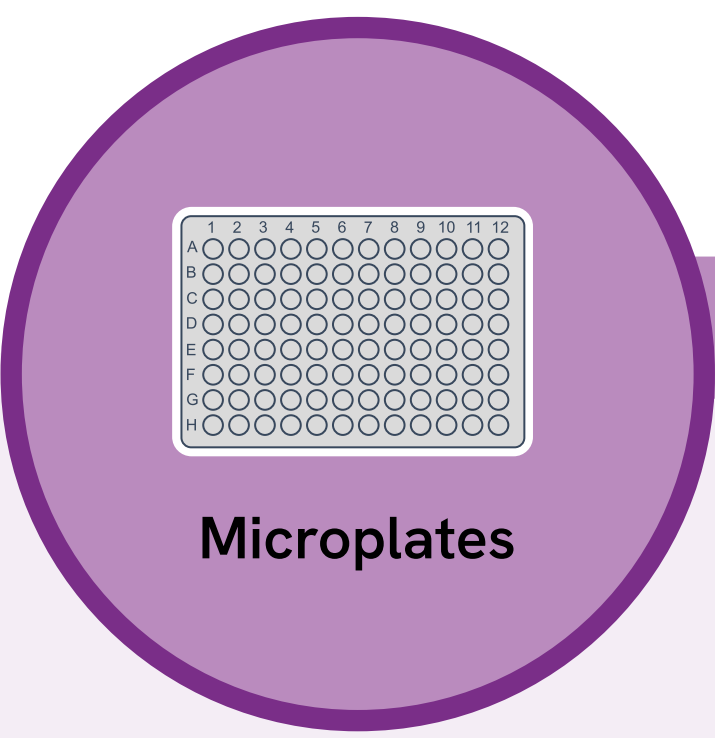
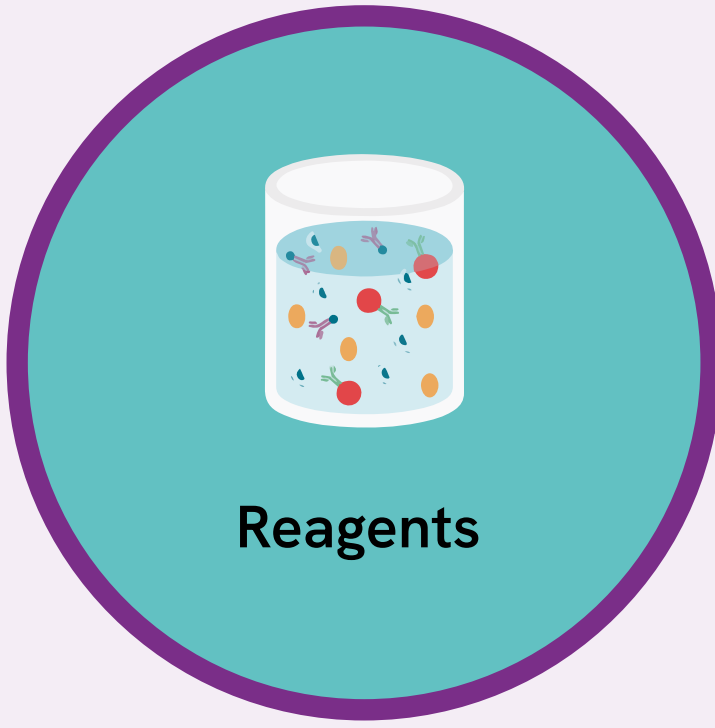


Accelerating drug discovery with walkaway automation.

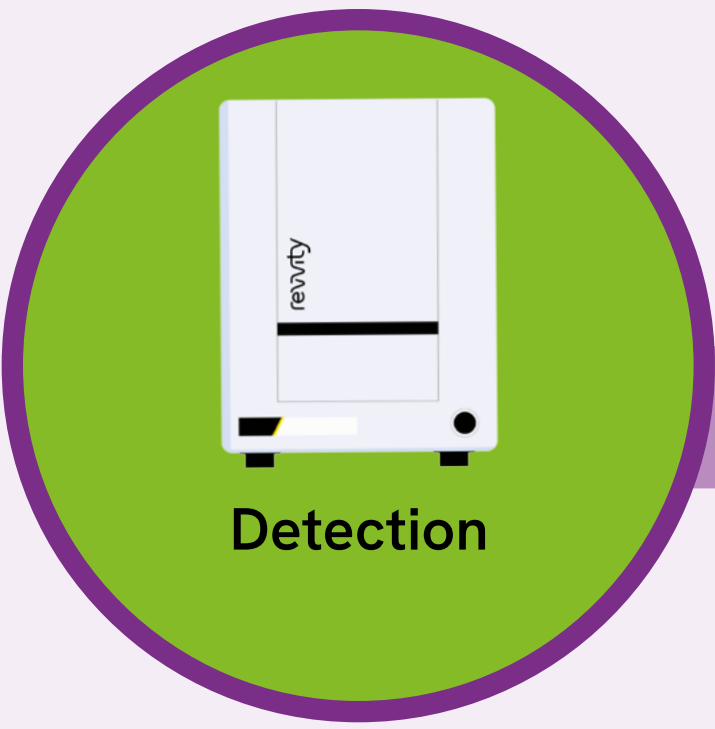
Optimize your workflow to set your lab up for success. Walkaway automation enables you to focus more on your research and less on plate management.



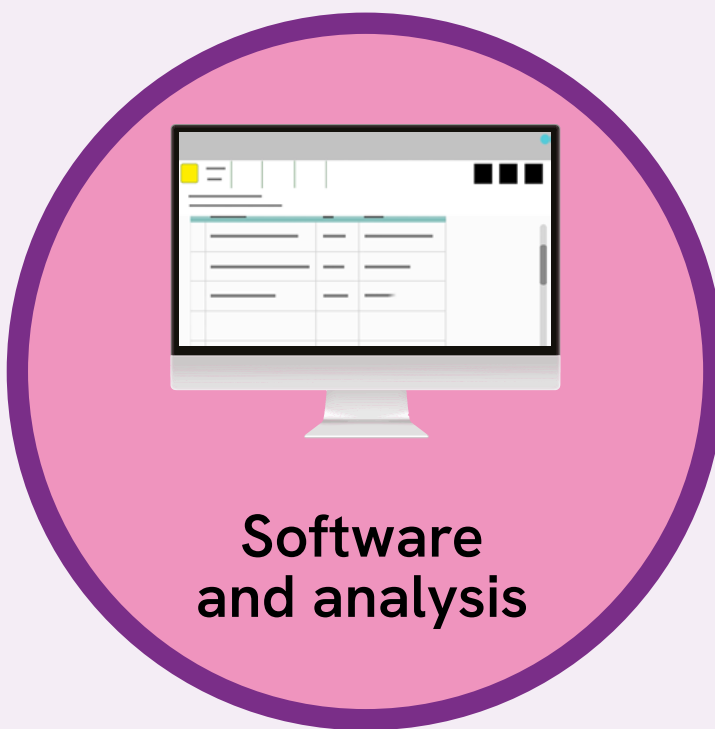
Microplates



Reagents



Detection



Software and analysis



QA/QC

Better plates for better results



Research

For fluorescence, luminescence, absorbance, Alpha, and HTRF assays



Cell culture

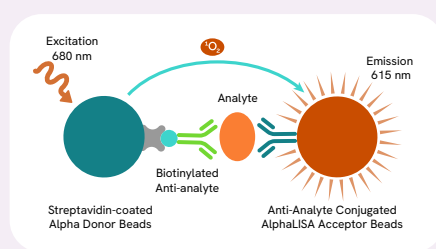
Improves cell growth consistency and reliability for general and 3D cell cultures



Next generation sequencing

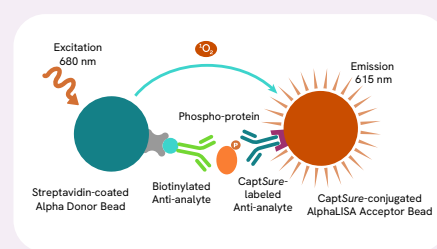
For next generation sequencing (NGS) used in sample preparation

Expand your research



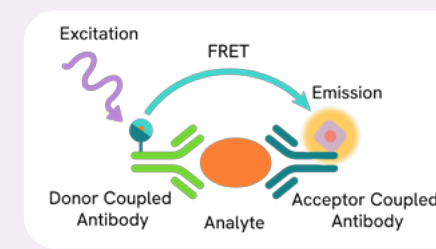
AlphaLISA™

Perfect for high sensitivity and compatibility with many complex samples



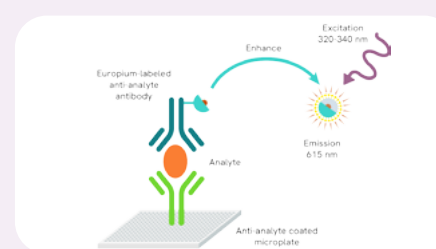
Alpha SureFire® Ultra™

Provides flexibility in using cultured cell lines, stem cells, and primary cells



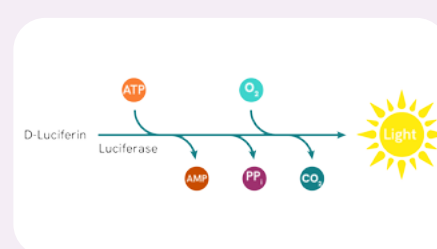
HTRF®

Delivers faster time-to-results and better reproduction than ELISA



DELFI A®

Detects and quantifies proteins in both simple and complex samples



Luminescence

Offers wide dynamic range and low interference with greater sensitivity than fluorescence technologies

Future-proof your lab



Absorbance

For ELISA, protein and nucleic acid quantification, and enzyme activity



Fluorescence Intensity

For DNA or protein quantitation, reporter-gene expression, and protein binding



Luminescence

Offers wide dynamic range, high sensitivity, and low interference



Time-Resolved Fluorescence

Improves sensitivity and dynamic range of immunoassays



Fluorescence Polarization

For analysis of molecular interactions to study binding or dissociation properties



Alpha

Assesses complex samples all-in-one well, no-wash steps, and quick readouts

Increase your productivity with the VICTOR Nivo™ and Stacker

Our VICTOR Nivo multimode plate reader is perfect for everyday biochemical and cell-based assays. When combined with the stacker, it delivers walkaway automation so you can focus more on your research and less on plate management.



[Watch video](#)

Research & discovery

Development & manufacturing

21 CFR PART 11 COMPLIANCE



Enhanced Security Software
Provides easy-to-use tools for regulated labs to simplify the journey to compliance

We perform validation and qualification services:

- Installation qualification (IQ)
- Software qualification
- Operation qualification (OQ)
- Computer system validation (CSV)
- Assay validation

For more information, please visit www.revvy.com

revvity

Revvity, Inc.
940 Winter Street
Waltham, MA 02451 USA
www.revvy.com

For a complete listing of our global offices, visit www.revvy.com
Copyright ©2024, Revvity. All rights reserved.

1394066 (1130911)