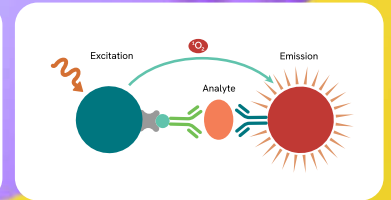
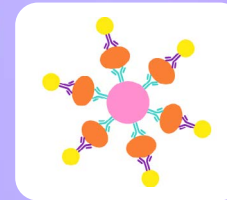


Simplify your  
immunoassays  
with an  
integrated  
resource.

revvity

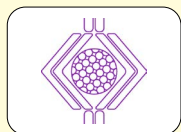


Immunoassay solutions

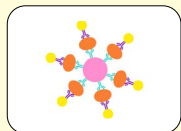
# Revolutionize your immunoassays with an integrated resource.

Advance your research with Revvity's comprehensive portfolio of immunoassays, including widely-used ELISA formats, innovative AlphaLISA™ and HTRF™ no-wash assays, and BioLegend LEGENDplex™ flow-based assays. Our diverse immunoassay technologies are designed to deliver reliable data and work effectively with our broader product portfolio to enhance the efficiency of your research workflows. Focus more of your time on your research objectives.

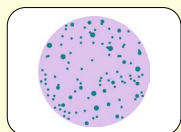
## Wash-based immunoassay technologies



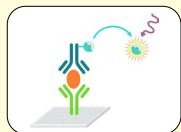
**ELISA** - Plate-based, colorimetric immunoassay



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

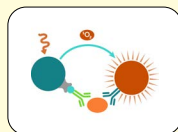


**ELISpot** - High-sensitivity, cell-based assay for single-cell level analyte detection

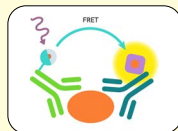


**DELFIA™** - Stable, time-resolved fluorescent signal to improve assay dynamic range

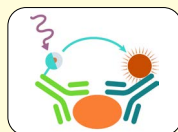
## No-wash immunoassay technologies



**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

## Determine the parameter most critical to your research to find the optimal immunoassay technology.

Ready-to-use assays to quantify biomarkers, cytokines, and other secreted proteins.

	ELISA	LEGENDplex <sup>1</sup>	HTRF	Alpha
Absorbance, luminescence, colorimetric	•			
TRF, TR-FRET, advanced fluorescence			•	
Advanced luminescence				•
Fluorescence intensity, standard fluorescence		•		
No automation	•	•	•	•
Automation			•	•
Wash steps	•	•		
No-wash steps			•	•
Testing < 1,000 data points per week	•		•	•
Testing > 1,000 data points per week		•	•	•
5 to 20µL sample volume			•	•
20 to 50µL sample volume	•	•	•	•
Assay dynamic range up to 2 logs (<100)	•		•	•
Assay dynamic range more than 2 logs (>100 fold)		•	•	•

Ready-to-use assays to measure second messenger detection, protein-protein interaction, cell signaling, and receptor ligand binding.

	HTRF	Alpha
TRF, TR-FRET, advanced fluorescence <sup>2</sup>	•	
Advanced luminescence		•
No automation	•	•
Automation	•	•
Wash steps		
No-wash steps	•	•
Testing < 1,000 data points per week	•	•
Testing > 1,000 data points per week	•	•
5 to 20µL sample volume	•	•
20 to 50µL sample volume	•	•
Assay dynamic range up to 2 logs (<100) <sup>3</sup>	•	•
Assay dynamic range more than 2 logs (>100 fold) <sup>3</sup>	•	•

1 - Requires flow cytometer capable of reading APC and PE or, in the case of instruments without a red laser, PE and PerCP.

2 - Advanced fluorescence does not apply to cell signaling applications.

3 - The assay dynamic range is only applicable to second messenger detection and receptor ligand binding.

## Immunoassay workflow

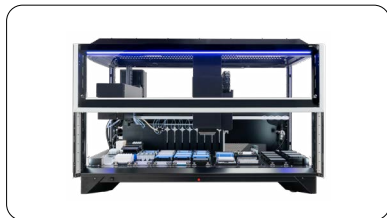
### Sample prep



#### Omni Ruptor Elite™ bead mill homogenizer and lysing kits

Up to 8 m/s processing speeds for analyte extraction from biological samples using optimized lysing bead kits.

### Liquid handling



#### Fontus™ workstations

Liquid handling powerhouse with easy-to-use software, optimized deck access, and more to simplify your workflows.

### Microplates



#### Assay 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.

### Reagents



#### Assay kits

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

### Readout



#### EnVision Nexus™ multimode plate reader

Next generation microplate reader for high-throughput screening and demanding applications.



#### Celigo™ image cytometer

Plate-based brightfield and fluorescent imaging system providing cell level insights.

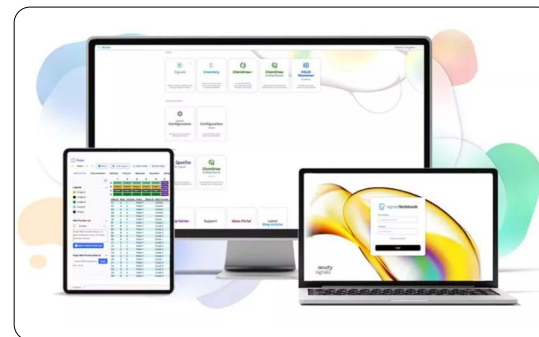
### Laboratory automation



#### explorer™ G3 workstation

From cellular testing and imaging applications to high-throughput screening and genomics-based applications, custom explorer™ G3 integrated workstations provide innovative application-focused laboratory automation solutions which help to simplify microplate handling, liquid handling, and detection.

### Software



#### Signals One™

Make your innovations a reality with a leading unified, cloud-native R&D SaaS solution that streamlines scientific complexity and promotes collaboration. Signals One is an end-to-end R&D workflow solution providing data capture & collaboration, data workflows and processing, and data-driven analytics. It sets the stage for AI/ML success by putting your data in context with AI-ready data management.

Click to learn more about our immunoassay solutions.

[www.revivity.com](http://www.revivity.com)

revvity

Copyright ©2025, Revvity, Inc. All rights reserved.  
Revvity is a trademark of Revvity, Inc. All other trademarks are the property of their respective owners.

1932554

For research use only. Not for use in diagnostic procedures.