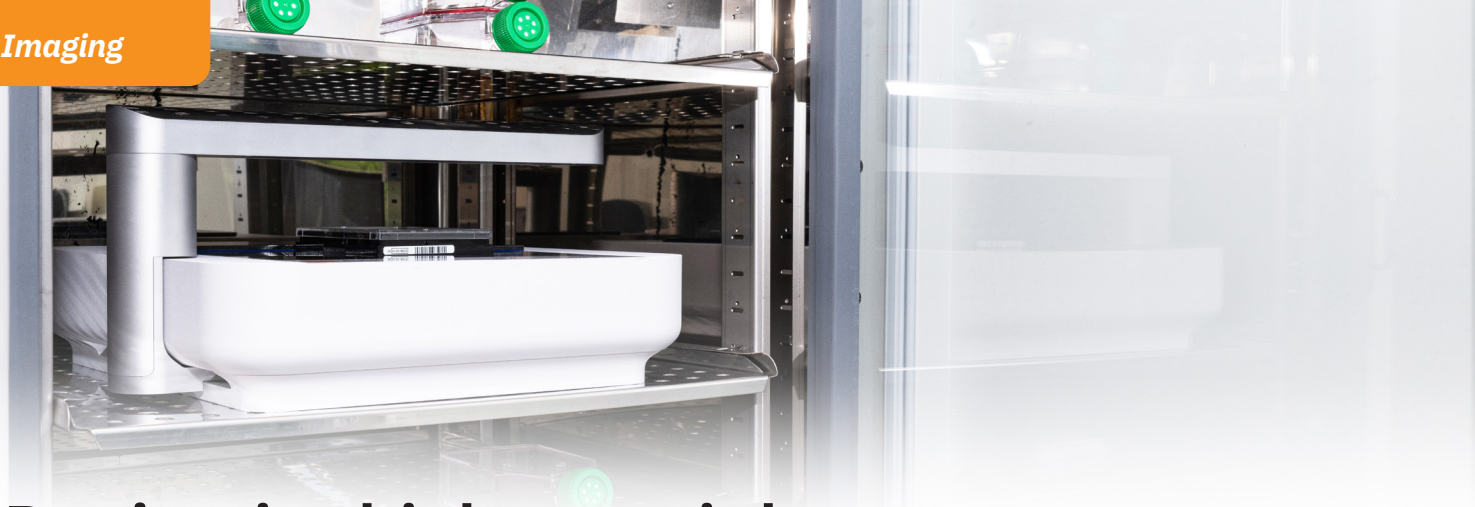


# Omni Platform

Powerful live-cell imaging.  
Seamless AI analysis.



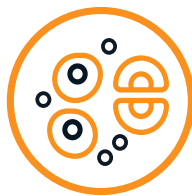


## Don't miss biology as it happens

Live-cell imaging has revolutionized cell biology research, making label-free, kinetic assays accessible. **With the Omni live-cell imaging platform you get:**



**Complete, realtime datasets**



**Physiologically relevant assays**



**Experimental flexibility**

## Omni: An essential tool for every lab

Visualize complex cellular biology with ease using Omni. Brightfield and fluorescence imaging combined with advanced software tools and intuitive design make the Omni an **essential tool for every laboratory.**

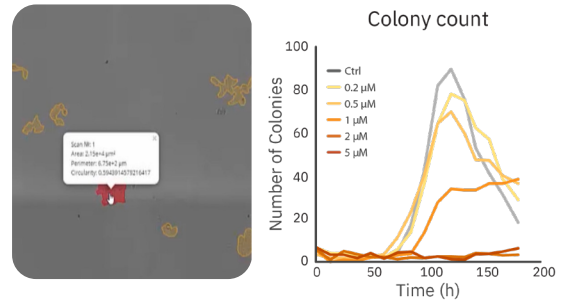
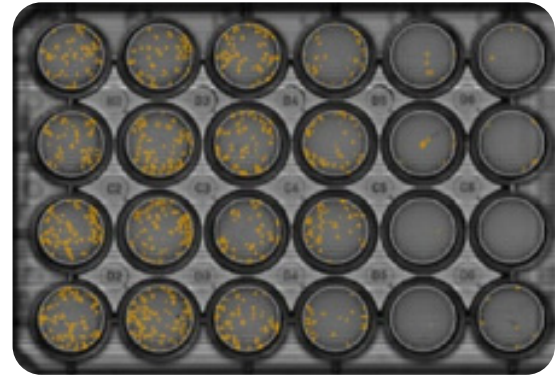
### *Key features*

- **Whole-vessel** brightfield imaging
- **Red & green** fluorescence
- **Powerful**, intuitive software
- **Incubator-friendly** design
- **AI analysis** with cloud-based computing



# The power of live-cell imaging + AI analysis

The Omni combines **real-time cellular imaging and fast, AI-powered analysis** into one imaging platform for labeled and label-free kinetic assays. Monitor growth across the entire plate with brightfield or evaluate specific mechanisms or markers with its fluorescence capabilities.

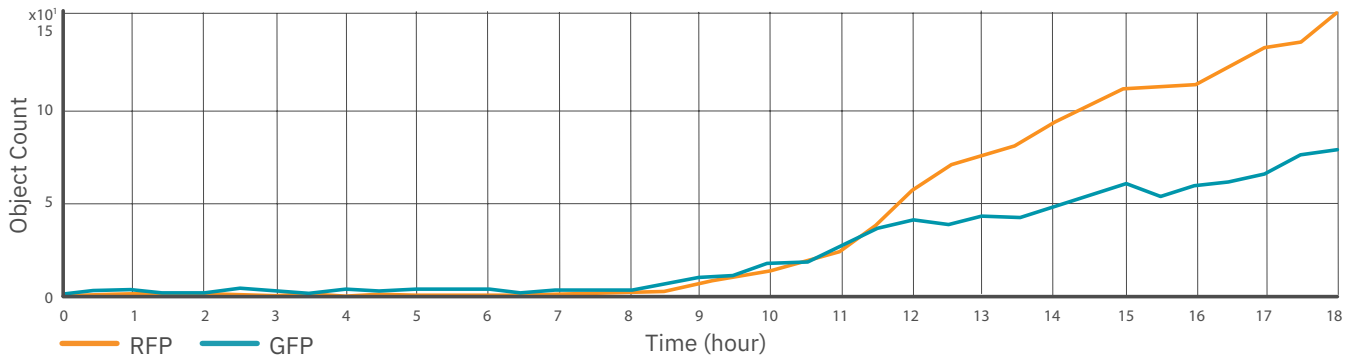
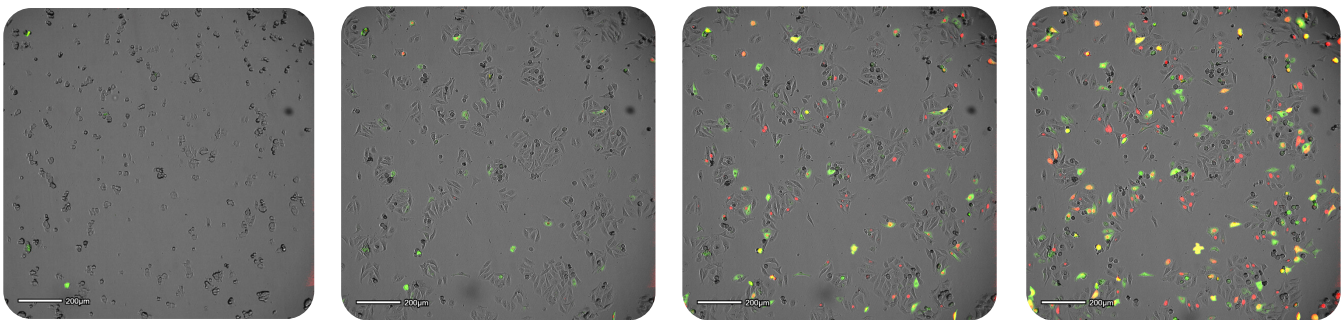


## Image and analyze the entire plate

**Whole-vessel brightfield imaging** gives unbiased results even in low density cultures found in clonogenic assays.

## Track cultures over time

**Advanced AI analysis easily quantifies every time point.** Here expression of cells transduced with actin GFP and nucleus RFP using BacMam is tracked over time using fluorescence.

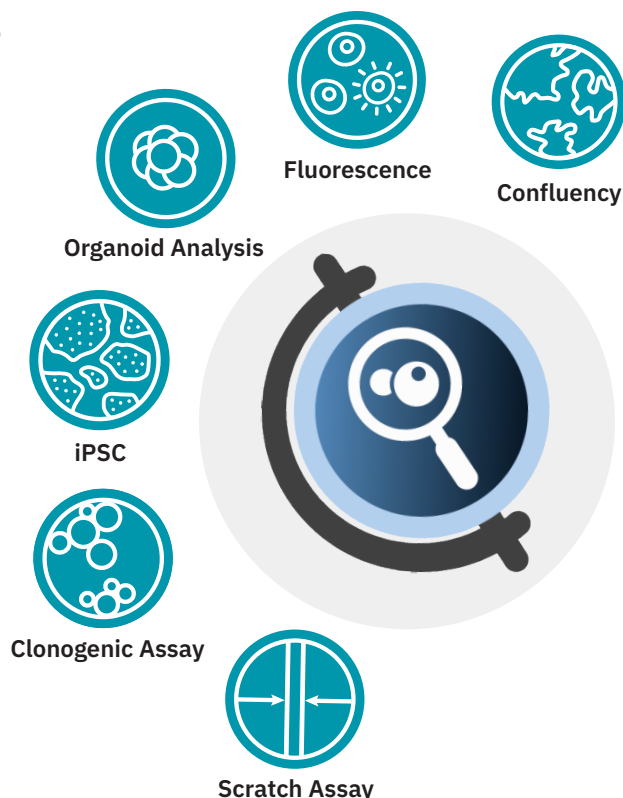


# Versatile and easy to use software

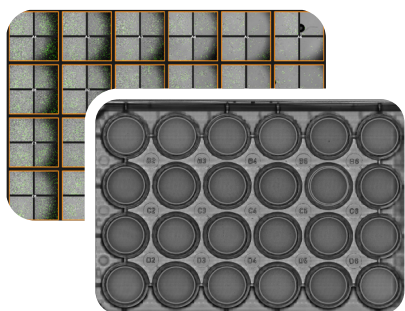
Omni's AxIS Vue software provides the flexibility to customize the platform to suit your research needs. Simply choose the modules you want to enable a wide range of applications:

## Applications

- Cell proliferation
- Cell migration
- Colony formation
- Cytotoxicity and cell viability
- Transfection and transduction
- Microfluidic chips
- Stem cell monitoring
- Spheroids & organoids

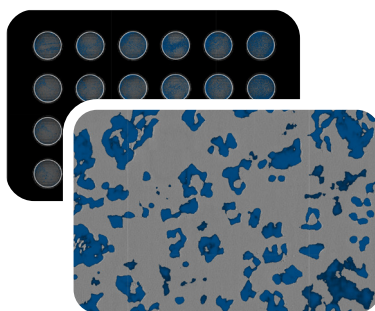


# Leveraging the cloud for fast, easy analysis



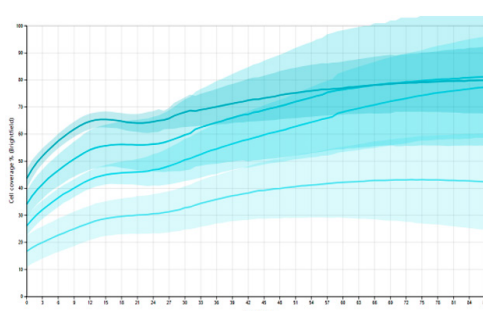
### >> View every moment

Rapidly takes 7850 images and combines them so you can **view the whole plate or zoom to any point of interest** at any time.



### >> Analyze with AI

**Processing the data is fast**, leveraging state-of-the-art AI detection and the computation power of the cloud.



### >> Get instant results

**Conveniently add plate maps and instantly view your results.** Export images and videos with a few clicks.



# The Axion Portal: A gateway to discovery

Imaging can generate large amounts of data requiring large amounts of computing power. The Axion Portal is a cloud-based computing environment for all Axion BioSystems imaging platforms. **Discover how it can accelerate your research:**



## Speed

Faster analysis with the processing power of the cloud



## Accessibility

Access your data remotely and easily collaborate



## Scalability

Save on infrastructure costs with unlimited data storage



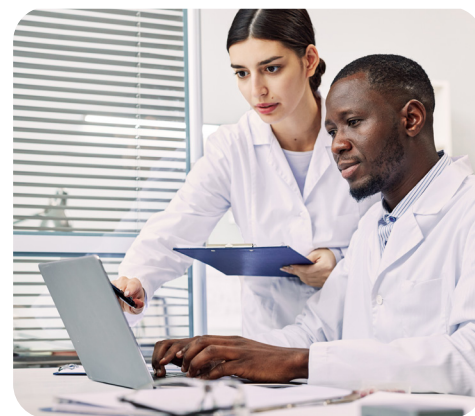
## Security

Keep your data safe with secure login and data encryption

## Working remotely, while working together

The Axion Portal makes it simple to analyze your data anywhere, anytime. Stay informed and make teamwork easy with convenient notifications and collaboration features.

- >> **Check your cells** *while away from the lab*
- >> **Analyze data** *with just a laptop*
- >> **Keep your data organized** *in one location*



## Keep your data safe

The Axion Portal was built to be secure, simple, and powerful. To ensure data security and integrity, the Axion Portal uses the Microsoft Azure cloud environment:



- >> **All data is fully encrypted.**
- >> **Data is never lost or accidentally deleted** *thanks to triple redundant storage.*
- >> **Your data is only accessible to you.** *You control who has access to your data with secure login.*

# Choose the *Imaging system for you*

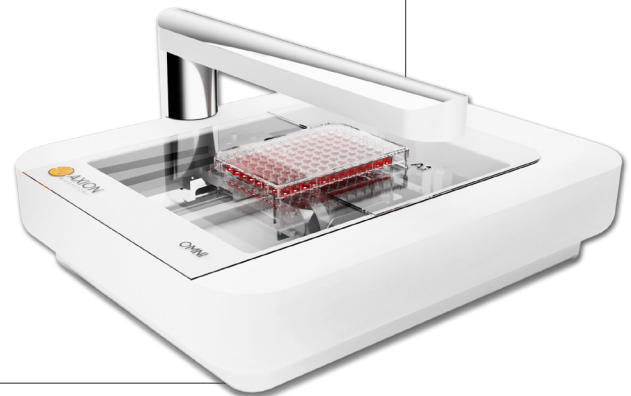
## **Omni Pro 12:** *Bringing imaging automation to your incubator*

The Omni Pro 12 is for labs that want more, featuring integrated robotics and **12-plate capacity for enhanced flexibility and efficiency**. Designed to provide an automated solution to accommodate multiple users and live-cell imaging applications.



## **Omni:** *Essential live-cell imaging for every lab*

The Omni imager adds dynamic visual results to any experiment. Featuring AI-supported analysis, compatibility with all analysis modules, and the versatility to scan all types of culture vessels, **do it all with Omni**.



## **Lux3:** *Convenient and affordable live-cell imaging*

**A compact live-cell analysis platform** with a single field of view designed to help scientists explore the kinetics of cell proliferation, cell migration, cell morphology, and more, while on a budget.



# Imaging system features comparison

There's a live-cell imaging solution for every lab. Choose the device that works best for you.



<i>Features</i>	<i>Lux3 (BR/FL)</i>	<i>Omni</i>	<i>Omni Pro 12</i>
Incubator compatible	✓	✓	✓
Whole-plate brightfield		✓	✓
Fluorescence Green Ex. 452/45 nm Em.512/23 nm Red: Ex. 549/15 nm Em. 630/90nm	(FL only)	✓	✓
Magnification	10X (20X digital)	10X	10X
Plate Handling	Manual	Manual	12-plate automated
# of Plates	Single view	1	12
Software	Confluency, Scratch, Fluorescence	All	All
Dimensions & weight	166 x 140 x 135 mm 1.3 kg	396 x 345 x 171mm 9.7 kg	460 x 417 x 439 mm 40.2 kg



Want to learn more?

Visit **axionbio.com** for more data, applications, and resources.

## Our commitment to our customers

With over **15 years of experience** bringing innovative new products to our customers, we strive to accelerate your research by making live, functional biology more accessible. Our design philosophy is to ensure all of our products are:



### Flexible

Hardware designed for broad, integrated functionality in one instrument



### Easy to use

Intuitive instruments, consumables, and software for fast, easy adoption



### Smart technology

Easy to run with no complicated steps, saves time and money

Contact our scientists to discuss how we can help your research with live-cell imaging:  
[axionbiosystems.com/contact](https://axionbiosystems.com/contact)