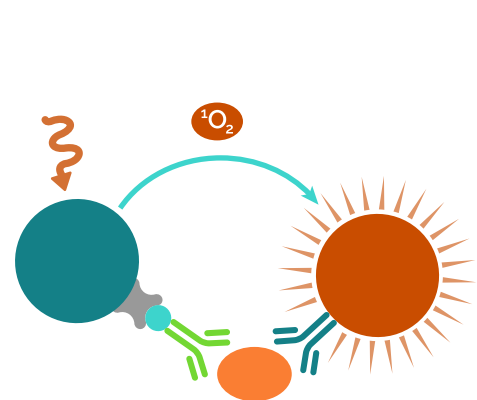
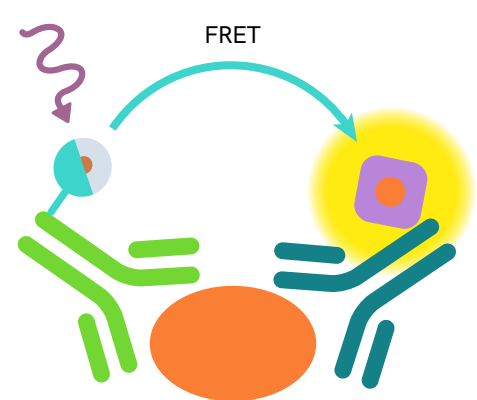


Simplify your immunoassays with a single-source solution.

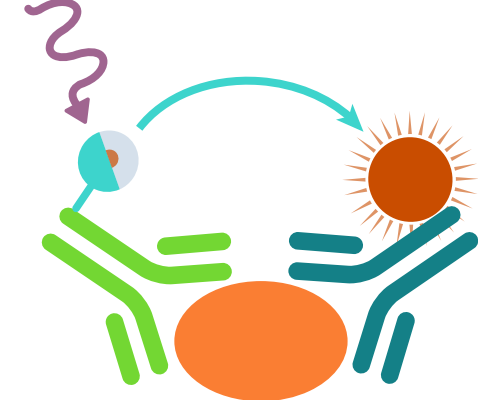
No-wash technologies



Alpha
Versatile bead-based, no-wash, luminescent technology.

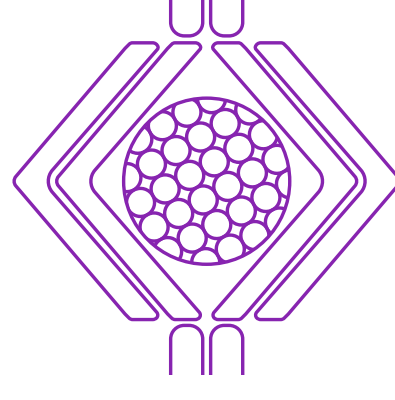


HTRF™
No-wash technology, gold reference for time-resolved FRET.

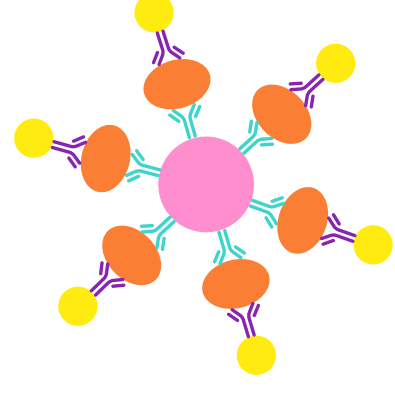


LANCE™
No-wash proximity assay technology that combines the benefits of TR with FRET.

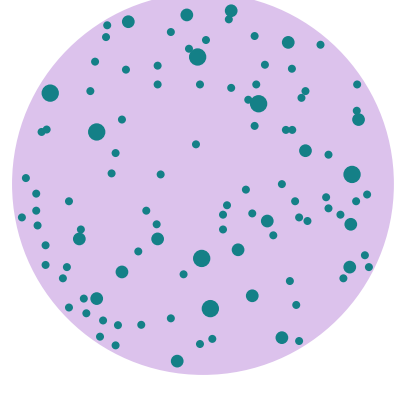
Wash-based technologies



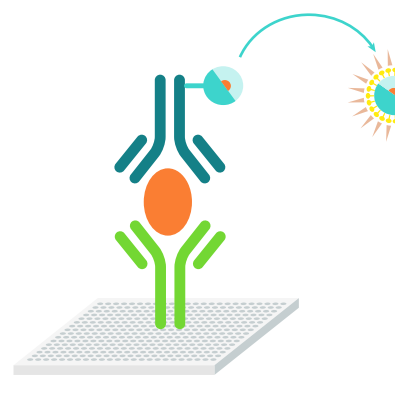
ELISA
Plate-based, colorimetric immunoassay.



LEGENDplex™
Fluorescence bead-based assay for simultaneous quantification of up to 14 analytes.



ELISpot
Cell-based assay for single-cell level analyte detection.



DELFI™
Stable, time-resolved fluorescent signal to improve assay dynamic range.

*For wash-based assays, our automation platform requires integration with a third-party washer.



Omni Bead Ruptor Elite™
bead mill homogenizer

Sample prep

Leverages the power of mechanical shearing to maximize tissue homogenization and cell lysis. Supports multiple tube sizes and bead lysis kits, with processing speeds up to 8 m/s.



No-wash and wash-based assay kits*

Immunoassays

Vast selection of immunoassay technology platforms, pre-configured kits, and customizable reagents, providing the flexibility to meet the demands of your specific research.



Fontus™ workstations

Liquid handling

Up to five configurations to accommodate workflow and space requirements. Both washable and disposable tips can be used, without user intervention, providing flexibility and time savings.

Assay microplates

Microplates

Our microplates are engineered to provide high signal integrity and well-to-well consistency for your research applications, and are available in both standard and custom-configured formats.

explorer™ G3 integrated workstations

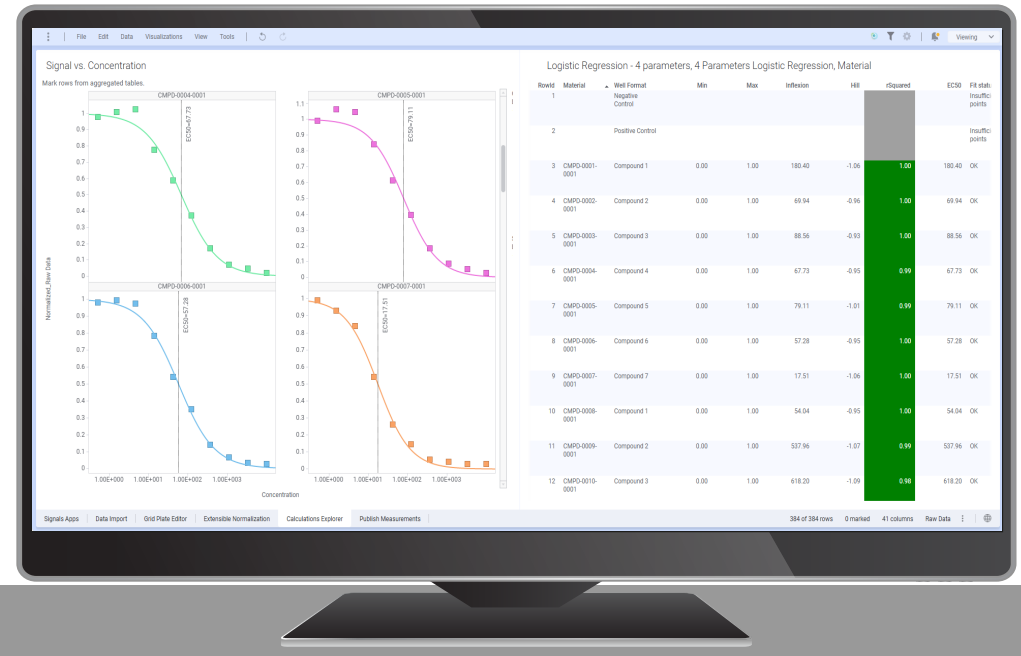
Automation

Custom integrated workstations, providing innovative automation solutions to simplify lab workflow.

EnVision Nexus™ multimode plate reader

Read-out

Configurable multimode plate readers equipped with the latest technologies for labs today.



Signals One™

Software

Software solutions that leverage the latest technologies to help simplify science and accelerate innovation.

For more information, visit: www.revvy.com/immunoassays-solutions

[Learn more](#)