

ProQinase™ RAF1 Y340D/Y341D (GST-HIS-tag)

Raf-1 proto-oncogene, serine/threonine kinase

Recombinant Human Active Protein Kinase

HGNC Symbol: RAF1

Synonyms: cRAF, CRAF, NS5

Product No.: 0792-0000-1

Lot: 001

Description: Human RAF1 C-terminal fragment, amino acids S₃₀₆-F₆₄₈ (as in [NCBI/Protein](#) entry NP_002871.1), Y340D and Y341D mutant, N-terminal GST-HIS₆ fusion protein with a 3C cleavage site, expressed in Sf9 insect cells

Product identity: RAF1 Y340D/Y341D Lot 001 was confirmed as RAF1 by mass spectroscopy LC-ESI-MS/MS

Theoretical MW_{Fusion Protein}: 67,409 Da

Expression host: Sf9 insect cells

Purification: GST-Affinity Chromatography

Activation: This kinase was not activated by special procedures

Storage buffer: 50 mM HEPES pH 7.5, 100 mM NaCl, 5 mM DTT, 15 mM reduced glutathione, 20 % glycerol

Storage temperature: -80°C

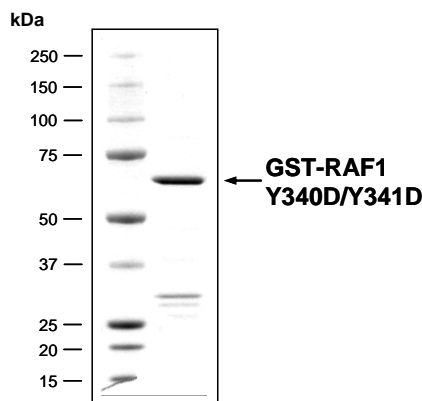
For complete recovery, mix well and spin before use. Product must not be stored in diluted solutions, aliquots below 10µl are not advisable. Avoid repeated freeze-thaw cycles!

Protein concentration: 0.301 µg/µl
(Bradford method using BSA [Sigma, cat# A-7638, Lot 79H7641] as standard protein)

Biochemical Parameters:
Specific kinase activity (P_i transfer): 124 pmol/µg × min
ATP-K_M: 0.32 µM

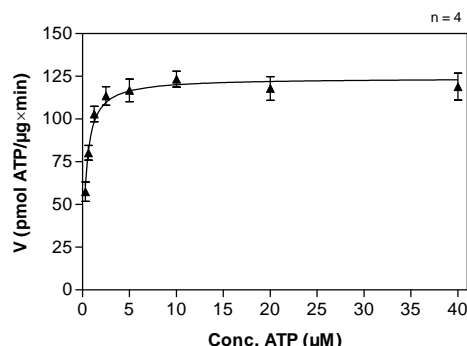
Additional assay technology:
RAF1 YDYD Lot 001 was also successfully tested by Reaction Biology for the use with the ADP-Glo™ Kinase assay from Promega ADP-Glo assay conditions may vary from radiometric assay conditions, please inquire for assay details

RAF1 YDYD Lot 001: Coomassie stain



2.0 µg GST-RAF1 Y340D/Y341D

RAF1 YDYD Lot 001: Determination of V_{max} and K_M value for ATP



Determination of K_M value & Specific activity:

- Assay conditions:
 - 60 mM HEPES-NaOH, pH 7.5
 - 3 mM MgCl₂
 - 3 mM MnCl₂
 - 3 µM Na-orthovanadate
 - 1.2 mM DTT
 - 50 µg/ml PEG_{20,000}
 - ATP (variable)
 - Substrate: MEK1-KM, 40 µg/ml
 - Kinase: 400 ng/ml
- Filter binding assay
- MSFC membrane (Millipore)

ProQinase™ RAF1 Y340D/Y341D (GST-HIS-tag)

Product No.: 0792-0000-1

GST-RAF1 YDYD Recombinant Fusion Protein Amino Acid Sequence							
1	MSPILGYWKI	KGLVQPTRL	LEYLEEKYEE	HLYERDEGDK	WRNKKFELGL	EFPNLPYYID	60
61	GDVKLTQSMA	IIRYIADKHN	MLGGCPKERA	EISMLEGAVL	DIRYGVSRIA	YSKDFETLKV	120
121	DFLSKLPPEML	KMFEDRLCHK	TYLNGDHVTH	PDFMLYDALD	VVLYMDPMCL	DAFPKLVCFK	180
181	KRIEAIPOID	KYLKSSKYIA	WPLQGWAQTF	GGGDHPPKSD	PMGHHHHHG	RDSLEVLFGG	240
241	PLAMGSQPKT	PVPAQRERAP	VSGTQEKNKI	RPRGQRDSSD	DWEIEASEVM	LSTRIGSGSF	300
301	GTVYKQKWHG	DVAVKILKVV	DPTPEQFQAF	RNEVAVLRKT	RHVNILLFMG	YMTKDNLAIV	360
361	TQWCEGSSLY	KHLHVQETKF	QMFQLIDIAR	QTAQGMDYLH	AKNIIHRDMK	SNNIFLHEGL	420
421	TVKIGDFGLA	TVKSRWGSQ	QVEQPTGSVL	WMAPEVIRMQ	DNNPFSFQSD	VYSYGIVLYE	480
481	LMTGELPYSH	INNDRQIIFM	VGRGYASPD	SKLYKNCPKA	MKRLVADCVK	KVKEERPLFP	540
541	QILSSIELQ	HSLPKINRSA	SEPSLHRAAH	TEDINACTLT	TSPRLPVF		600

1-218: GST **Red**: HIS6-tag **Green**: 3C cleavage site **blue**: RAF1 fragment **DD**: YDYD mutation

RAF1 wt ¹ Amino Acid Sequence							
1	MEHIQGAWKT	ISNGFGFKDA	VFDGSSCISP	TIVQQFGYQR	RASDDGKLT	PSKTSNTIRV	60
61	FLPNKQRTVV	NVRNGMSLHD	CLMKALKVRG	LQPECCAVER	LLHEHKGKKA	RLDWNTDAAS	120
121	LIGEELQVDF	LDHVPLTTHN	FARKTFLKLA	FCDICQKFL	NGFRCQTCGY	KFHEHCSTKV	180
181	PTMCDVWSNI	RQLLLFPNST	IGDSGVPALP	SLTMRRMRES	VSRMPVSSQH	RYSTPHAFTF	240
241	NTSSPSSEGS	LSQRQRSTST	PNVHVMVSTL	PVDSRMIEDA	IRSHSESASP	SALSSSPNNL	300
301	SPTGWSQPKT	PVPAQRERAP	VSGTQEKNKI	RPRGQRDSSY	YWEIEASEVM	LSTRIGSGSF	360
361	GTVYKQKWHG	DVAVKILKVV	DPTPEQFQAF	RNEVAVLRKT	RHVNILLFMG	YMTKDNLAIV	420
421	TQWCEGSSLY	KHLHVQETKF	QMFQLIDIAR	QTAQGMDYLH	AKNIIHRDMK	SNNIFLHEGL	480
481	TVKIGDFGLA	TVKSRWGSQ	QVEQPTGSVL	WMAPEVIRMQ	DNNPFSFQSD	VYSYGIVLYE	540
541	LMTGELPYSH	INNDRQIIFM	VGRGYASPD	SKLYKNCPKA	MKRLVADCVK	KVKEERPLFP	600
600	QILSSIELQ	HSLPKINRSA	SEPSLHRAAH	TEDINACTLT	TSPRLPVF		660

blue: RAF1 sequence expressed in recombinant protein **Red**: variant in recombinant protein

¹[NCBI/Protein](https://www.ncbi.nlm.nih.gov/Protein) accession number NP_002871.1