

GREATER CONFIDENCE IN YOUR RESULTS WITH MULTIPLE POWERFUL DETECTION TECHNOLOGIES



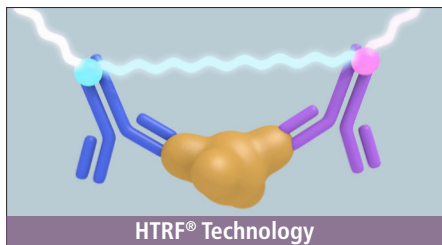
VICTOR® Nivo™ and EnVision® with HTRF® Certification and Alpha Technology



Alpha Technology

Now certified for use with HTRF® technology, PerkinElmer's VICTOR Nivo and EnVision with multiple detection technologies including Alpha empowers you to take an orthogonal approach to your research by gathering insights from different perspectives to give you greater confidence in your results.

With HTRF® technology, you will benefit from its ease of use and no wash assays. HTRF kits, with a portfolio that is constantly growing, can enhance your research capabilities on major indications, including but not limited to immuno-oncology, neuroscience, and cardiology.



HTRF® Technology

As for Alpha Technology, it is a versatile, bead-based platform that gives you the power to assay even the most complex samples – all in one well, with no wash steps. Alpha Technology has a solution for nearly every research application.

PerkinElmer has more than 30 years of experience in instrument development so you can expect excellent performance and reliability from our plate readers. With such robust detection capabilities, our microplate readers can help research, screening, and assay development labs achieve scientific discoveries faster and more efficient.

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

PerkinElmer, Inc.
940 Winter Street
Waltham, MA 02451 USA
P: (800) 762-4000 or
(+1) 203-925-4602
www.perkinelmer.com



For a complete listing of our global offices, visit www.perkinelmer.com/ContactUs

Copyright ©2019, PerkinElmer, Inc. All rights reserved. PerkinElmer® is a registered trademark of PerkinElmer, Inc. All other trademarks are the property of their respective owners.