

List of Target Proteins Validated for NanoBRET™ Target Engagement Assay

Updated August 2023 (Targets in blue are available via Early Access)

E3 Ubiquitin Ligases	OPRK1	PARP1
CRBN	OPRM1	PARP2
VHL	ADRA1A	PARP3
cIAP1	ADRA1B	PARP4
cIAP2	ADRA2A	PARP5a
XIAP	ADRA2B	PARP5b
MDM2	ADRA2C	PARP6
Bromodomains (BRDs)	DRD1	PARP7
BRD2	DRD2	PARP8
BRD2 BD1	DRD3	PARP10
BRD2 BD2	DRD4	PARP11
BRD3	DRD5	PARP12
BRD4	HRH1	RAF (Protomer & Homodimer)
BRD4 BD1	HTR1A	ARAF Protomer
BRD4 BD2	HTR1B	BRAF Protomer
BRDT	HTR1F	CRAF Protomer
BRD3 BD1	HTR2A	BRAF Homodimer
BRD3 BD2	HTR2C	CRAF Homodimer
BRD7 BD	HTR6	RAS
BRD9	Heat Shock Protein (HSP)	KRAS2B WT
BRD9 BD	HSP90AB1	KRAS2B G12C
BRPF1	Histone Deacetylase (HDAC)	KRAS2B G12A
BRPF1 BD	HDAC1	KRAS2B G12D
CECR2	HDAC2	KRAS2B G12S
CECR2 BD	HDAC3	KRAS2B G12V
TAF1	HDAC6 (FL, CD2)	KRAS2B G12R
TAF1 BD1-BD2	HDAC10	KRAS2B G13D
TAF1 BD2	Kinases	KRAS2B Q61H
TAF1L	>340 targets in catalog (see full list here)	KRAS2B Q61L
G-protein Coupled Receptors (GPCR)	Additional target available in Early Access (see next page)	KRAS2B Q61R
ADORA1	NLR	HRAS WT
ADORA2A	NLRP3	HRAS G12C
ADORA2B	Poly (ADP-Ribose) Polymerase (PARP)	HRAS G12V
ADORA3		
OPRD1		

NanoBRET™ Target Engagement

Kinase Targets in Early Access

Kinase/Kinase Complex		
AKT3 (isoform 1)*	EML4-ALK	NPM1-ALK
AKT3 (isoform 2)*	EPHA10	PDPK1
ALK	EPHB6	PI4KA
ALK(C1156Y)	FGFR2(V564F)	PI4KB
ALK(L1196M)	FGFR2(N549K)	PIK3CG*
AXL(R492C)	FGFR2(N549H)	PKN1
BRAF	FGFR2(L617F)	PKN2
BTK(L528W)	FGFR2(L617V)	PKN3*
BTK(M437R)	FGFR2(K659M)	PLK1
BTK(V416L)	FGFR2(M537I)	PNCK*
BTK(T474I)	FGFR3(K650E)	PRKCA + PMA stim.
CAMK2B	FGFR3(L608V)	PRKCB + PMA stim.*
CASK	FGFR3(N540K)*	PRKCD + PMA stim.
CDK1/Cyclin A2	FGFR3(V553M)	PRKCG + PMA stim.
CDK2/Cyclin A2	FGFR3(V555F)	PRKCH + PMA stim.
CDK3/Cyclin E2	FGFR3(V555M)	PRKCQ + PMA stim.
CDK6/Cyclin D3	GRK1	RIPK3
CDK12/Cyclin K*	GRK7	ROCK1
CDK13/Cyclin K*	HIPK1*	ROCK2
CIT*	HUNK	ROS1*
CRAF	JAK1*	ROS1(1883-2347)
DAPK3*	JAK1 (JH1 domain)	RPS6KB1*
DCLK1	KDR*	RPS6KB1 + KRAS(G12C)
DCLK1 (isoform 2)	MAP2K1 + KRAS(G12C)*	SGK3
DSTYK*	MAP2K2 + KRAS(G12C)*	TGFBFR1
DYRK3*	MAP2K5	TNIK*
EGFR(L858R,T790M)	MAPK15	TNIK + KRAS(G12C)
	MAPK7(1-713)	UHMK1*

* = Targets with low assay window. For information on the capabilities of low assay window assays, see Technical Manual TM598 or contact your local Promega representative.

Do not see your target of interest listed?

Explore the Do-It-Yourself option using the NanoBRET™ 590 dyes to design novel NanoBRET™ tracers against your target of interest. Or work with Promega's Tailored R&D Solutions team for new assay development.

[NanoBRET™ 590 Dyes](#)

[Custom Assay Development](#)