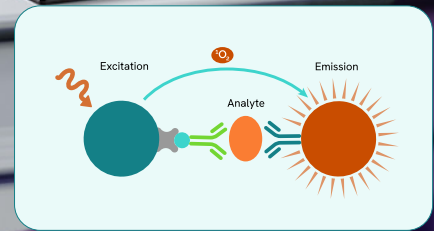


Research
at the speed
of now.



revvity

Multimode detection solutions

Detection technologies for today – and tomorrow

All our microplate readers are equipped with multiple detection modes so you can meet various assay requirements. We also offer optimized reagents and microplates that can expedite your workflows, reduce costs, and provide reproducible results.

In addition to our portfolio of products and services, our deep-seated scientific expertise provides you with a complete research solution.



VICTOR Nivo™ multimode plate reader

For low-throughput assays

- Compact, lightweight, and versatile
- For your everyday biochemical and cell-based assays
- Available with all major detection modes
- Intuitive browser-based software for remote access



EnVision™ XCite multimode plate reader

For medium to high throughput assays

- Hybrid system combining filters for sensitivity and monochromators for assay flexibility
- Freely configurable and upgradable



EnVision Nexus™ multimode plate reader

For demanding high-throughput assays

- Provides ultimate speed and sensitivity
- Freely configurable and upgradable
- Fully automatable for walkaway convenience
- Options for both high sensitivity and speed for Alpha, Luminescence, and TRF

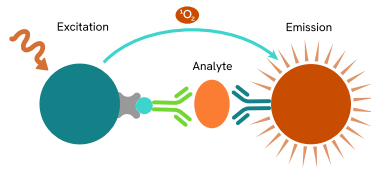
Choose the right plate reader for your research

Plate Reader	Absorbance	Fluorescence Intensity	Luminescence	Ultrasensitive Luminescence	TRF and TR-FRET	Fluorescence Polarization	Alpha (Laser-Based)	AlphaPlex (Laser-Based)	Dual PMT Detector
VICTOR Nivo	Filter or spectrometer	Filter	Including BRET/BRET2 assays		Lamp based, HTRF® certified	●	Alpha Standard		
EnVision XCite	Filter and quad monochromators	Filters and quad monochromators	Including BRET/BRET2 assays	●	Lamp based, HTRF® certified	●	Standard or HTS Alpha	●	
EnVision Nexus	Filter	Filter	Including BRET/BRET2 assays	●	Lamp or laser based, HTRF® certified	●	Enhanced Alpha and/or HTS Alpha	●	●

Choose the right technology for your research

Alpha

- Versatile, no-wash, intuitive platform
- Compatible with complex samples
- Capable of measuring virtually any target



Brands:

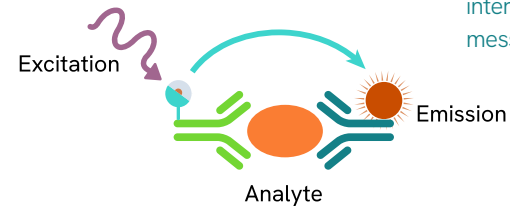
AlphaLISA®
AlphaLISA SureFire® Ultra™
Alpha Screen®
AlphaPlex®

Applications:

Proteins and phosphoproteins, protein: protein interactions, receptor-ligand interactions, DNA interactions, post-translational protein modifications

TR-FRET

- Easy-to-use, no-wash, and ultrasensitive
- Highly reproducible assay platform
- Used for detection, quantitation, and screening



Brands:

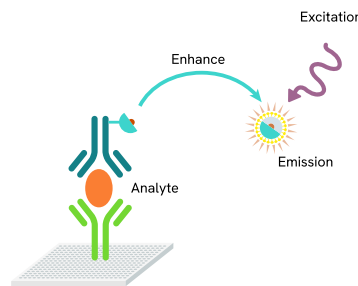
HTRF®
LANCE®

Applications:

Proteins and phosphoproteins, protein: protein interactions, receptor-ligand interactions, second messenger levels

TRF

- Long signal stability
- Multiplexing capability
- Broad sample compatibility



Brands:

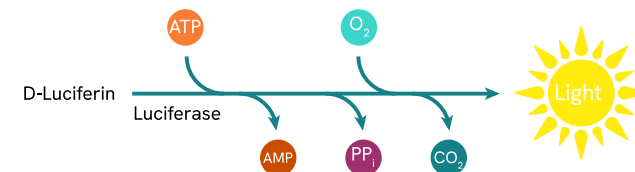
DELFI A®

Applications:

Proteins and phosphoproteins, protein: protein interactions, cytotoxicity, cell proliferation

Luminescence

- Significant for drug development
- For research in transcriptional gene expression, virus life cycles, and cell viability



Brands:

britelite™ plus, steadylite™ plus, neolite™, sensilite™, twinlite™, ATPlite™

Applications:

Cytotoxicity and cell proliferation, reporter gene, kinase activity



Discover better results with the right microplate.

Engineered to deliver the highest quality data, our microplates are optimized for our readers and reagents to provide the results you need - fast. Choosing the right microplate can help you reduce costs and meet project deadlines.

Explore our variety of microplates:

- Storage
- Cell culture
- Cell imaging
- Next generation sequencing
- Research

www.revivity.com/platereaders



revvity