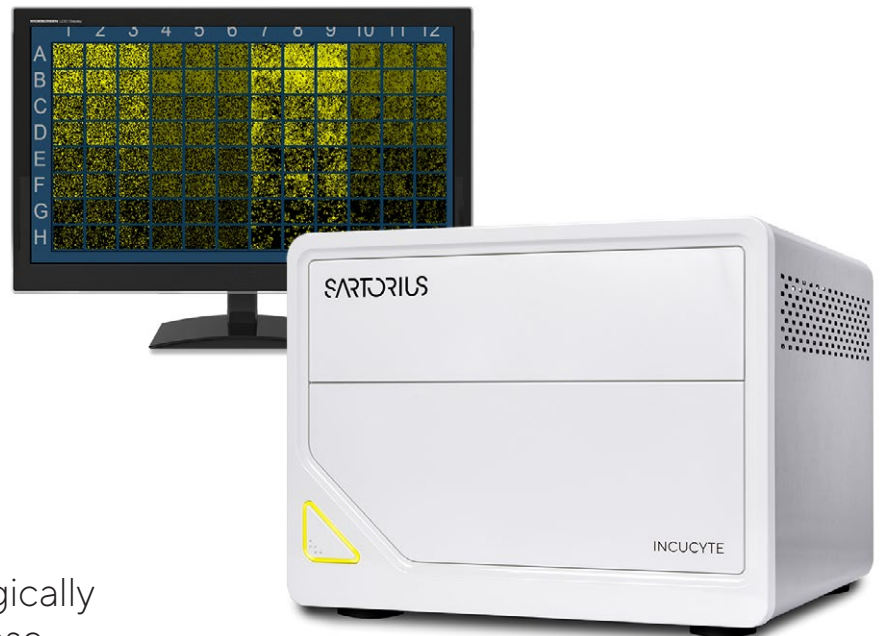


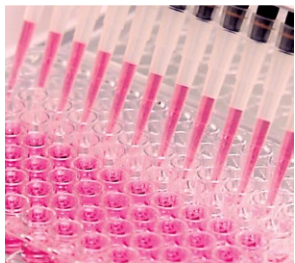
Incucyte® ATP Assay Gain Unprecedented Access to Metabolic Changes

With the Incucyte ATP Assay,
You Can...

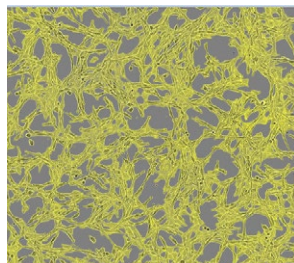
- Express a novel, genetically encoded fluorescent ATP indicator
- Analyze cell-type specific ATP changes in advanced cell models
- Conduct temporal studies of ATP for greater biological insight
- Inspect changes in cell morphology with HD phase images at every time point



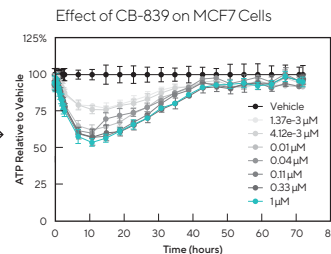
Measure and Visualize Dynamic
Metabolic Changes in a Physiologically
Relevant Environment with Purpose-
Built Software Tools, Novel Reagents,
and Guided Workflow.



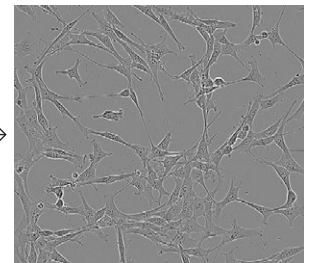
Infect cells with Incucyte
CytoATP Lentivirus



Capture images
over time



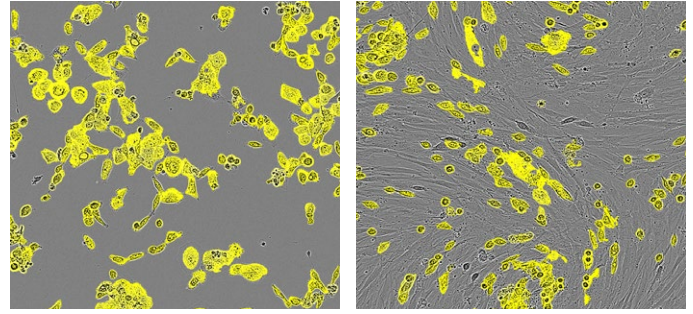
Quantify changes with
integrated kinetic analysis



Associate changes in
ATP metabolism with
cell morphology

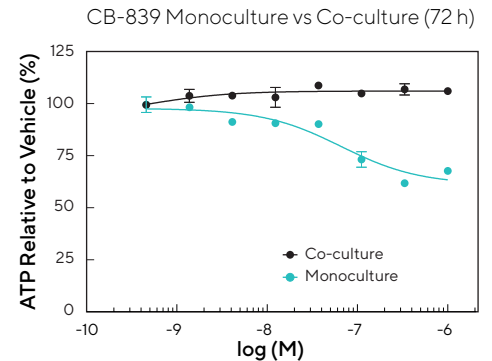
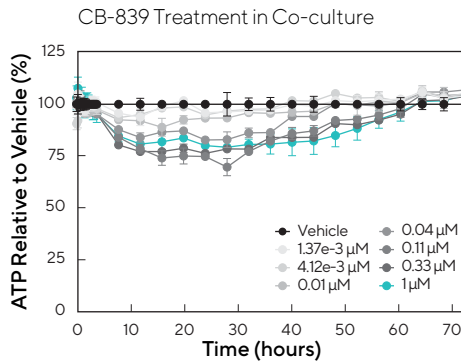
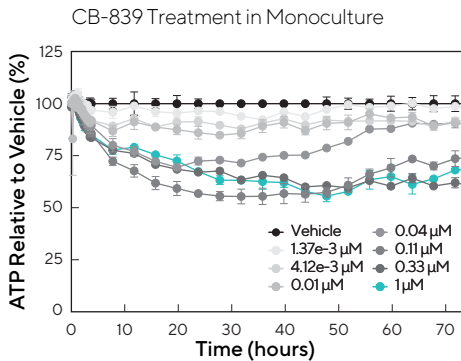
Distinguish Cell Type-Specific Metabolic Changes With Direct, Kinetic Measurements of ATP

- Reveal dynamic changes of tumor environment through analysis of cytosolic ATP and integrated normalization
- Maximize your metabolic insight with fully automated kinetic analysis of the same population of cells in a physiologically relevant environment



Monoculture

Co-culture



The triple negative breast cancer (TNBC) cell line HCC1806 stably expressing CytoATP was seeded in the presence or absence of CCD14086SK fibroblasts. Identification of cellular changes in ATP (color scaled masking) is performed using the integrated Incucyte® ATP Analysis Software Module. Masked images provide visualization of ATP in HCC1806 cells treated with 1 μM CB-839 in monoculture and co-culture with CCD14086SK cells. Data shows a sustained depletion of ATP in monoculture following CB-839 treatment, however, co-culture with stromal cells mediated resistance to CB-839.

Find out more:

www.sartorius.com/incucyte

Contact us: orders.US07@sartorius.com

North America: +1 734 769 1600, ext. 3

Europe: +44 7515 947101

Japan: +81 3 5826 4795

China: +86 21 6878 2300

Rest of APAC and other countries
around the world: +65 6872 3966