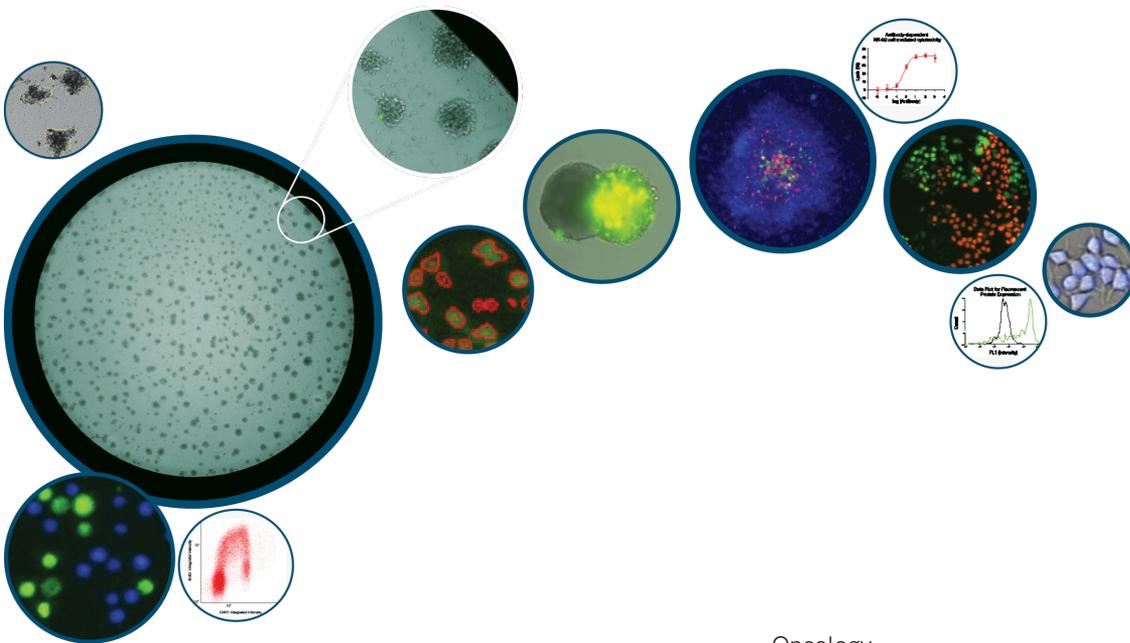
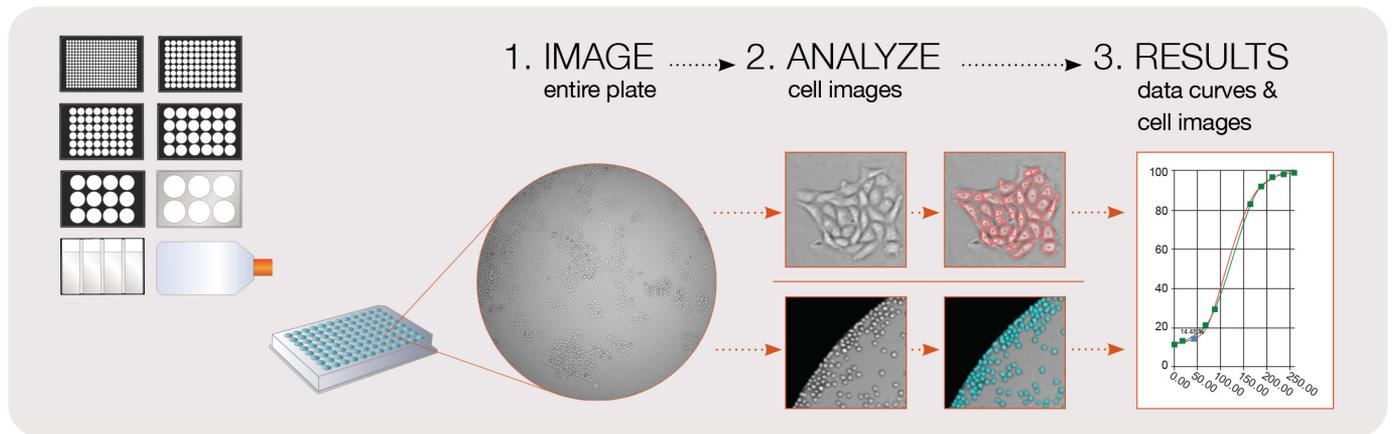


revvity

# Integrated cell imaging and analysis.



- Oncology
- Immuno-Oncology
- Cell Line Development and Biologics Production
- 3D Tumor Spheroid Modeling
- Stem Cell Research
- Virology Assays



## Oncology

- Apoptosis
- Cell cycle
- Cell viability
- DNA synthesis
- Cytotoxicity
- Cell migration
- Cell morphology
- Immunophenotyping
- Metabolism
- Wound healing
- Chemotaxis

## Stem cell research

- iPS cell line generation
- Embryoid body morphology
- Stem cell marker analysis

## Cell line development and biologics production

- Single cell detection
- Transfection/Transduction efficiency
- CRISPR fluorescent reporter monitoring
- Single cell to single colony monitoring
- Growth tracking
- Cell secretion
- Hybridoma screening
- Routine quality control

## Virology assays

- Foci and plaque counting
- Antibody neutralization
- Hemagglutination assay

## 3D tumor spheroid modeling

- Growth tracking
- Growth inhibition
- Apoptosis/Viability
- Invasion/Migration
- Patient Derived Organoids (PDOs)

## Immuno-oncology

- Direct cell-mediated cytotoxicity
- Antibody-Dependent Cell-mediated Cytotoxicity (ADCC)
- Complement-dependent cytotoxicity
- Antibody-mediated cytotoxicity

Plate Compatibility	6, 12, 24, 48, 96, 384, 1536 well plates (black, white and clear well plates) / T-25 and T-75 flasks. Slides and cell array plate profiles available upon request. Holders/adapters for 60 mm dish, and slide holders that fit up to 4 slides also available.				
Software	Proprietary image acquisition and processing software Powerful analysis software/Computer workstation Microsoft Windows				
Illumination/Optics	<ul style="list-style-type: none"> <li>1 LED-based enhanced bright field imaging channel with uniform well illumination</li> <li>4 LED-based fluorescent channels</li> <li>Proprietary F-theta lens with superior well edge-to-edge contrast</li> <li>Galvanometric mirrors for fast imaging of large areas</li> <li>Large chip CCD camera (2024x2024 pixels)</li> <li>1, 2, 4 or 8 um/pixel resolution</li> </ul>				
Fluorescent Channels	Channel	Excitation	Dichroic	Emission	Example Dyes
	Blue	377/50	409	470/22	Hoechst, DAPI
	Green	483/32	506	536/40	FITC, Calcein, GFP, Alexa Fluor® 488
	Red	531/40	593	629/53	R-PE, PI, Texas Red, Alexa Fluor® 568
	Far-Red	628/40	660	688/31	DRAQ5®, Alexa Fluor® 647
High-Speed Imaging	Less than 2 minutes per 384-well plate				
Weight and Dimensions	Dimensions: 19.5" W x 16"H x 24"D (49.5cm x40cm x 61cm) Weight: 117 lbs. (53 kg)				
Focusing Modes	Hardware-Based Auto Focus, Image-Based Auto Focus, Manual Focus				

