



Accelerate drug discovery: end-to-end imaging solutions to help support regulatory confidence.

From organoids to living models, Revvity's imaging solutions help provide researchers with a comprehensive picture to support the path toward safer, more effective therapeutics.

Modern drug discovery requires more than isolated data points. Researchers need integrated insights across biological complexity, from advanced 3D organoid models to whole-animal *in vivo* imaging.

Revvity delivers both high-content cellular imaging and *in vivo* imaging solutions within one connected ecosystem, helping researchers improve translational relevance, accelerate toxicity assessment, and make more confident decisions.

Why Revvity

- Integrated imaging portfolio from *in vitro* to *in vivo*
- Improved translational confidence
- Scalable workflows for drug discovery
- Trusted by leading pharma and biotech researchers
- Advanced analytics and automation capabilities



Organoid and cellular analysis

- High-content imaging
- Phenotypic screening
- Multiparametric analysis
- AI-enabled image analysis

Translational validation

- 3D disease models
- Functional assays
- Longitudinal studies
- Predictive toxicity insights

In vivo imaging

- Real-time monitoring
- Whole-body imaging
- Non-invasive analysis
- Faster evaluation of drug candidates

Organoid and cellular analysis



Opera Phenix OptIQ™ and Operetta CLS™ high-content screening and analysis platforms allow understanding the pathogenesis of viral infection and the host immune response, and support compound screens for drug discovery or repurposing as well as toxicity studies for drug development.



Phenologic.AI

The Phenologic.AI™ module in Harmony™ and Image Artist™ software harnesses the power of pretrained deep neural networks (DNNs) to provide an efficient and reliable method for identifying cells and cellular nuclei within fluorescent and brightfield images.

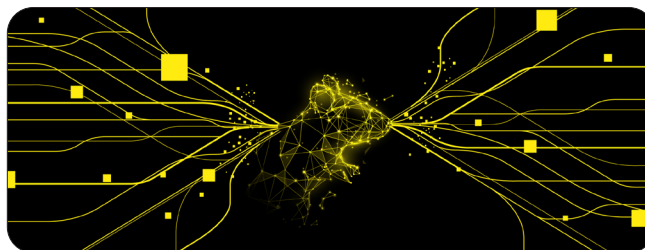
In vivo imaging solutions



Optical Imaging

Small animal *in vivo* imaging on the IVIS™ system allows studying tumor biology, infectious disease, drug biodistribution, safety, and efficacy *in vivo*. Animals can be studied repeatedly over time, thus limiting the number of animals needed and allowing to understand therapy or vaccination effectively over time.

In vivo imaging reagents - *In vivo* imaging reagents enable visualization of important biomarkers *in vivo*, such as those involved in lung inflammation and others.



Living image Synergy AI

Living Image™ Synergy AI multimodality software is our next generation platform for *in vivo* imaging analysis.

As a complement to existing Living Image, SonoEQ, or microCT software, Living Image Synergy AI offers imaging researchers a unified platform with AI capability to support seamless data analysis across optical, microCT, and ultrasound imaging modalities.

Revvity imaging ecosystem

Research stage	Solutions	Key features
Cellular analysis	Operetta CLS system	Fast, reliable phenotypic analysis
Advanced 3D imaging	Opera Phenix OptIQ system	Deep organoid imaging with high sensitivity
AI image analysis	Harmony, Image Artist, Phenologic.AI software, Living Imaging Synergy AI	Accelerated image interpretation
<i>In vivo</i> validation	IVIS imaging systems	Non-invasive longitudinal monitoring

Reach out to your local sales representative or visit www.revvity.com for more information.