

ProQinase™ ARAF YDYD

A-Raf proto-oncogene, serine/threonine kinase

Recombinant Human Active Protein Kinase

HGNC Symbol: ARAF

Synonyms: ARAF1

Product No.: 1768-0000-1

Lot: 004

Description: Human ARAF, C-terminal fragment, amino acids K₂₈₂-P₆₀₆ (as in [NCBI/Protein](#) entry NP_001645.1), point mutations Y₃₀₁D/Y₃₀₂D, N-terminal GST-HIS₆ fusion protein with a 3C cleavage site, expressed in Sf9 insect cells

Product identity: ARAF YDYD LOT004, has been verified by mass spectrometry LC-ESI-MS/MS

Theoretical MW_{Fusion Protein}: 65578 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 10 µl are not advisable. Avoid repeated freeze-thaw cycles!

Protein concentration: 0.09 µg/µl

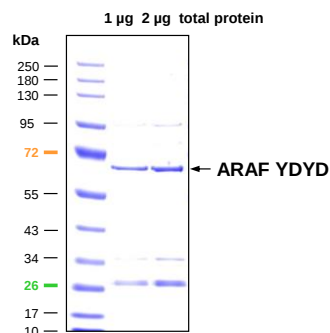
(Bradford method using BSA [Sigma, cat# A-7638, Lot 79H7641] as standard protein)

Biochemical Parameters:

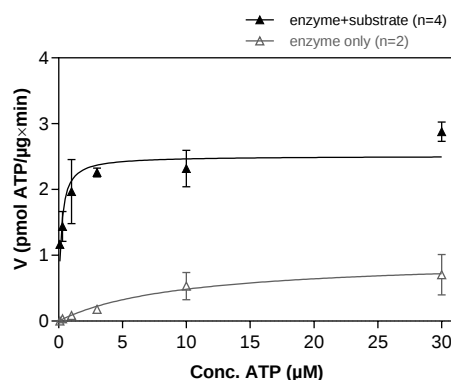
Specific kinase activity (P_i transfer): 2.5 pmol/µg*min

ATP-K_M: 0.2 µM

ARAF YDYD LOT004 Coomassie stain



ARAF YDYD LOT004 Determination of V_{max} and K_M value for ATP



- 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: 4 µg/ml

Assay technology:
Radiometric filter binding assay
MSFC membrane (96 well plate, Millipore)

Sequence information

GST-ARAF YDYD Recombinant Fusion Protein Amino Acid Sequence

1	MSPILGYWKI	KGLVQPTRLL	LEYLEEKYEE	HLIERDEGDK	WRNKKFELGL	EFPNLPYYID	60
61	GDVKLTQSMA	IIRYIADKHN	MLGGCPKERA	EISMLEGAVL	DIRYGVSRIA	YSKDFETLKV	120
121	DFLSKLPEML	KMFKDRLCHK	TYLNGDHVTH	PDFMLYDALD	VVLYMDPMCL	DAFPKLVCFK	180
181	KRIEAIPIQID	KYLKSSKYIA	WPLQGWQATF	GGGDHPPKSD	PMGHHHHHHG	RDSLEVLFGG	240
241	PLAMGARGRK	SLADDDKKVK	NLGYRDSGDD	WEVPPSEVQL	LKRIGTGSFG	TVFRGRWHGD	300
301	VAVKVLKVSQ	PTAEQAQAFK	NEMQVLRKTR	HVNILLFMGF	MTRPGFAIIT	QWCEGSSLYH	360
361	HLHVADTRFD	MVQLIDVARQ	TAQGMDYLHA	KNIIHRDLKS	NNIFLHEGLT	VKIGDFGLAT	420
421	VKTRWGAQAP	LEQPSGSVLW	MAAEVIRMQD	PNPYSFQSDV	YAYGVVLYEL	MTGSLPYSHI	480
481	GCRDQIIFMV	GRGYLSPDLS	KISSNCPKAM	RRLSDCLKF	QREERPLFPQ	ILATIELLQR	540
541	SLPKIERSAS	EPSLHRTQAD	ELPACLLSAA	RLVP			600

1-218: GST Red: HIS6-tag Green: 3C cleavage site blue: ARAF fragment boxed: Y₃₀₁DY₃₀₂D double mutation

ARAF wt¹ Amino Acid Sequence

1	MEPPRGPPAN	GAEPSRAVGT	VKVYLPNKQR	TVVTVRDGMS	VYDSLDKALK	VRGLNQDCCV	60
61	VYRLIKGRKT	VTAWDTAIAP	LDGEELIVEV	LEDVPLTMHN	FVRKTFFSLA	FCDFCLKFLF	120
121	HGFRCQTCGY	KFHQHCSSKV	PTVCVDMSTN	RQQFYHSVQD	LSGGSQRQHEA	PSNRPLNELL	180
181	TPQGSPSRTQ	HCDPEHFPPF	APANAPLQRI	RSTSTPNVHM	VSTTAPMDSN	LIQLTGQSFS	240
241	TDAAGSRGGS	DGTPRGSPSP	ASVSSGRKSP	HSKSPAEQRE	RKSLADDDKKK	VKNLGYRDSG	300
301	YYWEVPPSEV	QLLKRIGHTS	FGTVFRGRWH	GDVAVKVLKV	SQPTAEQAQA	FKNEMQVLRK	360
361	TRHVNILLFM	GFMTRPGFAI	ITQWCEGSSL	YHHLHVADTR	FDMVQLIDVA	RQTAQGMDYL	420
421	HAKNIIHRDL	KSNNIFLHEG	LTVKIGDFGL	ATVKTRWSGA	QPLEQPSGSV	LWMAAEVIRM	480
481	QDPNPYSFQS	DVYAYGVVLY	ELMTGSLPYS	HIGCRDQIIF	MVGRGYLSPD	LSKISSNCPK	540
541	AMRRLSDCL	KFQREERPLF	PQILATIELL	QRLPKIERS	ASEPSLHRTQ	ADELPACLLS	600
601	AARLVP						660

blue: ARAF sequence expressed in recombinant protein Red: variant in recombinant protein

¹[NCBI/Protein](https://www.ncbi.nlm.nih.gov/protein/1001645.1) accession number NP_001645.1