



Precision meets practicality: One delivers.

EnVision Nexus™ One

Fast, sensitive detection on an innovative platform

Operating on the advanced EnVision Nexus platform, the EnVision Nexus One brings flagship detection performance within reach of more labs, from high-throughput screening centers to emerging research teams.

With enhanced sensitivity in TRF, Alpha, and Luminescence, its single-detector configuration delivers the performance to support your demanding applications. From hit identification to lead optimization, the system helps you accelerate your drug discovery workflow.

Field-upgradeable and automation-ready, the EnVision Nexus One integrates smoothly into your existing Revvity ecosystem, backed by our comprehensive portfolio of reagents, microplates, and expert application support.



Features and benefits at a glance

	EnVision Nexus One
Detector configuration	Single detector for sequential measurements
Standard detection technologies	Absorbance, fluorescence intensity*, time-resolved fluorescence*, fluorescence polarization, luminescence*, enhanced luminescence
Advanced detection technologies (optional)	Enhanced Alpha, HTS-Alpha, ultrasensitive-luminescence, TRF Laser
Standard accessories	Shaker with 3 modes, plate ID barcode reader, desktop PC and monitor, temperature control
Automation options	Stacker (20 or 50 plates), RCI for automation integration
Software	Kaleido™ Software for reader control, Enhanced Security software option supporting 21 CFR Part 11 compliance, optional MyAssays® Desktop for data analysis
Throughput	Suitable for medium-throughput labs
Suited for:	Labs requiring enhanced sensitivity and practical scalability

* Top and bottom read mode available

Enhance your lab with the new EnVision Nexus One

- Built on an advanced platform with more than 20 years of expertise, the EnVision Nexus One represents the next generation of multimode detection, supporting your evolving demands in high-throughput screening and drug discovery. Here’s what it means for your lab:
- **Detect weaker signals with reliability** – Enhanced sensitivity across detection technologies helps you capture the subtle biological responses
 - **Optimize your Alpha screening workflow** – Choose between high-throughput screening (HTS) module for high speed or Enhanced module for multiplexed AlphaPlex™ assays
 - **Speed up time-resolved fluorescence assays** – Optional high power TRF laser supports fast, high-performance measurements with improved signal-to-noise ratio
 - **Streamline sample tracking** – Integrated barcode reading helps eliminate manual entry errors and supports complete traceability
 - **Simplify method development** – Preconfigured filter modules with optimized optics reduce setup time and deliver consistent results
 - **Support regulatory compliance** – Enhanced Security software package provides audit trails, electronic signatures, and controls to help satisfy 21 CFR Part 11 compliance
 - **Scale with your growing needs** – Plate stackers and robotic integration help enable walkaway automation for high-throughput laboratories

For more information, visit www.revvity.com/EnVisionNexus or contact your sales representative!