



# Concerned about potency? Trust Revvity!

## Hit key requirement guidelines with biochemical and cell-based assays

A biological drug's stability and efficacy are critical quality attributes that need evaluation with potency estimation being the most important parameter to confirm lot-to-lot consistency and equivalence.

To assess relative potency of a biological drug product bioassays are needed to compare biological response with a control via determination of mode of action.

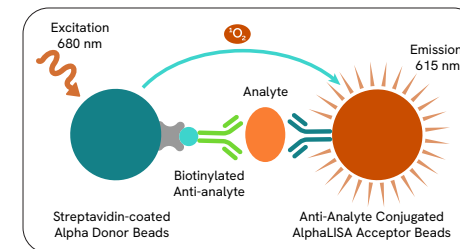
## Enable compliance with Alpha and HTRF technology

Our assay platforms are designed to provide an accurate and precise response that can evaluate the impact of drug concentrations for optimum potency ranges and help assess lot-to-lot consistency and equivalence of biotherapeutics.

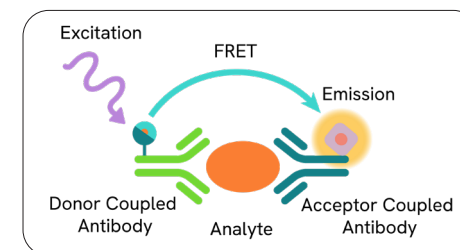
## Fit the solution to your needs with custom capabilities

With more than 30 years of experience in custom laboratory services, let's extend your lab capabilities in assay design and development to cater to your specific development and profiling requirements. Based on trust and collaboration, achieving efficiency has never been easier.

### Potency cell-based assays



| Alpha technology



| HTRF technology

### Detection



| VICTOR Nivo™ multimode plate reader



| EnVision™ Nexus™ multimode plate reader

## The platforms to detect and evaluate potency levels

### HTRF™ and AlphaLISA™ Kits for reliable potency assessment

Our no-wash HTRF and AlphaLISA cytokine and other biomarker detection assays aid potency assessments with enhanced precision, sensitivity, and efficacy.

- User-friendly with no-wash steps
- Easily automatable
- Fewer touchpoints for reproducible handling
- Assays display exceptional dilutional linearity
- Greater sensitivity and broader dynamic range
- Compatible with broad range of buffers
- Reliable antigen spike recovery
- Robust reproducibility
- Host cell protein detection kits also available

[View HTRF assay](#)

[View AlphaLISA assay](#)

### Customization services

Gain productivity and expert advice for solutions to meet your research challenges. Our commitment to you includes accurate timelines, minimized disruption of your workflow, and complete delivery of our agreed service to you.

- Assays and reagents
- Microplate services
- Custom antibody conjugation and labeling

[View services](#)

## Other solutions for your adjacent workflow

### Cell counting and image cytometry instruments

Offering a leading portfolio of automated cell counters, image cytometry systems, and cell counting reagents and consumables to perform a range of cell based functional assays. Providing viability and cell density testing with GMP and 21 CFR Part 11 compliant solutions.

[View platforms](#)



### Microfluidic protein characterization

LabChip GXII Touch is an excellent tool for high throughput titer and quality assessment of antibodies and analysis of Critical Quality Attributes such as size, purity, titer, charge variance, and glycosylation in one platform.

[View platforms](#)



### Microplate readers

Providing a portfolio of multimode plate readers equipped with the latest detection technologies for biochemical and cell-based assays. Software capabilities to support 21 CFR Part 11 compliance.

[View platforms](#)



### Automated liquid handling workstations

With workstations designed for a diverse variety of applications, along with the ability to customize, our innovative automated liquid handling solutions help minimize errors, reduce hands-on time, and increase throughput and reproducibility.

[View platforms](#)



revvity