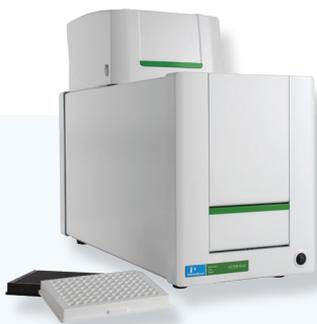


Multimode Microplate Readers for Every Application and Throughput Level



VICTOR® Nivo™ Industry's Smallest Multimode Plate Reader

- Compact and lightweight
- High performance for routine applications, including biochemical and cell-based assays
- Intuitive, browser-based software allows for remote operation
- All major detection modes including Alpha Technology
- Best for research labs with multiple users and relatively low throughput



EnVision® Our Fastest Benchtop Multimode Plate Reader

- Proven performance when the highest levels of speed and sensitivity are required
- Uses two detectors to enable simultaneous, dual-wavelength reading
- Well established for laser-based Alpha and TRF technologies, as well as ultrasensitive luminescence
- Best for screening and assay development labs needing high throughput and high performance

Summary of Modalities¹

	VICTOR Nivo	EnVision
UV-Vis Absorbance	Filter or spectrometer	Filter and/or quad monochromators
Fluorescence Intensity	Filter	Filter and/or quad monochromators
Luminescence	• ²	• ²
Ultrasensitive Luminescence		•
TRF and TR-FRET	Lamp based ³	Lamp or laser based ³
Fluorescence Polarization	•	•
Alpha (Laser-Based)	Alpha Standard	Alpha Standard or Alpha HTS
AlphaPlex (Laser-Based)		Alpha Standard
Dual PMT Detector		•

1 - All instruments are configurable. 2 - Capable of BRET/BRET2 assays. 3 - HTRF® Certified.

For more information, visit PerkinElmer.com/category/microplate-readers or Cisbio.com/readers