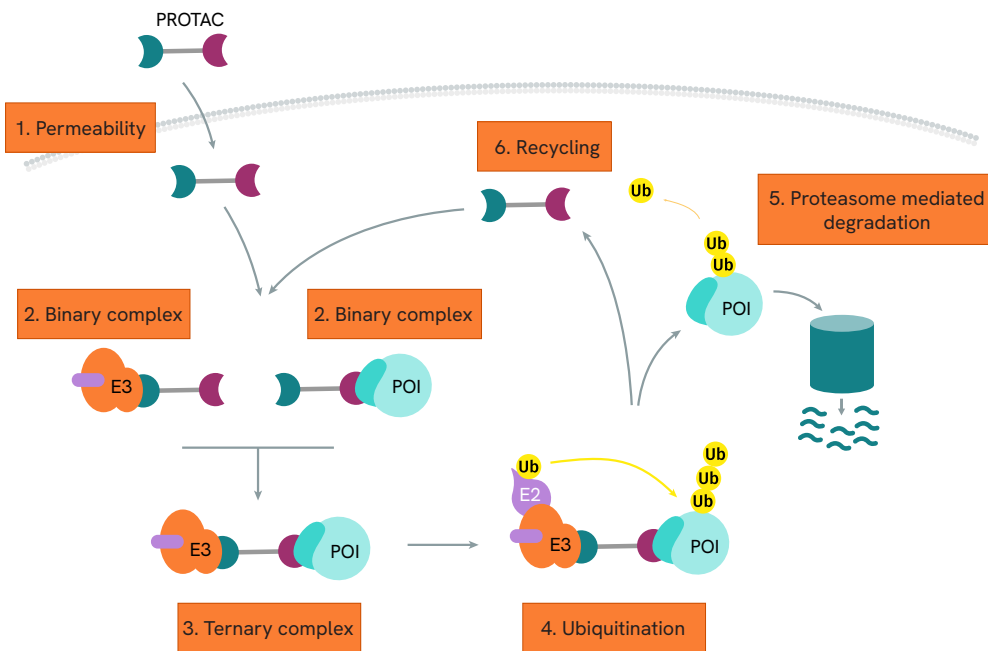




# Reagent solutions for targeted protein degradation.



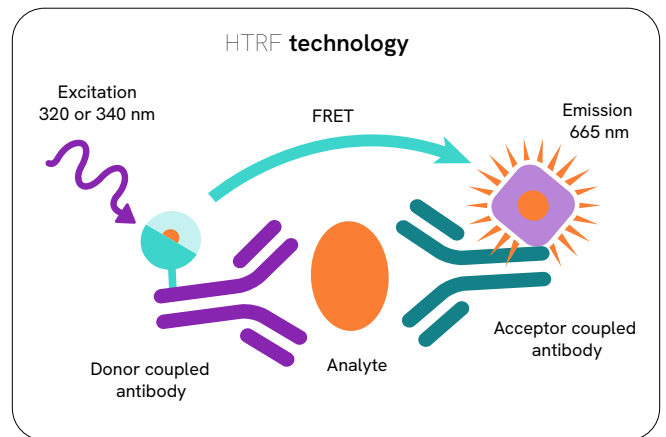
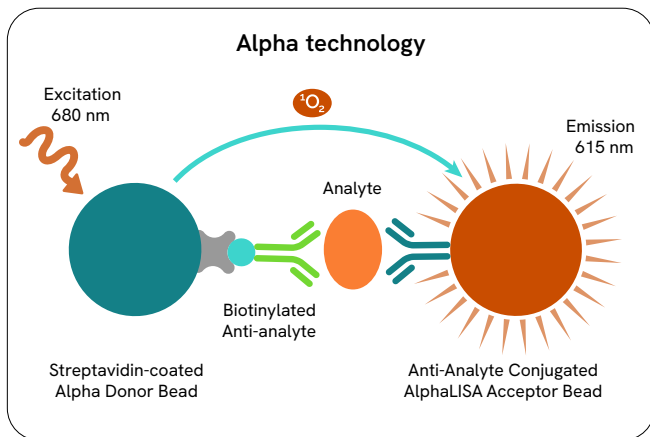
Proteolysis Targeting Chimaeras (PROTACs) is a game changing mechanism for treating the untreatable in the area of drug discovery.

A PROTAC is a small molecule with a binding region on each side. One recognizes an E3 ubiquitin ligase and the other a protein of interest. Both are joined by a linker. When the linker recognizes both E3 ligase and POI, it forms a ternary complex. The protein of interest is ubiquitinated and is then degraded by the proteasome.

Contrary to current drugs which inhibit the proteasome, this mode of action leads to the destruction of the target. Consequently, it could be more effective than current cancer treatments.

## Advance your PROTAC<sup>®</sup> research and benefit from Revvity's broad offering of reagent technologies:

- A large biochemical assay portfolio, including GPCR research, cell signaling, biomarkers, or protein-protein interaction
- Homogeneous easy-to-use immunoassay kits for cell-based detection of cytokines, biomarkers, and cell signaling proteins which provide fast and accurate results utilizing Alpha and HTRF assay technologies
- Alpha CETSA<sup>®</sup> technology to monitor binding of your compound of interest in the natural living cellular environment by bringing together the Cellular Thermal Shift Assays (CETSA<sup>®</sup>) from Pelago Bioscience with Revvity Alpha technology



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.

### Does my PROTAC bind to E3 ligase?

Streptavidin-coated Alpha Donor Bead

Anti-TAG Conjugated AlphaLISA Acceptor Bead

Anti-GST Eu3+ antibody

GST

Cereblon protein GST tagged

Thaldomide-Red

- E3 ligase binding kits
- HTRF PPI reagents
- Alpha PPI kit

### Is a ternary complex formed?

Streptavidin-coated Alpha Donor Bead

Anti-Analyte Conjugated AlphaLISA Acceptor Bead

HTRF donor and acceptor labeled anti-TAG antibodies

FRET

- HTRF PPI reagents
- Alpha PPI kit

### Does my PROTAC compound bind and degrade my target?

Ub

- Alpha CETSA<sup>®</sup> kits
- HTRF total protein kits
- HTRF total Cereblon kit
- AlphaLISA<sup>®</sup> SureFire<sup>®</sup> Ultra<sup>™</sup> total protein kit
- Alpha total Cereblon kit

## Custom labeling and custom assay development services

Don't see the assay you need listed here? Revvity has partnered with industry and academia professionals to develop custom solutions to dramatically shorten the drug development process using our powerful, accurate HTRF®, DELFIA® and AlphaLISA® technologies.

An experienced Project Manager will evaluate your project's chance of success, then work with you to create a customized assay for the target you need to advance your research.

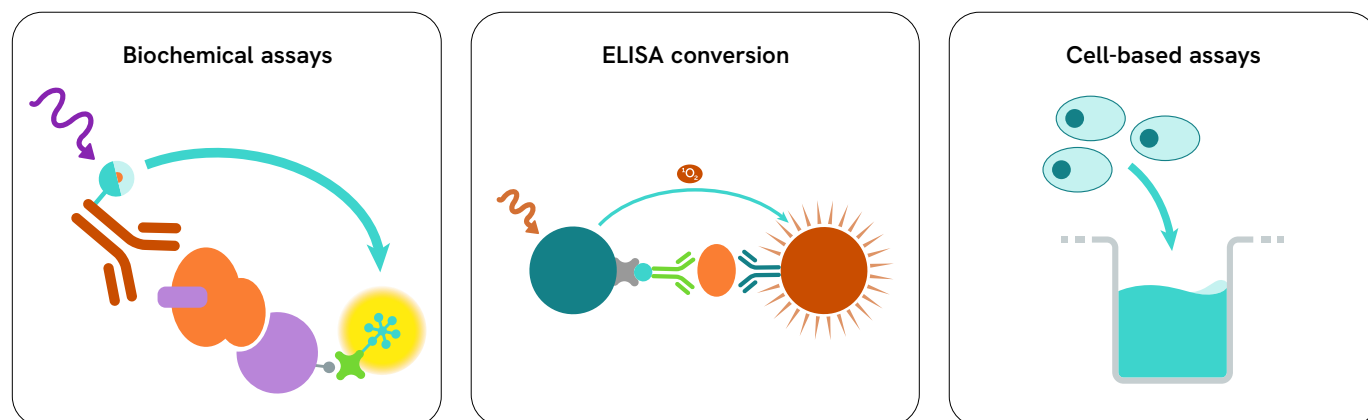


Table 1. HTRF and AlphaLISA E3 ligase binding kits.

Product	Size	Part number	
		HTRF	AlphaLISA
Cereblon Binding kit	500 tests	64BDCRBNPEG	AL3147C
VHL Binding kit	500 tests	64BDVHLPEG	
cIAP1 BIR2 Binding Kit	500 tests	64BDIAPB2PEG	AL3166C
cIAP1 BIR3 Binding Kit	500 tests	64BDIAPB3PEG	AL3167C
XIAP1 BIR2 Binding Kit	500 tests	64XIAPB2PEG	AL3172C
XIAP1 BIR3 Binding Kit	500 tests	64XIAPB3PEG	AL3173C
MDM2 Binding Kit	500 tests	64BDMDM2PEG	AL3168C
Thalidomide - Red Ligand	10 tests	64BDCRBNRED	
VHL - Red Ligand	10 tests	64BDVHLRED	
LCL161 - Red Ligand	10 tests	64BDIAPB3RED	
MI-1061 - Red Ligand	10 tests	64BDMDM2RED	
XIAP BIR3 Red Ligand	10 tests	64XIAPFLRED	
PROTAC Binding Buffer 1	200 mL	64BDE31RDF	
PROTAC Binding Buffer 2	220 mL	64BDE32RDF	
PROTAC Binding Buffer 3	220 mL	64BDE33RDF	

Table 2. Cell based kits.

Product	HTRF	AlphaLISA
ACC Total	64NACPEG	ALSU-TACC-A500
AKT1 Total	63ADK079PEG	ALSU-TAKT-B-HV
AKT2 Total	63ADK081PEG	ALSU-TAKT-B-HV
AKT3 Total	63ADK083PEG	ALSU-TAKT-B-HV
AMPK/2 Total	63ADK060PEG	ALSU-TAMPK-A500
Androgen Receptor Total	64ANDRPEG	ALSU-TANDR-A500
ATF-4 Total		ALSU-TATF4-A500
ATM Total	64ATMTPEG	ALSU-TATM-A500
Aurora A Total		ALSU-TAURA-A500
Aurora B Total		ALSU-TAURB-A500
BCL-2		AL384C
Beta-Catenin Total	64NCAPEG	ALSU-TBCAT-A500
BRD4 Total	64BRD4TPEG	ALSU-TBRD4-A500
BRD2 Total	64BRD2TPEG	ALSU-TBRD2-A500
BRD3 Total	64BRD3TPEG	ALSU-TBRD3-A500
BRD7 Total	64BRD7TPEG	ALSU-TBRD7-A500
BRD9 Total	64BRD9TPEG	ALSU-TBRD9-A500
BTK Total	63ADK064PEG	ALSU-TBTK-A-HV
CaMKII Total		ALSU-TCAM-A500
CD3 Zeta Total		ALSU-TCD3-A500
CDK1 Total	64CDK1TPEG	ALSU-TCDK1-A500
CDK2 Total	64CDK2TPEG	ALSU-TCDK2-A500
CDK4 Total	64CDK4TPEG	ALSU-TCDK4-A500
CDK5 Total	64CDK5TPEG	ALSU-TCDK5-A500
CDK6 Total	64CDK6TPEG	ALSU-TCDK6-A500
CDK7 Total	64CDK7TPEG	ALSU-TCDK7-A500
CDK8 Total	64CDK8TPEG	
CDK9 Total	64CDK9TPEG	ALSU-TCDK9-A500
CDK12 Total	64CDK12TPEG	ALSU-TCDK12-A500
Cereblon total	64CRBNPEG	AL3139C
CHK1 Total	64CHK1TPEG	ALSU-TCHK1-A500
CHK2 Total	64CHK2TPEG	ALSU-TCHK2-A500
CREB Total	63ADK052PEG	ALSU-TCREB-A-HV
CREBBP Total		ALSU-TCRBP-A500
Cyclin B1 Total		ALSU-TCYCB1-A500
Cyclin D1 Total	64CYCD1TPEG	ALSU-TCYCD1-A500
C-Jun Total		ALSU-TCJUN-A500
c-Met Total		ALSU-TCMET-A500
c-Myb Total		ALSU-TCMYB-A500
c-Src Total	64NSRPEG	
DNA PKcs Total		ALSU-TDNA-A500
EGFR Total	64NG1PEG	ALSU-TEGFR-A-HV
eIF2a Total	64NEFPEG	ALSU-TEIF2-A-HV
ErbB2 (HER2) Total	64NR2PEG	ALSU-TEB2-A-HV
ErbB3 (HER3) Total	64NR3PEG	ALSU-TERB3-A500
ERK1/2 Total	64NRKPET	ALSU-TERK-A-HV

Table 2. Cell based kits. (continued).

Product	HTRF	AlphaLISA
ERα Total	64ERAPEG	ALSU-TERA-A500
FAK Total		ALSU-TFAK-A500
FGFR2 Total	63ADK057PEG	ALSU-TFGFR2-A500
FLT3 Total	64FLT3TPEG	ALSU-TFLT3-A500
FOXM1 Total		ALSU-TFOXM1-A500
FoxO1 Total	64NFOPEG	
GSK3α Total	64GTAPEG	
GSK3β Total	64GTBPEG	ALSU-TGS3B-A500
HDAC4 Total	64HDAC4TPEG	
HER3 Total	64NR3PEG	
Histone H3 Total	64NH3PEG	
HSP27 Total		ALSU-THSP-A500
H2AX Total	64H2AXTPEG	
IκBα Total		ALSU-TIKBA-A500
IKKA Total	64KKATPEG	
IKKβ Total	63ADK097PEG	ALSU-TIKKB-A500
InsR beta Total		ALSU-TINR-A500
IRF5 Total	64HIRF5PEG	
IRS1 Total	64NRSPEG	
IRS2 Total	64IRTPEG	
JAK1, Total	64JAK1TPEG	ALSU-TJAK1-A500
JAK2 Total	64JAK2TPEG	ALSU-TJAK2-A500
JAK3 Total	64JAK3TPEG	
LRRK2 Total	6FNRKPEG	
MDM2 Total		ALSU-TMDM2-A500
MEK1 Total	64NE1PEG	ALSU-TMEK1-A-HV
MKK4 Total	64NK4PEG	ALSU-TMKK4-A-HV
MLKL Total	64MLKLTPEG	ALSU-TMLKL-A500
NDRG1 Total	63ADK094PEG	
NFκB Total	64NFTPEG	ALSU-TNFKB-A500
NRF2 Total	64NRF2TPEG	ALSU-TNRF2-A500
PDHA1 Total		ALSU-TPDHA1-A500
PERK Total Kit		ALSU-TPERK-A500
Phospholamban Total	63ADK075PEG	
PLK1 Total		ALSU-TPLK1-A500
Progesterone Receptor Total	64HPRTPEG	
P21 Total	64HP21PEG	
p27 KIP1 Total		ALSU-TP27K-A500
P38 MAPK Total	64P38PET	ALSU-TP38-B500
p53 Total	64T53PET	ALSU-TP53-A-HV
p300 Total		ALSU-TP300-A500
PINK1 Total	64PINKTPEG	ALSU-TPINK-A500
RAB 8A Total		ALSU-TRAB8-A500
RAB10 Total	64RAB10TPEG	
RBM39 Total	64RBM39TPEG	ALSU-TRBM39-A500
RNA POL II Total		ALSU-TRNAP-A500

Table 2. Cell based kits. (continued).

Product	HTRF	AlphaLISA
SHP1 Total	64NH1PEG	ALSU-TSHP1-A-HV
SHP2 Total	64NH2PEG	ALSU-TSHP2-A-HV
SLP76 Total	63ADK077PEG	ALSU-TSLP-A-HV
SMAD3 Total	64ND3PEG	ALSU-TSM3-A500
SMAD4 Total	64SMAD4TPEG	ALSU-TSM4-A500
SMARCA2 Total	64SMARCA2TPEG	
SMARCA4 Total	64SMARCA4TPEG	ALSU-TBRG1-A500
SQSTM1 p62 Total		ALSU-TSQSTM-A500
STAT1 Total	63ADK096PEG	ALSU-TST1-A500
STAT3 Total	64NT3PEG	ALSU-TST3-A-HV
STAT6 Total	64STAT6TPEG	
SYK Total	64SYKTPEG	ALSU-TSYK-A-HV
Tau Total	64NTAUPEG	ALSU-TTAU-A500
TBK1 Total	64NTBPEG	ALSU-TTBK1-A500
TDP43 Aggr Total		ALSU-ATDP43-A500
TDP43 Total	64TDP4TPEG	ALSU-TTDP43-A500
TPD43 (Mouse) Total		ALSU-TTDP43-B500
TFEB Total	64TFEBTPEG	ALSU-TTFEB-A500
TYK2 Total	64TYK2TPEG	
VASP Total	63ADK067PEG	
ULK1 Total	64ULK1TPEG	ALSU-TULK1-A500
Wee1 Total	64WEE1TPEG	
XBP1 Total	64XBPTPEG	
YAP Total	64YATPEG	ALSU-TYAP-A500
ZAP70 Total	64ZATPEG	ALSU-TZAP-A-HV

Table 3. HTRF kinase binding kits.

Product	Part number
Kinase-6HIS Binding Discovery Kit	62KBD02PEA
Kinase-Biotin Binding Discovery Kit	62KBD03PEA
Kinase-GST Binding Discovery Kit	62KBD01PEA

Table 4. Alpha CETSA.

CETSA® is a registered trademark of Pelago Bioscience.

Product	Part number
	<b>500 Assay points</b>
Akt1	CETSA-AKT1-A500
BCL-2	CETSA-BCL2-A500
CDK2	CETSA-CDK2-A500
CDK4	CETSA-CDK4-A500
Cereblon Kit	CETSA-CRBN-A500
EGFR	CETSA-EGFR-A500
ERK1/2	CETSA-ERK-A500
ErbB2	CETSA-ERBB2-A500
MEK1	CETSA-MEK1-A500
NFκB p65	CETSA-NFKB-A500
p38	CETSA-p38-A500
Raf-1	CETSA-RAF1-A500
STING	CESTA-STING-A500

Table 5. HTRF labeled antibodies.

Product	Size	Part number
MAb Anti 6HIS-d2	1 000 tests	61HISDLF
MAb Anti 6HIS-Eu cryptate	5 000 tests	61HISKLA
MAb Anti 6HIS-Eu cryptate Gold	5 000 tests	61HI2KLA
MAb Anti 6HIS-Tb cryptate	1 000 tests	61HISTLF
MAb Anti 6HIS-Tb cryptate Gold	1 000 tests	61HI2TLF
MAb Anti 6HIS-XL665	1 000 tests	61HISXLF
MAb Anti cmc-d2	1 000 tests	61MYCDAF
MAb Anti cmc-Eu cryptate	5 000 tests	61MYCKLA
MAb Anti cmc-Tb cryptate	1 000 tests	61MYCTAF
MAb Anti cmc-XL665	1 000 tests	61MYCXLF
MAb Anti DNP-Eu cryptate	5 000 tests	61DNPKLA
MAb Anti DNP-XL665	5 000 tests	61DNPXLA
MAb Anti FLAG M2-d2	1 000 tests	61FG2DLF
MAb Anti FLAG M2-Eu cryptate	5 000 tests	61FG2KLA
MAb Anti FLAG M2-Tb cryptate	1 000 tests	61FG2TLF
MAb Anti FLAG M2-XL665	1 000 tests	61FG2XLF
MAb Anti GST-d2	1 000 tests	61GSTDLF
MAb Anti GST-Eu cryptate	5 000 tests	61GSTKLA

Table 5. HTRF labeled antibodies. (continued).

Product	Size	Part number
MAB Anti GST-Tb cryptate	1 000 tests	61GSTTLF
MAB Anti GST-XL665	1 000 tests	61GSTXLF
MAB Anti HA-d2	1 000 tests	610HADAF
MAB Anti HA-Eu cryptate	5 000 tests	610HAKLA
MAB Anti HA-Tb cryptate	1 000 tests	610HATAF
MAB Anti HA-XL665	1 000 tests	610HAXLF
MAB Anti MBP-d2	5 000 tests	61MBPDAA
MAB Anti MBP-Eu cryptate	5 000 tests	61MBPKAA
MAB Anti MBP-Tb cryptate	5 000 tests	61MBPTAA
PAb Anti Human IgG-Eu cryptate	5 000 tests	61HFCKLA
PAb Anti Human IgG-Tb cryptate	1 000 tests	61HFCTAF
PAb Anti Human IgG-d2	1 000 tests	61HFCDAF
PAb Anti Human IgG-XL665	1 000 tests	61HFCXLF
MAB Anti Human Kappa-Eu cryptate	5 000 tests	61KAPKAA
MAB Anti Human Kappa-d2	5 000 tests	61KAPDAA
Streptavidin-d2	1 000 tests	610SADLF
Streptavidin-XL665	1 000 tests	610SAXLF
Ubiquitin-Eu cryptate	5 000 tests	61UBIKLA

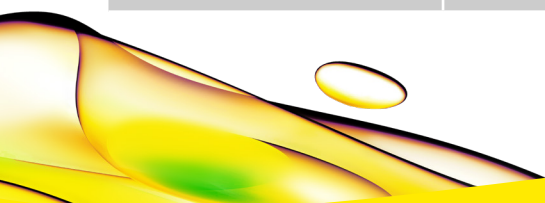
## Available products

Table 6. Alpha toolkit. (continued).

Alpha	Donor beads	Acceptor beads
Anti-6x His		AL178
Anti-chicken IgY		AL131
Anti-c-myc		AL111
Anti-DIG	AS108	AL113
Anti-DNP	AS111	AL173
Anti-FITC		AL127
Anti-FLAG®	AS103	AL112
Anti-GFP	AS112	AL133

Table 6. Alpha toolkit. (continued).

Alpha	Donor beads	Acceptor beads
Anti-GST		AL110
Anti-HA		AL170
Anti-HRP	AS109	AL171
Human IgG (Fc-fragment)	AS113	AL174
Anti-human IgG (Fc-specific)	AS114	AL103
Anti-human IgG1		AL141
Anti-human IgG2 (isotyping)		AL154
Anti-human IgG4		AL142
Anti-Maltose-Binding Protein (MBP)		AL134
Anti-mouse IgG	AS104	AL105
Anti-Protein C tag		AL172
Anti-rabbit IgG	AS105	AL104
Anti-rat IgG	AS110	AL106
Anti-sheep IgG		AL132
Anti-V5		AL129
Glutathione (GSH)	6765300/1/2/3	AL109
LCA (Lens Culinaris Agglutinin)		AL140
Nickel Chelate	AS101	AL108
Protein A	AS102	AL101
Protein G		AL102
Protein L		AL126
Strep Tactin®	AS106	AL136
Streptavidin	6760002	AL125



revvity