



# Living biologics.

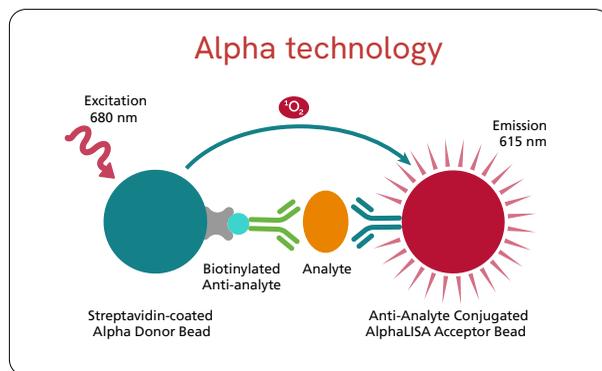
Biotherapeutics research is by far the fastest-growing sector in the pharmaceutical industry and it's no surprise that scientists around the world are constantly seeking ever more efficient tools.

Developing these therapeutic proteins requires extensive analytical characterization at every step, from lead selection to lead optimization and characterization, including functional activity testing, right through to bioprocessing qualification.

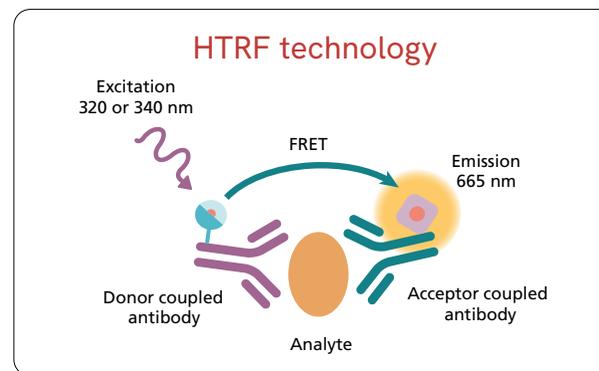
To meet this need, Revvity offers a comprehensive line of tools for biologics screening, mechanism-of-action studies, and biomanufacturing, using the robustness and convenience of a no-wash technology.

## Advance your biologics research with Revvity solutions

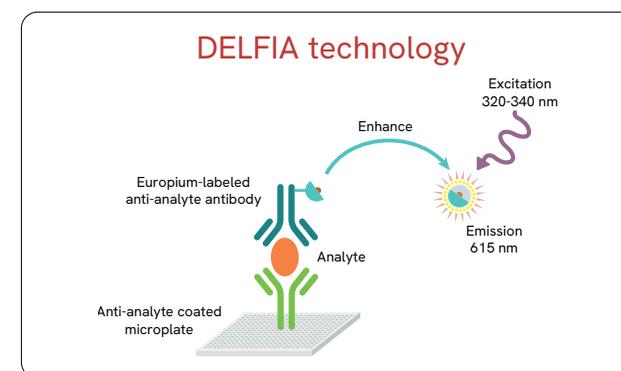
From discovery through manufacturing, harnessing living cells to create safe and effective therapies is complex and expensive. Revvity helps scientists develop and streamline their entire biologics workflow so they can overcome these challenges and bring consistent, high-quality, biological medicines to market faster.



Alpha is a versatile, bead-based platform that enables you to assay the most complex samples in one well and with no wash steps.



HTRF is a fast, sensitive, homogeneous, and ready to use assay platform with no wash steps.



DELFI has a similar assay principle and workflow to that of a traditional ELISA, but with the added benefits of a stable, time-resolved fluorescent signal improving assay dynamic range.



VICTOR™ Nivo™ multimode plate reader



EnSight™ multimode plate reader



EnVision™ Nexus™ multimode plate reader

## Ig detection

Ready-to-use kits to quantify IgGs and biologics fragments, offering a homogeneous, easily miniaturized alternative to conventional ELISAs.

	AlphaLISA™	HTRF™
Human IgG	AL205C	6FHIGPEG
Human IgG1	AL307C	
Human IgG4	AL310C	
Pan Mouse IgG	AL584C	
Mouse IgG	AL529C	
Mouse IgG1	AL522C	
Mouse IgG2a	AL523C	
Mouse IgG2b	AL524C	
Mouse IgG3	AL525C	
Human Fc	AL3129C	62HFCPEG
Mouse Fc		6FMIGPEG
Rabbit Fc		6FRIGPEG
Kappa MAb	AL3023C	64KAPPEG
Lambda MAb		64LAMPEG
Human IgM	AL263C	
Mouse IgM	AL527C	
Mouse IgE	AL526C	
Hybridoma (Mouse) Binding - GST Antigen		64HYGSTPEG
Hybridoma (Mouse) Binding - HIS Antigen		64HYHISPEG

## Cytokine receptor/ligand binding assays

Ready-to-use kits to discover blocking biologics.

	AlphaLISA™	HTRF™
Human TNFa/TNFR1 Binding Kit	AL3120C	64BDTNFPEG
Human TNFa/TNFR2 Binding Kit	AL3121C	
Human IL6/IL6R Binding Kit		64BDIL6PEG
Human IL12/IL12Rβ1 Binding Kit		64BDIL12PEG
Human IL17A/IL17RA Binding Kit		64BDIL17PEG
Human IL23/IL12Rβ1 Binding Kit		64BDPIL12PEG
Human IL23/IL23R Binding Kit		64BDPIL23PEG
Human 4-1BB/4-1BBL Binding Kit	AL3113C	

## ADCC

DELFIATM EuTDA Cytotoxicity Reagents facilitate the sensitive detection of ADCC using the same workflow as the Chromium-51 release assay but in a convenient-to-use, non-radioactive format.

	DELFIATM
DELFIATM EuTDA Cytotoxicity Reagents	AD0116

## Other binding assays

	AlphaLISA™	HTRF™
Human EGF/EGFR Binding Kit	AL366C	
Human PCSK9/LDLR Binding Kit	AL3132C	
SARS-CoV-2 S1/ACE2 Binding kit	AL3146C	64BDS1PEG

## Fc gamma receptor (FcγR) and neonatal Fc (FcRn) binding assays

Tag-lite™-based cellular assays or biochemical binding assays for measuring FcγR-specific antibody binding. These assays are ideal for predicting antibodies with strong ADCC activity and longer circulating half-lives. They can also be used to assess the degree of antibody afucosylation.

	AlphaLISA™	HTRF™
CD16a (FcγRIIIa, F158)	AL347C	6FR3APAG
CD16a (FcγRIIIa, V158)	AL348C	62C16PAG
CD16b (FcγRIIIb)	AL3094C	
CD32a (FcγRIIa, H131/167)	AL3086C	6FRAHPAG
CD32a (FcγRIIa, R131/167)	AL3087C	6FRARPAG
CD32b (FcγRIIb)	AL3080C	6FR2BPAG
CD64 (FcγRI)	AL3081C	6FC64PAG
FcRn	AL3095C	64FCRNPET

## ADCC and CDC related readouts

	AlphaLISA™	HTRF™
Human Granzyme B	AL3170C	63ADK095PEH
Human Perforin-1	AL3170C	63ADK100PEG
Human C1Q binding		64C1QPET
Serum Albumin (human) - HSA	AL294C	
Serum Albumin (human) Biotin-Free	AL363C	
Serum Albumin (mouse)	AL508C	
Anti-Peg IgM	AL313C	
Host Cell Residual DNA	AL331C	

## Bioprocess

	AlphaLISA™	HTRF™
Anti-Peg IgM	AL313C	
CHO HCP <b>NEW</b>	AL3176C	64CHOPEG
Host Cell Residual DNA	AL331C	
Serum Albumin (Human) Biotin-Free	AL363C	
Serum Albumin (Mouse)	AL508C	
Serum Albumin (Human) - HSA	AL294C	

## Immune checkpoint binding assays

Ready-to-use kits to screen for interaction inhibitors.

	AlphaLISA™	HTRF™
Human CD47 / SIRP alpha Binding Kit		64SIRPEEG
Human CTLA4/B7-1(CD80) Binding Kit	AL3046C	64CTLA80PEG
Human CTLA4/B7-2 (CD86) Binding Kit	AL3047C	64CTLA86PEG
Human LAG3/MHC II Binding Kit		64LAG3PEG
Human HLA-DRA/LAG3 Binding Kit	AL3066C	
Human PD1/PD-L1 Binding Kit	AL356C	64PD1PEG
Cynomolgus PD1/PD-L1 Binding Kit	AL583C	
Mouse PD1/PD-L1 Binding Kit	AL580C	
Human PD1/PD-L2 Binding Kit	AL3128C	
Human CD28/CD86 Binding Kit	AL3131C	
Human PDGF-BB/PDGFR $\alpha$ Binding kit	AL3064C	
Human PDGF-BB/PDGFR $\beta$ Binding kit	AL3065C	

References indicated for 500 tests, other kit sizes are also available.

## Toolbox reagents

With the generation of complex molecules such as biotherapeutics drugs comes the need to better characterize their potential as drug candidates. One of the criteria for evaluating their therapeutic potential is the binding efficiency of an antigen present on the cell surface, or more generally, the protein-protein interactions.

Our reagents can help screen medium to large biologics libraries produced by hybridoma or in periplasmic crude extract or run affinity maturation experiments.

[View reagents here](#)

## Custom solutions for drug discovery

Revvity is proud to be able to share our more than 30 years of experience in custom laboratory services, and to accompany scientists each year in their drug discovery research in a personalized way based on trust, exchange and efficiency.

We offer a range of custom assay development, profiling and labeling, to custom radiosynthesis.

[Learn more here](#)

For more information, visit [www.revvity.com/category/protein-analysis](http://www.revvity.com/category/protein-analysis)

revvity