



revvity

The nexus of speed and sensitivity.

EnVision® Nexus™

Our fastest, most sensitive plate reader ever

Operating on a brand-new innovative platform, our EnVision Nexus multimode plate reader fast tracks your research, delivering the speed and accuracy you need for your most demanding applications.

Fully automatable for walkaway convenience, the EnVision Nexus system provides the speed and sensitivity you need to boost your lab's productivity. It's flexible, configurable, and upgradable to adapt to your changing needs.

Combined with a lineup of exciting enhancements, this system elevates your science and takes your research to the next level.



For research use only. Not for use in diagnostic procedures.

Features and benefits at a glance

Modalities	EnVision Nexus 2-Detector	Standard features	Optional extras
Absorbance	✓	<ul style="list-style-type: none"> 5 filter-based detection technologies 2 detectors enabling simultaneous dual wavelength reading Top and bottom reading for FI, TRF and LUM Shaker with 3 modes: linear, orbital, and double orbital Plate ID barcode reader (short side only) Kaleido™ 4.0 instrument control software Desktop PC, including monitor Temperature control 	<ul style="list-style-type: none"> Ultrasensitive Luminescence Enhanced Alpha (with aperture) High-throughput Alpha TRF Laser Preconfigured, assay ready filter modules Plate stacker (20 or 50 plates) Fully automatable Data analysis software from MyAssays® Desktop Enhanced Security software supporting 21 CFR Part 11 compliance in regulated environments, with secure export to MyAssays® Desktop
Luminescence (LUM)	✓ (standard and enhanced)		
Fluorescence Intensity (FI)	✓		
Time-resolved Fluorescence (TRF)	✓ (lamp)		
Fluorescence Polarization	✓		
Alpha	Optional		

Enhance your lab with the new EnVision® Nexus™ platform

For more than twenty years, our EnVision multimode plate reader has been the proven leader in high-throughput screening - and the EnVision Nexus is the next generation of superior detection. This new platform includes:

- Dual detectors for highest speed measurements
- Optional Ultrasensitive Luminescence detection mode boosts sensitivity up to 25x compared to Enhanced Luminescence Two Alpha detection modes: high-throughput screening (HTS) module for highest speed, or enhanced module to enable AlphaPlex technology
- Optional laser-based TRF module for highest speed
- Bottom read mode for all standard technologies
- Standard barcode reading (left side)
- Preconfigured filter modules for optimized optics and simplified workflows
- Enhanced Security software package for regulated environments provides technological controls and features to support 21 CFR Part 11 compliance
- Plate stackers and robotic integration for HTS

For more information, visit www.revvy.com or contact your sales representative!

