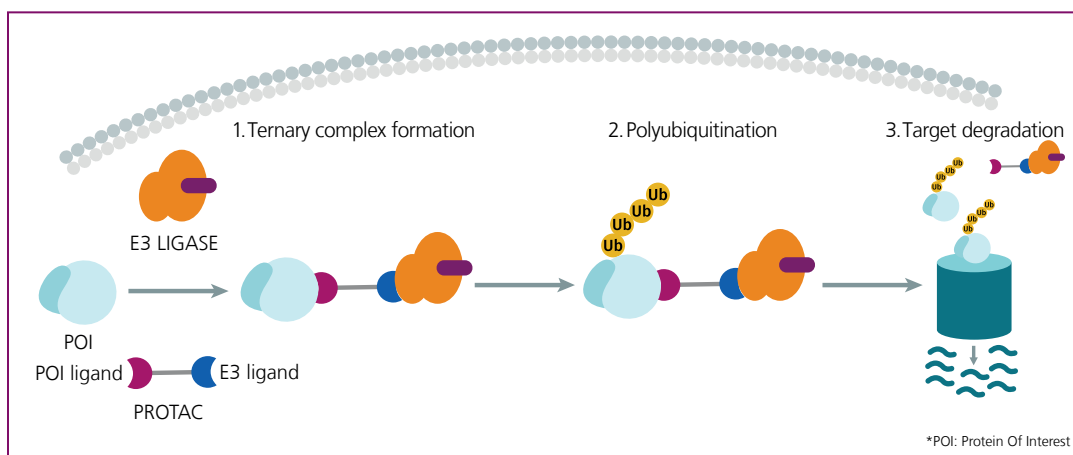


REAGENT SOLUTIONS FOR PROTAC RESEARCH



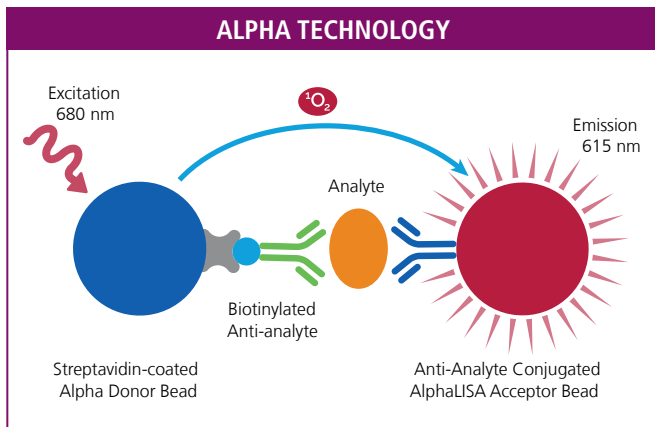
Proteolysis Targeting Chimaeras (PROTACs) is a game changing mechanism for treating the untreatable in the area of drug discovery.

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.

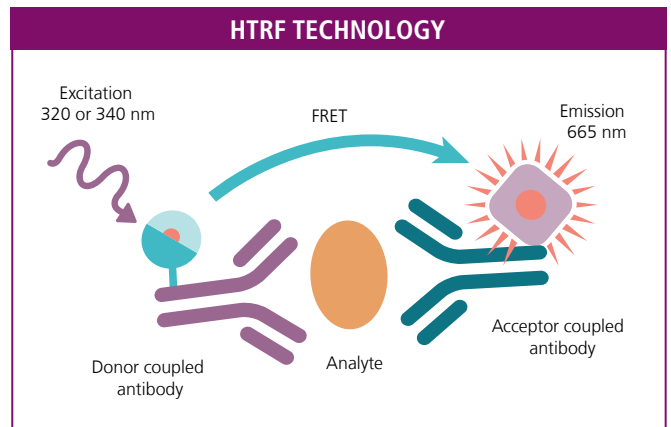
Contrary to current drugs which inhibit the proteasome, this mode of action leads to the destruction of the target. Consequently, it could be more effective than current cancer treatments.

Advance your PROTAC® research and benefit from PerkinElmer'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® technology to monitor binding of your compound of interest in the natural living cellular environment by bringing together the Cellular Thermal Shift assays (CETSA®) from Pelago Bioscience with PerkinElmer Alpha technology

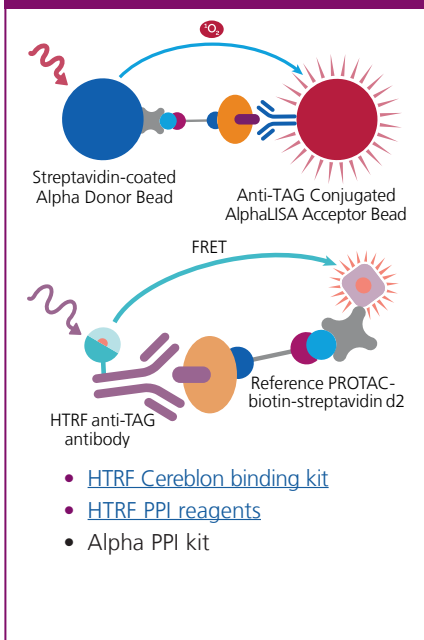


Alpha is a versatile, bead-based platform that enables you to assay the most complex samples in one well and with no wash steps.

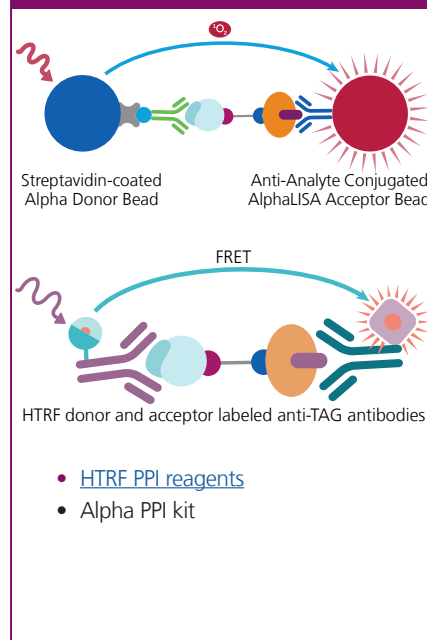


HTRF is a fast, sensitive, homogeneous, and ready to use assay platform with no wash steps.

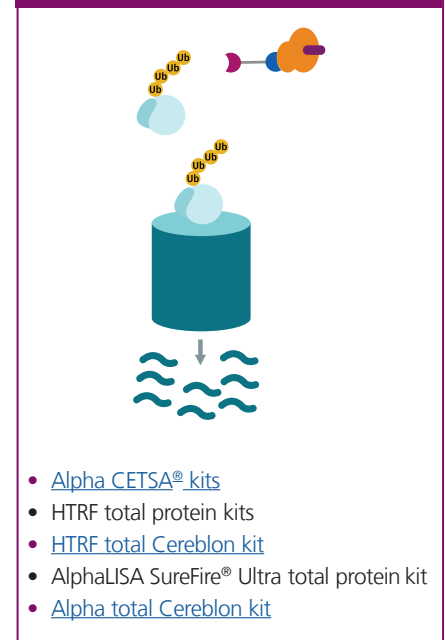
Does my PROTAC bind to E3 ligase?



Is a ternary complex formed?



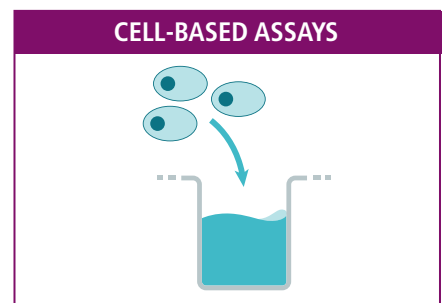
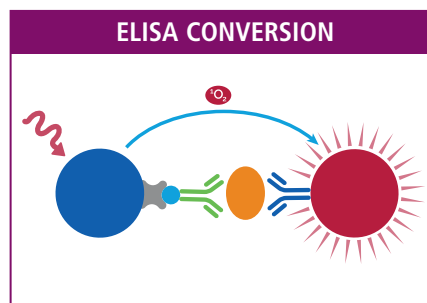
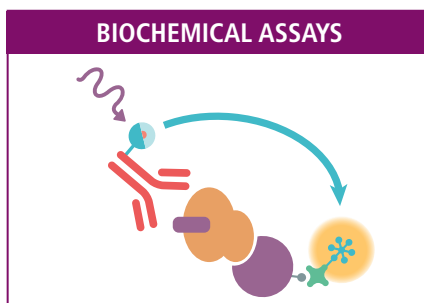
Does my PROTAC compound bind and degrade my target?



Custom Labeling and Custom Assay Development Services

Don't see the assay you need listed here? PerkinElmer 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.



Contact us today: www.cisbio.com/content/assay-development

Table 1. 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 cmyc-d2	1 000 tests	61MYCDAF
MAb Anti cmyc-Eu cryptate	5 000 tests	61MYCKLA
MAb Anti cmyc-Tb cryptate	1 000 tests	61MYCTAF
MAb Anti cmyc-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
MAb Anti GST-Tb cryptate	1 000 tests	61GSTTLF
MAb Anti GST-XL665	1 000 tests	61GSTXLF
MAb Anti HA-d2	1 000 tests	61OHADAF
MAb Anti HA-Eu cryptate	5 000 tests	61OHAKLA
MAb Anti HA-Tb cryptate	1 000 tests	61OHATAF
MAb Anti HA-XL665	1 000 tests	61OHAXLF
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

Table 2. Cell based kits.

Product	AlphaLISA	HTRF
ACC Total	ALSU-TACC-A500	64NACPEG
AKT1 Total	ALSU-TAKT-B-HV	63ADK079PEG
AKT2 Total	ALSU-TAKT-B-HV	63ADK081PEG
AKT3 Total	ALSU-TAKT-B-HV	63ADK083PEG
AMPK/2 Total	ALSU-TAMPK-A500	63ADK060PEG
BCL-2	AL384	
Beta-Catenin Total		64NCAPEG
BTK Total	ALSU-TBTK-A-HV	63ADK064PEG
CDK2 Total	ALSU-TCDK2-A500	64CDK2TPEG
CDK4 Total	ALSU-TCDK4-A500	64CDK4TPEG
Cereblon total	AL3139C	64CRBNPEG
CREB Total	ALSU-TCREB-A-HV	63ADK052PEG
c-Src Total		64NSRPEG
EGFR Total	ALSU-TEGFR-A-HV	64NG1PEG
eIF2a Total	ALSU-TEIF2-A-HV	64NEFPEG
ErbB2 (HER2) Total	ALSU-TEB2-A-HV	64NR2PEG
FGFR2 Total	ALSU-TFGFR2-A500	63ADK057PEG
FoxO1 Total		64NFOPEG
GSK3 α Total		64GTAPEG
GSK3 β Total		64GTBPEG
HER3 Total		64NR3PEG
Histone H3 Total		64NH3PEG
Human Tau Total	AL271C	64NTAUPEG
IKKB Total		63ADK097PEG
InsR beta Total	ALSU-TINR-A500	
IRS1 Total		64NRSPEG
IRS2 Total		64IRTPEG
LRRK2 Total		6FNRKPEG
MEK1 Total	ALSU-TMEK1-A-HV	64NE1PEG
MKK4 Total	ALSU-TMKK4-A-HV	64NK4PEG
NDRG1 Total		63ADK094PEG
NF κ B Total	ALSU-TNF κ B-A500	64NFTPEG
P38 MAPK Total	ALSU-TP38-B500	64P38PET
p53 Total	ALSU-TP53-A-HV	64T53PET
Phospholamban Total		63ADK075PEG
Raf1	ALSU-RAF1-A500	
Rb Total	ALSU-TRB-A500	63ADK106PEG
SHP1 Total	ALSU-TSHP1-A-HV	64NH1PEG
SHP2 Total	ALSU-TSHP2-A-HV	64NH2PEG
SLP76 Total	ALSU-TSLP-A-HV	63ADK077PEG
SMAD3 Total	ALSU-TSM3-A500	64ND3PEG
STAT1 Total	ALSU-TST1-A500	63ADK096PEG
STAT3 Total	ALSU-TST3-A-HV	64NT3PEG
SYK Total	ALSU-TSYK-A-HV	64SYKTPEG
TBK1 Total	ALSU-TTBK1-A500	64NTBPEG
Total-ERK1/2	ALSU-TERK-A-HV	64NRKPET
VASP Total		63ADK067PEG
YAP Total		64YATPEG
ZAP70 Total	ALSU-TZAP-A-HV	64ZATPEG

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.

Product	Part Number
500 Assay Points	
Akt1	CETSA-AKT1-A500
BCL-2	CETSA-BCL2-A500
CDK2	CETSA-CDK2-A500
CDK4	CETSA-CDK4-A500
EGFR	CETSA-EGFR-A500
ERK1/2	CETSA-ERK-A500
ErbB2	CETSA-ERBB2-A500
MEK1	CETSA-MEK1-A500
NF κ B p65	CETSA-NF κ B-A500
p38	CETSA-p38-A500
Raf-1	CETSA-RAF1-A500

CETSA® is a registered trademark of Pelago Bioscience.

Available products

Table 5. Alpha toolkit.

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
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

FOR MORE INFORMATION

Europe and other countries +33(0)466 796 705 U.S. and Canada 1-888-963-4567 China +86 21 5018 9880 Japan +81-(0)43-306-8712
Visit www.cisbio.com to find a list of our regional distributors

PerkinElmer, Inc.
940 Winter Street
Waltham, MA 02451 USA
P: (800) 762-4000 or
(+1) 203-925-4602
www.perkinelmer.com



cisbio

For a complete listing of our global offices, visit www.perkinelmer.com/ContactUs

Copyright ©2020, PerkinElmer, Inc. All rights reserved. PerkinElmer® is a registered trademark of PerkinElmer, Inc. All other trademarks are the property of their respective owners.

145833 (105765) PKI