

OMNI

Live-cell imaging and analysis
from your incubator



DISCOVER

Omni

The Omni live-cell imaging systems are designed for continuous multiwell imaging in your incubator. Through intuitive design and robust image acquisition and analysis, the systems accommodate multiple applications and users. With automated brightfield and fluorescent imaging, you can study **complex biological processes** with ease.



*Track every cell and every moment.
Do it all with **Omni.***

VERSATILITY

Single platform, many uses

From cell culture monitoring to complex, multistep assays, the Omni platform is developed to accelerate research and enhance your cell culture workflow.

Incubator-based, *whole-plate imaging*

See every cell without disturbing the biology.

Monitor one plate *or many*

Choose your throughput with single-plate or automated multiplate handling.

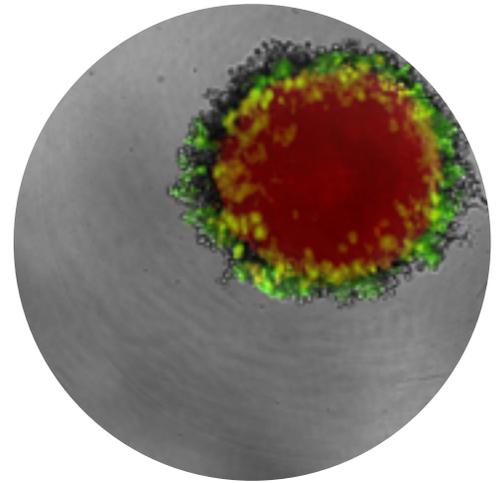
AI-driven *analysis*

Quantify dynamic cellular processes in real time.



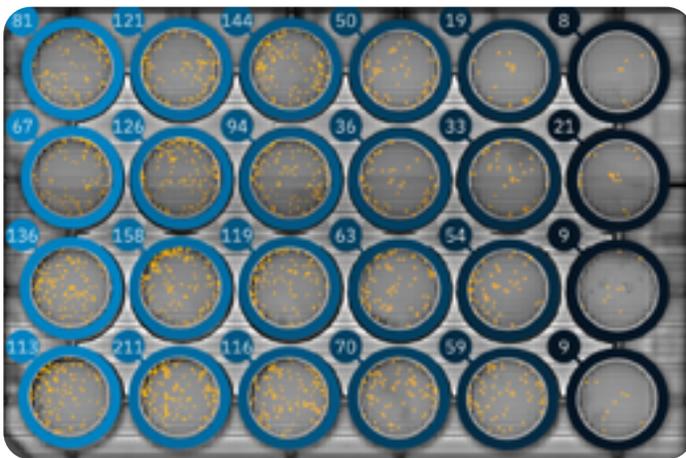
SEE MORE

Cellular biology up close

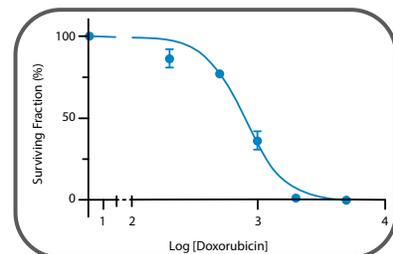
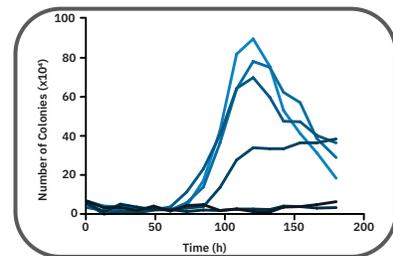


Live-cell imaging offers noninvasive analysis of cell health, morphology, and function. Generate more accurate and reproducible data using whole-plate brightfield imaging and fluorescence-based assays.

Image and analyze the whole plate



Colony formation of cancer cells exposed to increasing concentrations of doxorubicin (left to right). Omni tracks colony growth over time (right-top) and quantifies the dose response (right-bottom).

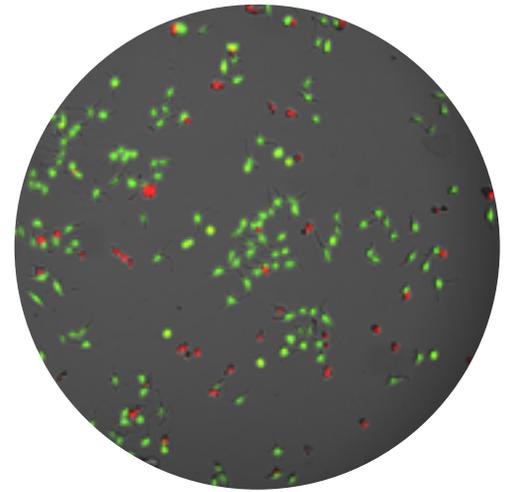


Measuring the whole well captures every cell, and AI-driven, machine-learning algorithms quickly analyze the results for easy, label-free brightfield assays. Even at low densities, like in this clonogenic assay, whole-well imaging ensures consistent, quantifiable results.

DO MORE

With simple assays

Red and green fluorescent channels make the Omni compatible with a wide range of assays. Here, the interactions between immune cells (red) and cancer cells (green) are monitored over time.



See what it can bring to your cell culture workflow:

- Cell health and proliferation
- Immune cell killing
- Cytotoxicity
- Tumor spheroids and organoids
- Cell migration
- Immune cell activation
- Colony detection

To find out more:

Visit axionbio.com/products/imaging/omni
or scan the QR code



Choose the **OMNI** for you

Omni BR

Omni Pro 12

Omni FL



<i>Features</i>	<i>Omni BR</i>	<i>Omni Pro 12</i>	<i>Omni FL</i>
Whole-well brightfield	✓	✓	✓
Automated acquisition	✓	✓	✓
Incubator compatible	✓	✓	✓
Fluorescence (Red)		✓	✓
Fluorescence (Green)		✓	✓
Number of plates	1	12	1
Plate handling	Manual	Automated	Manual

SOFTWARE MODULES

To meet your needs



Clonogenic assay *module*

Track colony formation for clonogenic assays



Confluency *module*

Track confluence to monitor cell proliferation or death



Fluorescent object *count*

Quantify and track fluorescent objects



Organoid analysis *module*

Identify, track, and analyze organoids



Scratch assay *module*

Measure wound healing and cell migration/invasion

Learn more:

axionbiosystems.com/products/imaging/omni

Contact us:

axionbiosystems.com/contact

Office locations:

North America - Europe - Asia Pacific

For pricing and ordering:

sales@axionbio.com or scan the QR code



For research use only. Not for use in diagnostic procedures.
This information is subject to change without notice. © Axion BioSystems, Inc, 2023.