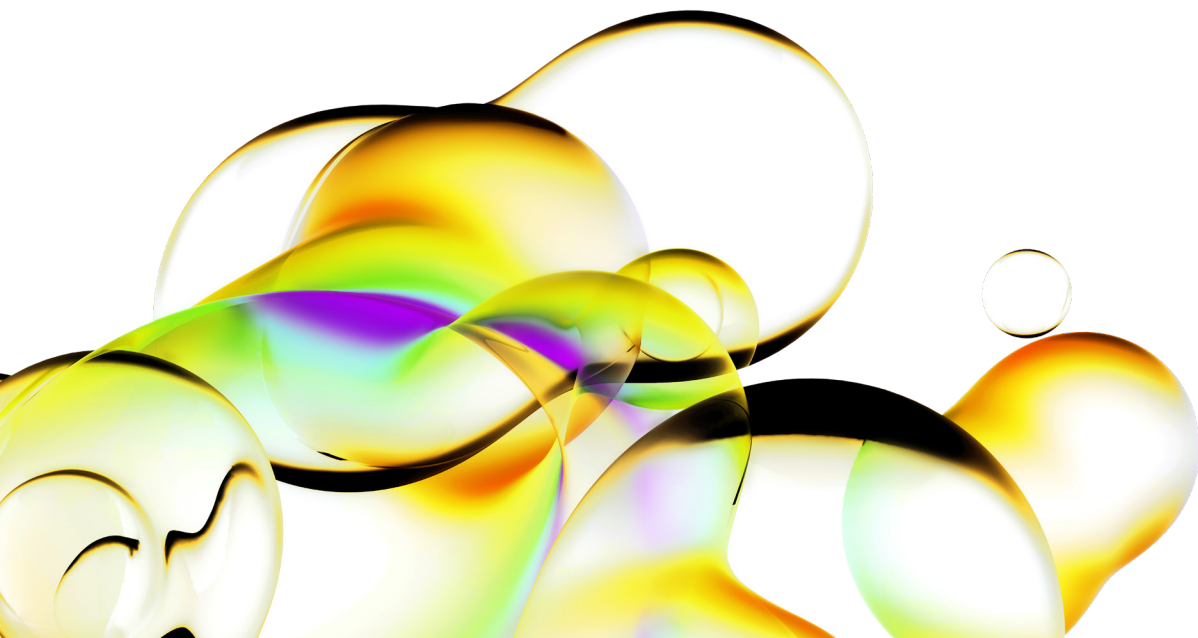




Extend your knowledge on protein-protein interaction.



HTRF® protein-protein interaction reagents

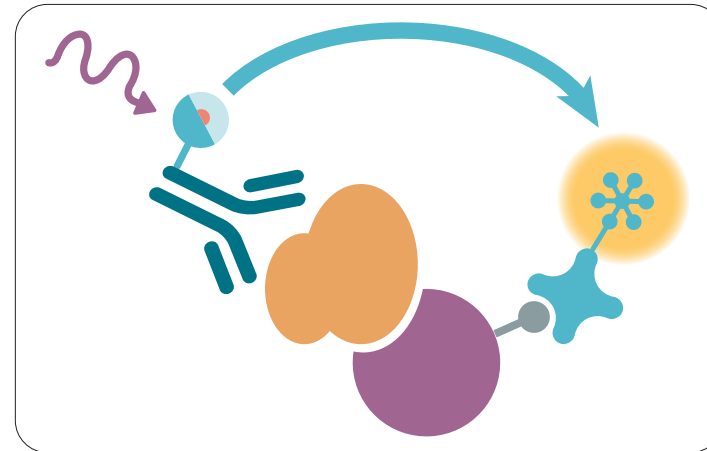
With their central role in biological processes, protein-protein interactions (PPIs) are compelling targets for drug discovery. But only the right tools will take you where you want to go.

Revvity's HTRF® PPI enables you to build robust assays that are precisely adapted to your own biochemical assay format. Your assays will be ready for screening with minimum development time.

Satisfying researchers for over a decade, HTRF PPI reagents help you investigate and screen many types of targets, from proteases and immunoassays to protein modification assessment and nucleic acid-related assays.

Benefit from unlimited flexibility for your assay development

- Tools to support any tag, any configuration
- Add-and-read format for all throughput needs
- Validated on numerous target configurations
- Labeling kits and custom labeling services available for coupling untagged proteins directly



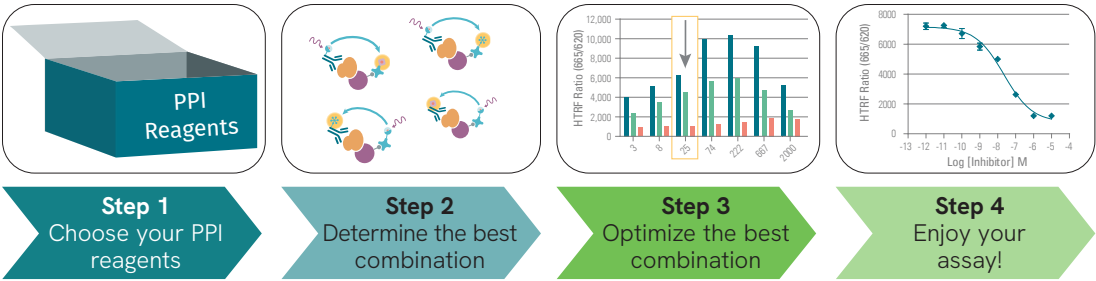
Studying protein-protein interactions

With support for multiple anti-tag probes, streptavidins, Strep-Tactin® and anti-species, HTRF PPI reagents adapt effectively to your own proteins and target materials.

How it works

Assay tools that let you explore all your best ideas

HTRF PPI reagents include Europium and Terbium cryptates, as well as XL665 and d2 labeled anti-tag antibodies, anti-species antibodies, and streptavidins. For any assay type, tag, or format, these reagents enable you to move forward with screening in record time by taking advantage of reliable and easy-to-use tools.



Ordering Information

Surrogates	Cryptate Donors					Acceptors					
	Eu3+		Tb3+			XL665			d2		
	5,000 Tests	20,000 Tests	1,000 Tests	5,000 Tests	20,000 Tests	1,000 Tests	5,000 Tests	20,000 Tests	1,000 Tests	5,000 Tests	20,000 Tests
Anti-GST	61GSTKLA	61GSTKLB	61GSTTLF	61GSTTLA	61GSTTLB	61GSTXLF	61GSTXLA	61GSTXLB	61GSTDLF	61GSTDLA	61GSTDLB
Anti-6HIS	61HISKLA	61HISKLB	61HISTLF	61HISTLA	61HISTLB	61HISXLF	61HISXLA	61HISXLB	61HISDLF	61HISDLA	61HISDLB
Anti-6HIS Gold	61HI2KLA	61HI2KLB	61HI2TLF	61HI2TLA	61HI2TLB	-	-	-	-	-	-
Anti-c-myc	61MYCKLA	61MYCKLB	61MYCTAF	61MYCTAA	61MYCTAB	61MYCXLF	61MYCXLA	61MYCXLB	61MYCDAF	61MYCDAA	61MYCDAB
Anti-FLAG®	61FG2KLA	61FG2KLB	61FG2TLF	61FG2TLA	61FG2TLB	61FG2XLF	61FG2XLA	61FG2XLB	61FG2DLF	61FG2DLA	61FG2DLB
Anti-HA	610HAKLA	610HAKLB	610HATAF	610HATAA	610HATAB	610HAXLF	610HAXLA	610HAXLB	610HADAF	610HADAA	610HADAB
Anti-MBP	61MBPKAA	61MBPKAB	-	61MBPTAA	61MBPTAB	-	-	-	-	61MBPDAA	61MBPDAB
Anti-DNP	61DNPKLA	61DNPKLB	-	-	-	-	61DNPXLA	61DNPXLB	-	-	-
Anti-mouse Ig	61PAMKLA	61PAMKLB	61PAMTAF	61PAMTAA	61PAMTAB	61PAMXLF	61PAMXLA	61PAMXLB	61PAMDAF	61PAMDAA	61PAMDAB
Anti-rabbit Ig	61PARKLA	61PARKLB	61PARTAF	61PARTAA	61PARTAB	61PARXLF	61PARXLA	61PARXLB	61PARDAF	61PARDAA	61PARDAB
Anti-human Ig	61HFCKLA	61HFCKLB	61HFCTAF	61HFCTAA	61HFCTAB	61HFCXLF	61HFCXLA	61HFCXLB	61HFCDAF	61HFCDAA	61PARDAB
Protein A	61PRAKLA	61PRAKLB	-	-	-	-	61PRAXLA	61PRAXLB	-	-	-
Streptavidin	610SAKLA	610SAKLB	610SATLF	610SATLA	610SATLB	610SAXLF	610SAXLA	610SAXLB	610SADLF	610SADLA	610SADLB
Streptavidin Xlent!	-	-	-	-	-	-	611SAXLA	611SAXLB	-	-	-
Strep-Tactin®	610STKLA	610STKLB	-	610STTLA	610STTLB	-	610STXLA	610STXLB	-	610STDLA	610STDLB

PPI detection buffers	
PPI - Terbium detection buffer - 200 mL	61DB10RDF
PPI - Europium detection buffer - 200 mL	61DB9RDF

