

# Powerful yet simple high-content image analysis.

### Harmony high-content imaging and analysis software

Harmony<sup>™</sup> high-content imaging and analysis software streamlines your entire workflow, from acquisition to results evaluation. It allows for quantification of complex 3D cellular models, reliable phenotype discrimination, and efficient data interpretation. With its intuitive interface, Harmony guides you through every step of your experiment on your Opera Phenix<sup>™</sup> Plus or Operetta CLS<sup>™</sup> experiment, providing instant feedback and comprehensive control, ultimately transforming your biological data into meaningful discoveries.

- Analyze common assays with more than 30 ready-made solutions, or create your own with simple image-analysis building blocks.
- Easily quantify complex cellular phenotypes based on changes in morphology, fluorescence intensity, intensity distribution, and texture parameters.
- Follow phenotypic changes over time and accurately quantify label-free live cell images.
- Visualize and analyze your samples in 3D for greater depth of information and insights in a more physiologically relevant context.

- Phenotypic screening
- Live cell imaging
- 3D Cell models
- Rare cell phenotypes





- A. Workflow-based interface with easy-to-read icons
- **B.** Analysis building blocks for easy protocol design
- **C.** Clear plate navigation and wizard for easy setup of new plate types
- **D.** Results summary with immediate numerical output for faster insights

Ready-made analysis templates

- Cell counting
- Cell painting
- CytotoxicityNuclear translocation
- Receptor internalization
- Morphology

- Spot analysis
- Neurite outgrowth
- Micronucleus
- Lipid droplet
- Colony formation
- Microtissue analysis
- Migration
- Confluency
- Cell tracking
- Fluorescence distribution
- Cell cycle classification
- ... plus you can build your own.

## New in Harmony 5.3

**Harmony 5.3** simplifies data management for faster image data handling and offers intuitive, powerful search functionality. Capabilities include:

- Faster data transfer: Speed up your data transfer processes, reducing wait times.
- Seamless experiment setup: Move data easily as part of your experiment setup, enhancing your workflow efficiency.
- **Scalable storage**: Distribute storage for enhanced scalability, accommodating growing data needs.
- **Intuitive search experience**: Enjoy an easy and efficient search experience that saves you time when identifying your data.
- Efficient multi-location search: Quickly search across multiple storage locations, improving data retrieval.
- **Open data format**: Organize data in files and folders with humanreadable names for easy access.
- **Rapid analysis with HPC**: Leverage high-performance computing (HPC) to analyze large datasets quickly, accelerating your research timelines through integration with Signals Image Artist.

# New Phenologic.AI 1.0

**Phenologic.AI™**, a module in Harmony and Signals Image Artist<sup>™</sup> software, harnesses the power of pre-trained deep neural networks (DNNs) to provide an efficient and reliable method for identifying cells and cellular nuclei within fluorescent and brightfield images.

- **Turnkey Al image analysis**: Utilizes pre-trained Al for easy and efficient analysis of both fluorescent and brightfield images.
- Label-free detection: Simplifies detection of cellular nuclei in brightfield images without nuclear staining, saving time in sample preparation.
- **Robust identification**: Trained on diverse cell lines, it reliably identifies cellular structures and provides phase contrast imaging.
- **Increase multiplexing capabilities**: Enhance your experimental flexibility and use more fluorescent dyes for other markers.
- **Improved viability**: Lower phototoxicity for live-cell experiments through avoiding stress from fluorescent dyes, improving cell viability.

### www.revvity.com Copyright ©2025, Revvity, Inc. All rights reserved. Revvity is a trademark of Revvity, Inc. All other trademarks are the property of their respective owners.