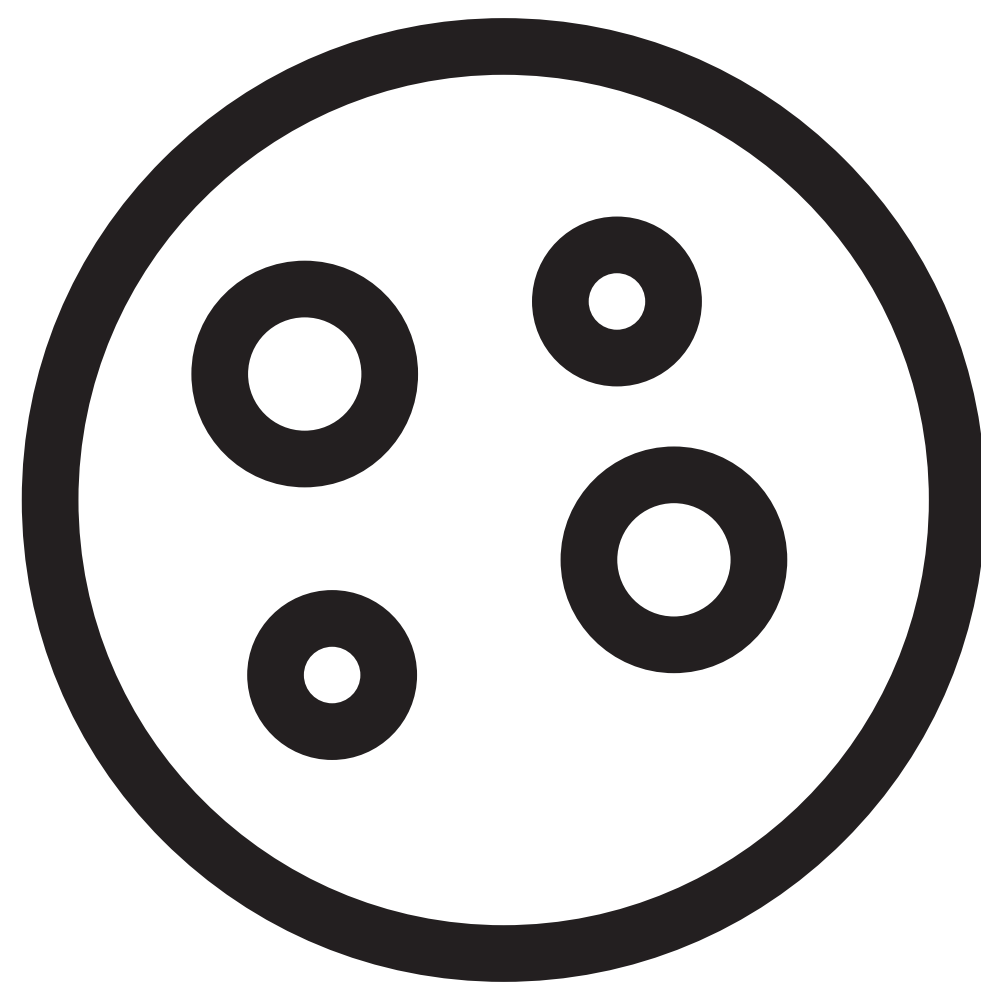
For further information, visit
www.sartorius.com

Specifications subject to change without notice.

© 2023 Copyright Sartorius Lab Instruments GmbH & Co. KG, Otto-Brenner-Strasse 20, 37079 Goettingen, Germany

Status: 08 | 2023

Live-Cell Imaging as the End-Point Assay



Organoid Growth

Any plates can be used, from Petri dishes to specially designed 6 well or 24 well round bottom Nanowell plates for Organoid growth and identification

Solutions include:

- [CellCelector Flex Capillaries and Consumables](#)

Sample Preparation

Live-Cell Imaging

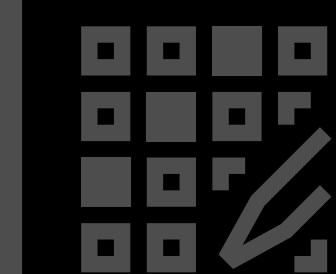
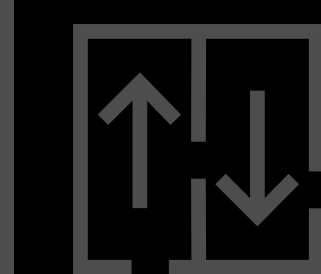
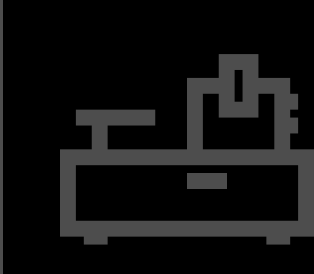
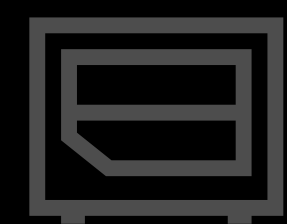
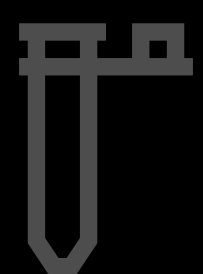
Growth and Development

Compound testing

Organoid Transfer

Molecular Analysis

Expression and sequencing



Live-Cell Imaging as the End-Point Assay



Automated Organoid Selection

Gently transfer organoids into individual wells without disturbing morphology and maintaining viability

Solutions include:

- [Cell Selection and Retrieval | CellCelector | Sartorius](#)

Sample Preparation

Live-Cell Imaging

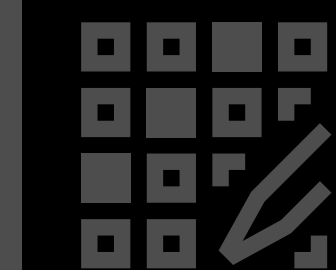
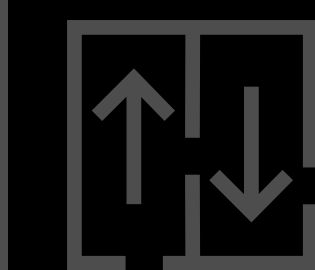
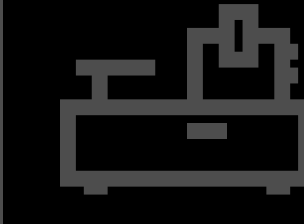
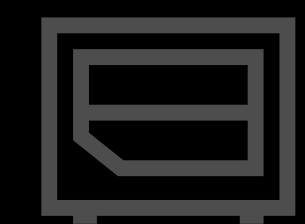
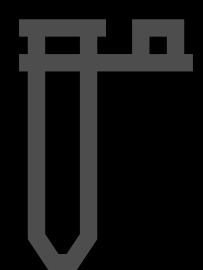
Growth and Development

Compound testing

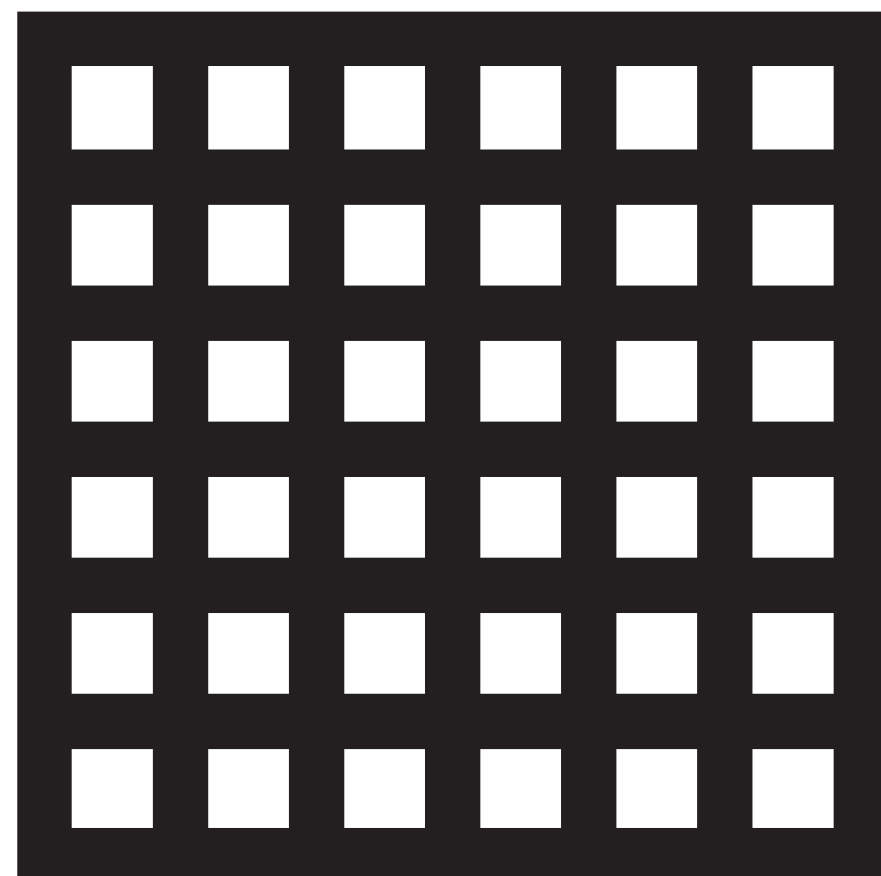
Organoid Transfer

Molecular Analysis

Expression and sequencing



Live-Cell Imaging as the End-Point Assay



High-Throughput Organoid Transfer

High-throughput automated transfer of organoids into 96 or 384 well plates with bubble-free Hydrogel deposition

Solutions include:

- [Organoid Technical Note | Sartorius](#)

Sample Preparation

Live-Cell Imaging

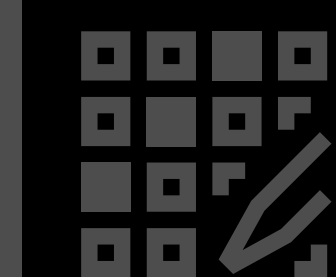
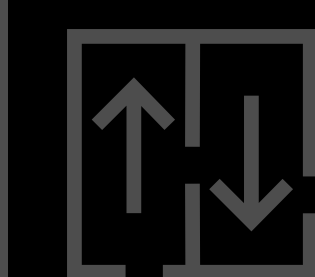
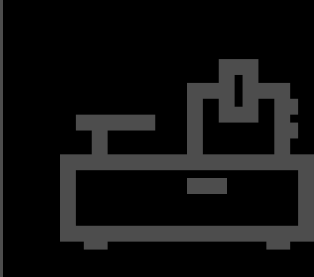
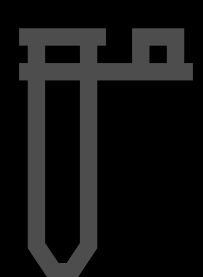
Growth and Development

Compound testing

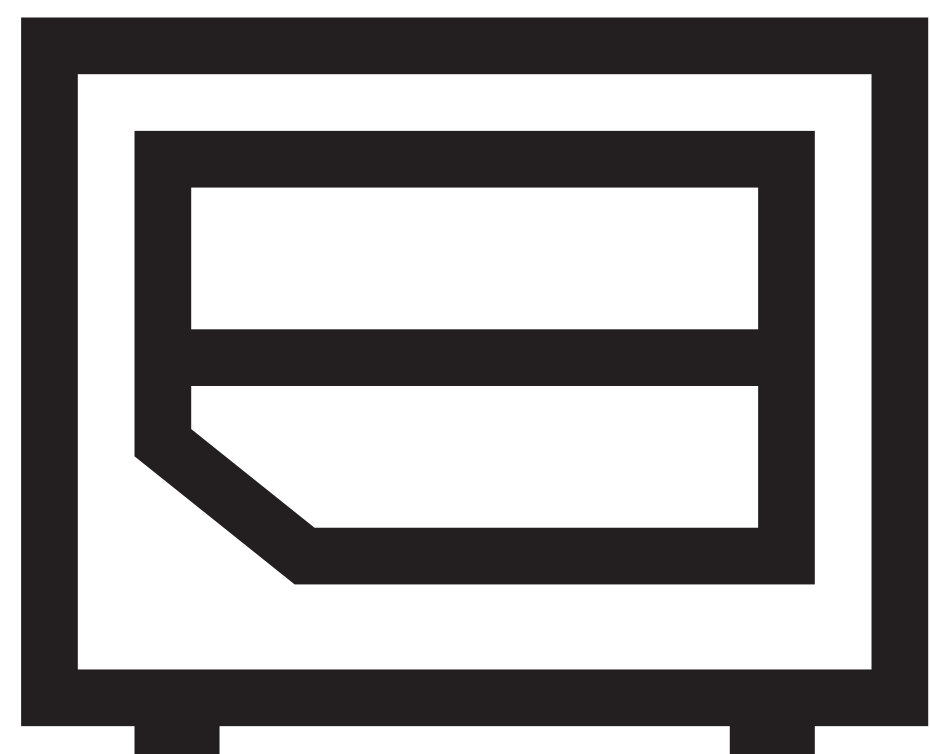
Organoid Transfer

Molecular Analysis

Expression and sequencing



Live-Cell Imaging as the End-Point Assay



Live-Cell Imaging

Analyze cells around the clock within the stable environment of the tissue culture incubator

Solutions include:

- [Organoid Assay | Sartorius](#)

Sample Preparation

Live-Cell Imaging

Growth and Development

Compound testing

Organoid Transfer

Molecular Analysis

Expression and sequencing

