

NEW PRODUCTS TO MEET YOUR RESEARCH AND DRUG DISCOVERY NEEDS

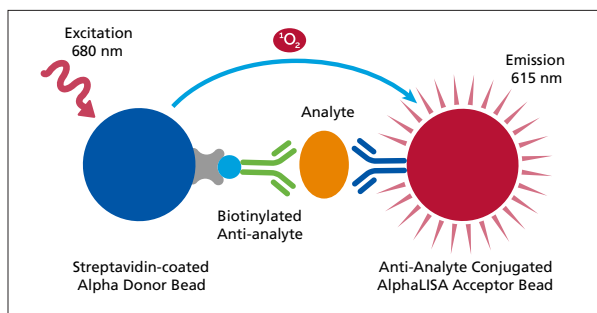
AlphaLISA and HTRF As your science evolves, our product offering expands!

Our R&D scientists have been working tirelessly to bring you fully validated and ready-to-use immunoassay Kits for the new targets you need most.

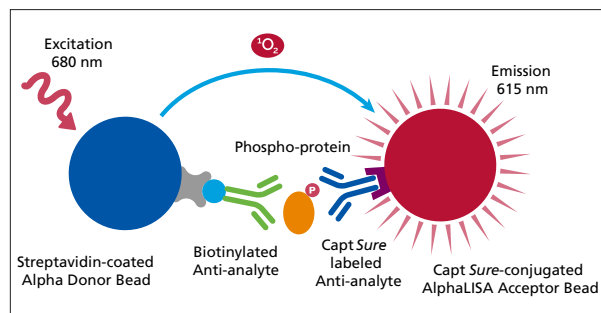
AlphaLISA and HTRF no-wash technologies provide a larger dynamic range, improved sensitivity, and higher signal-to-noise ratios than traditional wash-based ELISAs.

We are pleased to announce our newest AlphaLISA and HTRF assays. Benefit from PerkinElmer's broad offering of reagent technologies and ditch your ELISA wash steps for good!

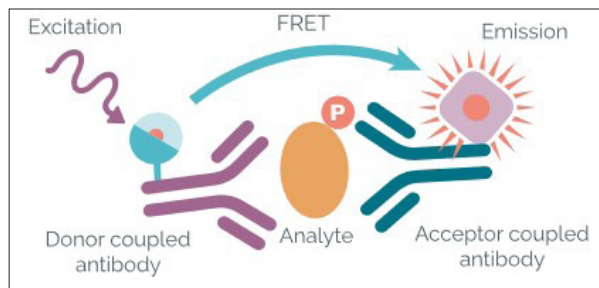
AlphaLISA® technology



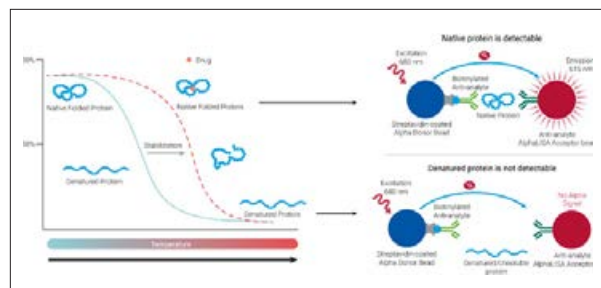
AlphaLISA® SureFire® Ultra™ technology



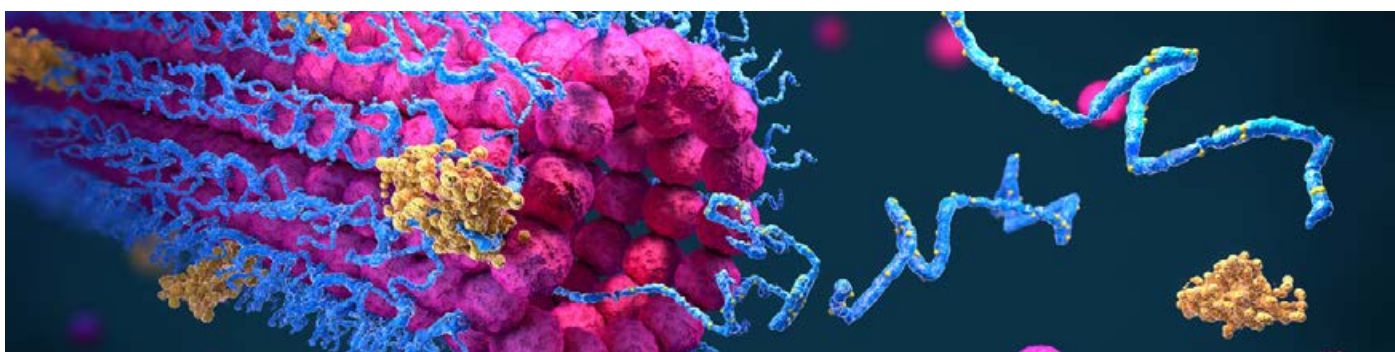
HTRF® technology



Alpha® CETSA



For more information please visit: www.perkinelmer.com/newestassays



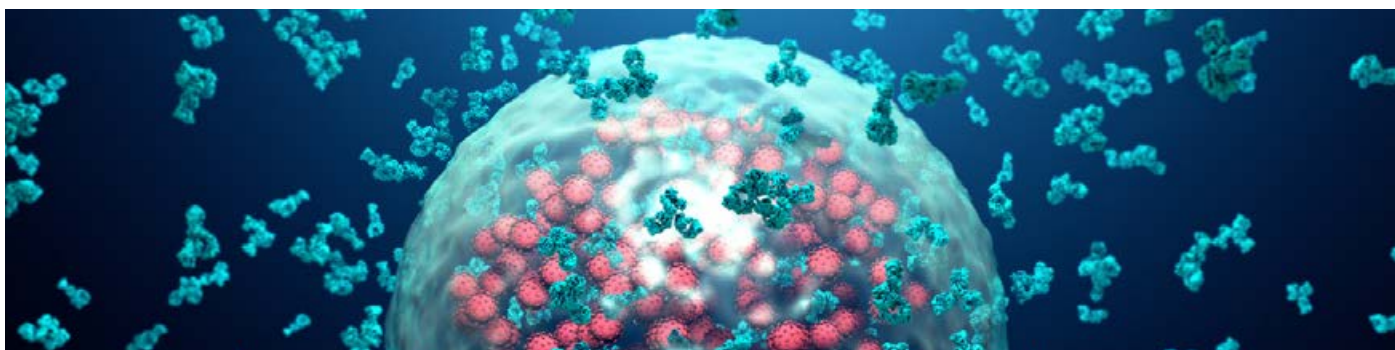
ORDERING INFORMATION – MAY 2022

Biomarkers Assays

AlphaLISA PRODUCTS	Part - 500 tests
AAV1 Capsid Kit	AL3181C
Insulin Human and Mouse Kit	AL3184C
TOM20 Kit	AL3178C

Cellular Phospho-Protein Assays

AlphaLISA SureFire Ultra PRODUCTS	Part - 500 tests
ErbB3, Phospho-Y1289 Kit	ALSU-PERB3-A500
ErbB3, Total Kit	ALSU-TERB3-A500
PLK1, Phospho-T210 Kit	ALSU-PPLK1-A500
PLK1, Total Kit	ALSU-TPLK1-A500
SMAD4, Total Kit	ALSU-TSM4-A500
SMARCA4, Total Kit	ALSU-TBRG1-A500
HTRF PRODUCTS	Part - 500 tests
BRD2, Total Kit	64BRD2TPEG
BRD3, Total Kit	64BRD3TPEG
CHK1, Phospho-S345 Kit	64CHK1S5PEG
CHK1, Total Kit	64CHK1TPEG
CHK2, Phospho-T68 Kit	64CHK2T8PEG
CHK2, Total Kit	64CHK2TPEG



ORDERING INFORMATION – MARCH 2022

Biomarkers Assays

HTRF PRODUCTS	Part - 500 tests
HTT, Total Kit	64HTTTPEG
HTT, Mutant Kit	64HTTMPEG

Cellular Phospho-Proteins Assays

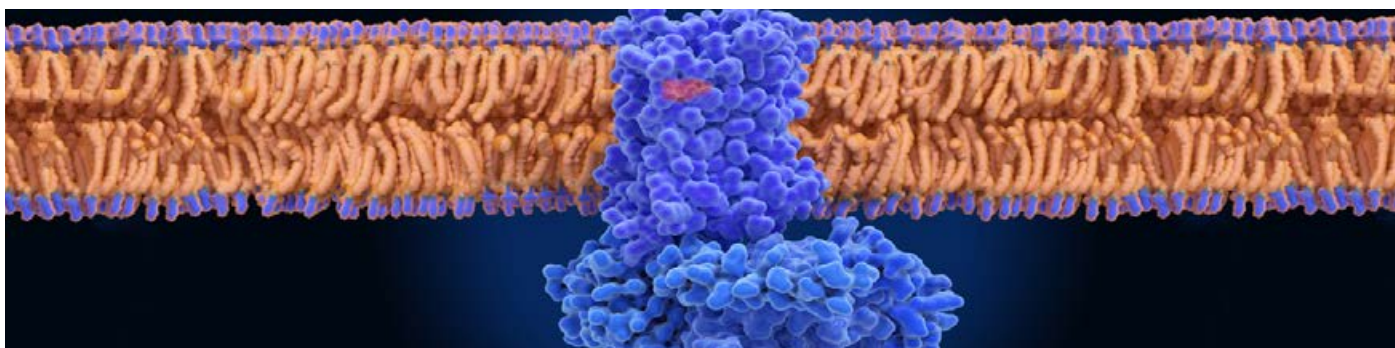
p21, Total Kit	64HP21PEG
SMAD4, Phospho-T277 Kit	64SMAD4T7PEG
SMAD4, Total Kit	64SMAD4TPEG

Cellular Phospho-Proteins Assays

AlphaLISA <i>SureFire Ultra</i> PRODUCTS	Part - 500 tests
BRD4, Total Kit	ALSU-TBRD4-A500
BRD9, Total Kit	ALSU-TBRD9-A500
CDK1, Phospho-T14 Kit	ALSU-PCDK1-A500
CDK1, Phospho-Y15 Kit	ALSU-PCDK1-B500
CDK1, Phospho-T161 Kit	ALSU-PCDK1-C500
CDK1, Total Kit	ALSU-TCDK1-A500
CDK7, Total Kit	ALSU-TCDK7-A500
CDK9, Total Kit	ALSU-TCDK9-A500
Cyclin D1, Phospho-T286 Kit	ALSU-PCYCD1-A500
Cyclin D1, Total Kit	ALSU-TCYCD1-A500
FLT3, Phospho-Y591 Kit	ALSU-PFLT3-A500
FLT3, Phospho-Y842 Kit	ALSU-PFLT3-B500
FLT3, Total Kit	ALSU-TFLT3-A500
PERK, Phospho-T980 Kit	ALSU-PPERK-A500
PERK, Total Kit	ALSU-TPERK-A500
YAP, Total Kit	ALSU-TYAP-A500
HTRF PRODUCTS	Part - 500 tests
BRD4, Total Kit	64BRD4TPEG
BRD7, Total Kit	64BRD7TPEG
BRD9, Total Kit	64BRD9TPEG
CDK1, Total Kit	64CDK1TPEG
CDK7, Total Kit	64CDK7TPEG
CDK9, Total Kit	64CDK9TPEG
CDK12, Total Kit	64CDK12TPEG
Cyclin D1, Phospho-T286 Kit	64CYCD1T6PEG
Cyclin D1, Total Kit	64CYCD1TPEG
IRF5, Total Kit	64HIRF5PEG
MEK1, Phospho-S298 Kit	64MEK1S98PEG

Bioprocess Assays

AlphaLISA PRODUCTS	Part - 500 tests
CHO HCP Kit	AL3176C
HTRF PRODUCTS	Part - 500 tests
CHO HCP Kit	64CHOPEG

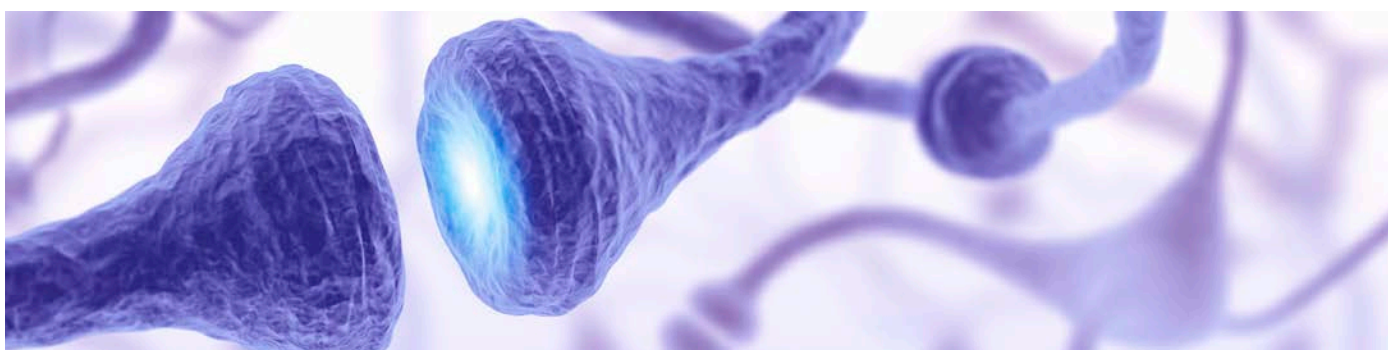


ORDERING INFORMATION – JANUARY 2022

Biomarkers Assays	
AlphaLISA PRODUCTS	Part - 500 tests
Fibronectin Human and Mouse Kit	AL3174C
Fibronectin-EDA Human Kit	AL3175C
HTRF PRODUCTS	Part - 500 tests
ApoE Human Kit	64APOEPEG
Estrogen Receptor Human Kit	64ERAPEG
Fibronectin Human and Mouse Kit	64HMFNPEG
Fibronectin-EDA Human Kit	64HFNEDAPEG
Flag Check Kit	63ADK036PEG
TYK2 Total Kit	64TYK2TPEG

Protein-Protein Interactions Assays	
AlphaLISA PRODUCTS	Part - 500 tests
C1Q Human Binding Kit	AL3169C
XIAP BIR2 Human Binding Kit	AL3172C
XIAP BIR3 Human Binding Kit	AL3173C
HTRF PRODUCTS	Part - 500 tests
VHL Human Binding Kit	64BDVHLEPEG
VHL Red Ligand	64BDVHLRED

Cellular Phospho-Protein Assays	
AlphaLISA SureFire Ultra PRODUCTS	Part - 500 tests
c-Jun, Phospho-S63 Kit	ALSU-PCJUN-A500
c-Jun, Phospho-S73 Kit	ALSU-PCJUN-B500
c-Jun, Total Kit	ALSU-TJUN-A500
DNA PKcs, Phospho-S2056 Kit	ALSU-PDNA-A500
DNA PKcs, Total Kit	ALSU-TDNA-A500
JAK1, Total Kit	ALSU-TJAK1-A500
ULK1, Phospho-S556 Kit	ALSU-PULK1-A500
ULK1, Phospho-S758 Kit	ALSU-PULK1-B500
ULK1, Total Kit	ALSU-TULK1-A500
HTRF PRODUCTS	Part - 500 tests
TYK2, Phospho-T1054/1055 Human Kit	64TYK2Y10PEG
TYK2, Total Kit	64TYK2TPEG



ORDERING INFORMATION – OCTOBER 2021

Cellular Phospho-Proteins Assays

AlphaLISA® SureFire® Ultra™	Part - 500 tests
AKT 1/2/3, Phospho-T450 Kit	ALSU-PAKT-C500
ALK, Phospho-Y1604 Kit	ALSU-PALK-C500
CaMKII, Phospho-T286 Kit	ALSU-PCAM-A500
CaMKII, Total Kit	ALSU-TCAM-A500
CD3 Zeta, Phospho-Y142 Kit	ALSU-PCD3-B500
CD3 Zeta, Phospho-Y83 Kit	ALSU-PCD3-A500
CD3 Zeta, Total Kit	ALSU-TCD3-A500
c-Myb, Phospho-S11 Kit	ALSU-PCMYB-A500
c-Myb, Total Kit	ALSU-TCMYB-A500
ERα, Phospho-S118 Kit	ALSU-PERA-A500
ERα, Phospho-S167 Kit	ALSU-PERA-B500
ERα, Total Kit	ALSU-TERA-A500
IκBα, Phospho-S32/36 Kit	ALSU-PIKBA-A500
IκBα, Total Kit	ALSU-TIKBA-A500
IKKβ, Phospho-S177/181 Kit	ALSU-PIKKB-A500
IKKβ, Total Kit	ALSU-TIKKB-A500
PDHA1, Phospho-S293 Kit	ALSU-PPDHA1-A500
PDHA1, Total Kit	ALSU-TPDHA1-A500
SQSTM1 p62, Phospho-S403 Kit	ALSU-PSQSTM-A500
SQSTM1 p62, Total Kit	ALSU-TSQSTM-A500
TDP43, Phospho-S409/410 Kit	ALSU-PTDP43-A500
TDP43, Total Kit	ALSU-TTDP43-A500

Biomarkers Assays

AlphaLISA PRODUCTS	Part - 500 tests
Granzyme B Kit	AL3170C

Cellular Phospho-Proteins Assays

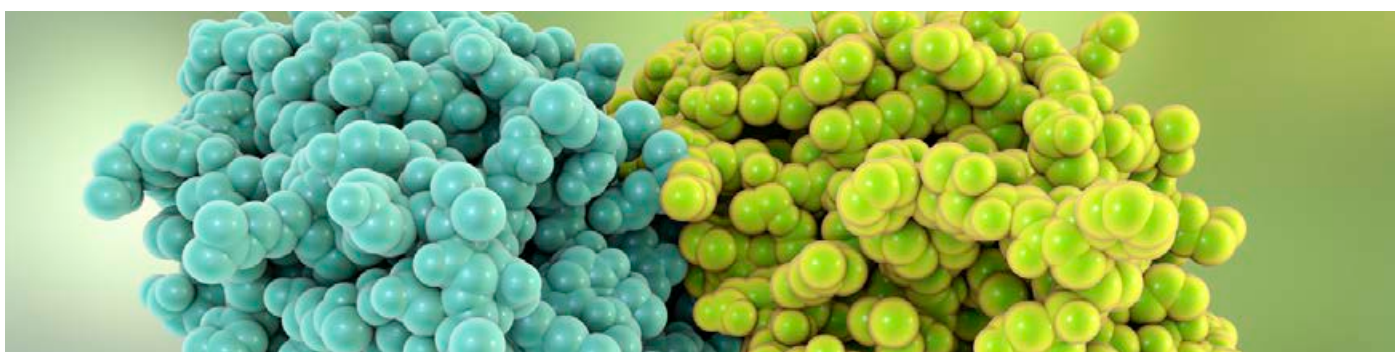
HTRF PRODUCTS	Part - 500 tests
Androgen Receptor Kit	64ANDRPEG
JAK1, Phospho-Y1034/1035 Kit	64JAK1Y10PEG
JAK1, Total Kit	64JAK1TPEG
JAK2, Total Kit	64JAK2TPEG
JAK2, Phospho-Y1007/1008 Kit	64JAK2Y10PEG
JAK3, Total Kit	64JAK3TPEG
AlphaLISA SureFire Ultra Multiplex PRODUCTS	Part - 500 tests
AKT 1/2/3, Phospho-T308 + AKT 1/2/3, Total Kit	MPSU-PTAKT-O500
AKT 1/2/3, Phospho-T450 + AKT 1/2/3, Total Kit	MPSU-PTAKT-N500
ALK, Phospho-Y1604 + ALK, Total Kit	MPSU-PTALK-K500
Ret, Phospho-panY + Ret, Total Kit	MPSU-PTRET-K500
SLP-76, Phospho-S376 + SLP-76, Total Kit	MPSU-PTSLP-K500

PROTAC biochemical Assays

AlphaLISA PRODUCTS	Part - 500 tests
cIAP-BIR2 Binding Kit	AL3166C
cIAP-BIR3 Binding Kit	AL3167C
MDM2 Binding Kit	AL3168C
HTRF PRODUCTS	Part - 500 tests
XIAP-Red Ligand	64APOEPEG
XIAP-BIR2 Binding Kit	64XIAPB2PEG
XIAP-BIR3 Binding Kit	64XIAPB3PEG

HTRF® products also available under 10 000 tests & bulk size
AlphaLISA® products also available under 100 – 10 000 – 50 000 tests

For more information please visit:
www.perkinelmer.com/newestassays



ORDERING INFORMATION – JULY 2021

Cellular Phospho-Proteins Assays

AlphaLISA SureFire Ultra PRODUCTS	Part - 500 tests
ATG16L1, Phospho-S278 Kit	ALSU-PATG-A500
ATG16L1, Total Kit	ALSU-TATG-A500
β -Catenin, Phospho-S37 Kit	ALSU-PBCAT-A500
β -Catenin, Phospho-T41/S45 Kit	ALSU-PBCAT-B500
β -Catenin, Phospho-S675 Kit	ALSU-PBCAT-C500
β -Catenin, Total Kit	ALSU-TBCAT-A500
Histone H3, Phospho-S10/T11 Kit	ALSU-PHISH3-A500
Histone H3, Total Kit	ALSU-THISH3-A500
RAB10, Phospho-T73 Kit	ALSU-PRAB-A500
RAB10, Total Kit	ALSU-TRAB-A500
SLP76, Phospho-S376 Mouse Kit	ALSU-PSLP-D500
SLP76, Total Mouse Kit	ALSU-TSLP-B500
HTRF PRODUCTS	Part - 500 tests
cRaf, Phospho-S43 Kit	64CRAFS43PEG
cRaf, Total Kit	64CRAFTPEG
MLKL, Phospho-S358 Kit	64MLKLS8PEG
MLKL, Total Kit	64MLKLTPEG
RET, Phospho-pan Kit	64RETYPEG
RET, Total Kit	64RETTPEG
RSK1, Phospho-T359/S363 Kit	64RSK1T9PEG
RSK1, Phospho-S380 Kit	64RSK1S0PEG
RSK1, Total Kit	64RSK1TPEG

Protein-Protein Interactions Assays

AlphaLISA PRODUCTS	Part - 500 tests
KRAS G12C GTP Kit	AL3149C
KRAS G12C/SOS1 Kit	AL3151C
KRAS WT GTP Kit	AL3148C
KRAS WT/SOS1 Kit	AL3150C

Biomarkers Assays

HTRF PRODUCTS	Part - 500 tests
TOM20 Kit	64TOM20PEG
AlphaLISA PRODUCTS	Part - 500 tests
TREM2 Human Kit	AL3152C

PROTAC biochemical Assays

HTRF PRODUCTS	Part - 500 tests
cIAP1 BIR2 Binding Kit	64BDIAPB2PEG
cIAP1 BIR3 Binding Kit	64BDIAPB3PEG
LCL161 - Red Ligand	64BDIAPB3RED
MDM2 Binding Kit	64BDMDM2PEG
MI-1061 - Red Ligand	64BDMDM2RED
PROTAC Binding Buffer 2	64BDE32RDF
PROTAC Binding Buffer 3	64BDE33RDF

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



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

Copyright © 2022, PerkinElmer, Inc. All rights reserved. PerkinElmer® is a registered trademark of PerkinElmer, Inc. All other trademarks are the property of their respective owners.

695753 (614272) PKI