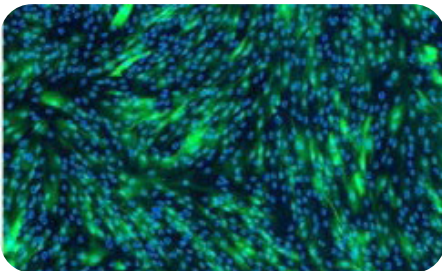


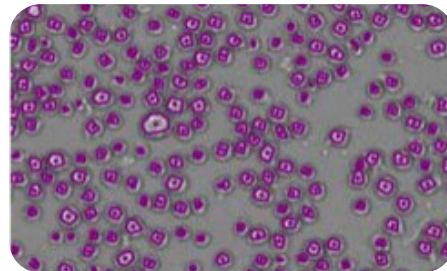


Cell-based screening with the Celigo image cytometer.

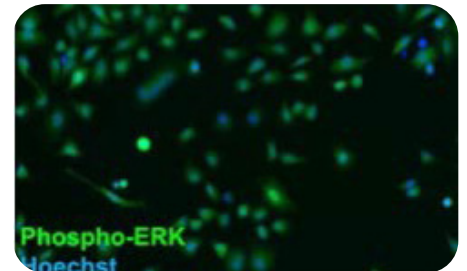
The Celigo™ image cytometer enables rapid, high quality, in situ, whole-well brightfield imaging for accurate label-free cell analysis. In addition, three channel fluorescence can be combined with brightfield for quantitative analysis of multiplexed assays. The Celigo is easy-to-use and offers a flow cytometry-like gating interface for optimal analysis of many different adherent and non-adherent cell types in multi-well plates (1536-well to 6-well) and T-flasks (T-25, T-75).



| High-throughput screening



| Cell line characterization



| Cell signaling

Assay development

- Straightforward cell segmentation software
- Flow cytometry-like gating interface with Boolean operators for cell classification
- Accurate normalization of wells using the actual number of cells in each well

Cell line characterization

- Determine growth characteristics of cells in situ, directly where they are grown
- Report growth curves, cell counts, confluence, doubling time, and doubling rate for each well
- Quickly analyze cells growing in T-flasks, T-75 flasks can be analyzed in ~15 min

Cell signaling

- Analysis of phosphorylated proteins associated with cell signaling responses
- Very little optimization required using “off the shelf” validated assays

High-throughput screening

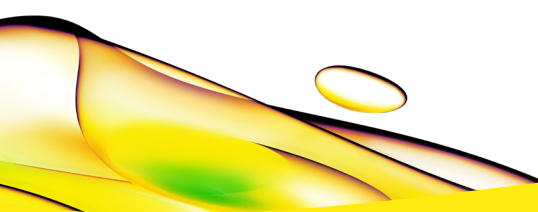
- Rapid scanning and analysis of multi-well plates (384-well plate can be analyzed in <5 min)
- Scan and analyze a 96-well plate using 2 fluorescent channels in <6 min
- Easy-to-use interface with specific applications that streamline image acquisition to data reporting

Proliferation assay

- Accurate, whole-well brightfield imaging and segmentation counts every cell in every well
- Label-free analysis allows multiple reads of the same sample, reducing cost and effort

Transfection & transduction optimization

- Quickly identify optimal parameters for high-efficiency transfection
- Determine transient and stable transfection rates and evaluate antibiotic induction using live imaging



revvity