

ProQinase™ NIK

mitogen-activated protein kinase kinase kinase 14

Recombinant Human Active Protein Kinase

HGNC Symbol: MAP3K14

Synonyms: FTDCR1B, HS, HSNIK

Product No.: 0344-0000-1

Lot: 041

Description: Human NIK, full length, amino acids M₁-P₉₄₇ (as in [NCBI/Protein](#) entry NP_003945.2), N-terminal GST-HIS₆ fusion protein with a Thrombin cleavage site, expressed in Sf9 insect cells

Product identity: NIK Lot 041, was confirmed as NIK by mass spectroscopy LC-ESI-MS/MS

Theoretical MW_{Fusion Protein}: 133,437 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

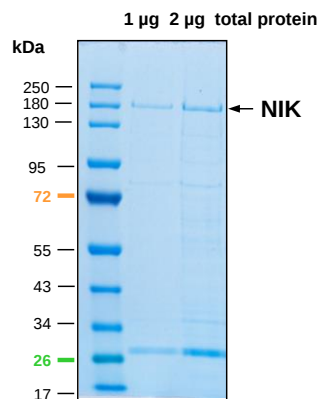
For complete recovery, mix well and spin before use. Product must not be stored in diluted solutions, aliquots below 10 µl are not advisable. Avoid repeated freeze-thaw cycles!

Protein concentration: 0.116 µg/µl
(Bradford method using BSA [Sigma, cat# A-7638, Lot 79H7641] as standard protein)

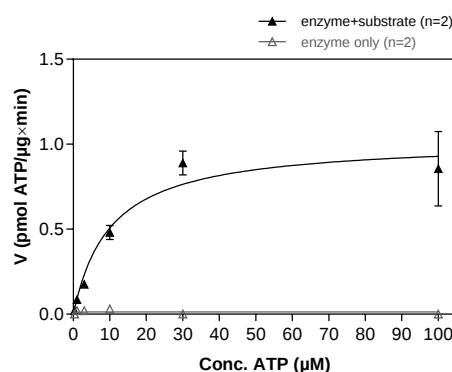
Biochemical Parameters:

Specific kinase activity (P_i transfer): 1 pmol/µg x min
ATP-K_M: 10 µM

NIK Lot 041:
Coomassie stain



NIK Lot 041:
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: CDC25C-derived peptide 160 µg/ml
 - Kinase: 2 µg/ml
- Filter binding assay
MSPH membrane (Millipore)

Sequence information

GST-NIK Recombinant Fusion Protein Amino Acid Sequence							
1	MSPILGYWKI	KGLVQPTRLL	LEYLEEKYEE	HLIERDEGDK	WRNKKFELGL	EFPNLPYYID	60
61	GDVKLTQSMA	IIRYIADKHN	MLGGCPKERA	EISMLEGAVL	DIRYGVSRIA	YSKDFETLKV	120
121	DFLSKLPEML	KMFEDRLCHK	TYLNGDHVTH	PDFMLYDALD	VVLYMDPMCL	DAFPKLVCFK	180
181	KRIEAIPIQID	KYLKSSKYIA	WPLQGWQATF	GGGDHPPKSD	PMGHHHHHHG	RRRASVAAGI	240
241	LVPKPGSPGLD	GICSRMAVME	MACPGAPGSA	VGQQKELPKA	KEKTPPLGKK	QSSVYKLEAV	300
301	EKSPVFCGKW	EILNDVITKG	TAKEGSEAGP	AAISIIAQAE	CENSQEFSP	FSEIFIFIAGS	360
361	KQYSQSESLD	QIPNNVAHAT	EGKMARVCWK	GKRRSKARKK	RKKKSSKSIA	HAGVALAKPL	420
421	PRTEQESCT	IPVQEDSPL	GAPYVRNTPQ	FTKPLKEPGL	GQLCFKQLGE	GLRPALPRSE	480
481	LHKLISPLQC	LNHVWKLHHP	QDGGPLPLPT	HPFPYSRLPH	PFPFHPLQPW	KPHPLESFLG	540
541	KLACVDSQKP	LPDPHLSKLA	CVDSPKPLPG	PHLEPSCLSR	GAHEKFSVEE	YLVHALQGSV	600
601	SSGQAHSLTS	LAKTWAARGS	RSREPSPKTE	DNEGVLLETK	LKPVDYREYRE	EVHWATHQLR	660
661	LGRGSFGEVH	RMEDKQGTGFQ	CAVKKVRLEV	FRAEELMACA	GLTSPRIVPL	YGAVREGPWV	720
721	NIFMELLEGG	SLGQLVKEQG	CLPEDRALYY	LGOALEGLEY	LHSRRILHGD	VKADNVLLSS	780
781	DGSHALCDF	GHAVCLQPDG	LKSLLLTGDI	IPGTETHMAP	EVVLGRSCDA	KVDVWSSCCM	840
841	MLHMLNGCHP	WTQFFRGPLC	LKIASEPPPV	REIPPSCAPL	TAQAIQEGLE	KEPIHRVSAA	900
901	ELGGKVNRL	QQVGGLKSPW	RGEYKEPRHP	PPNQANYHOT	LHAQPRELSP	RAPGPRPAEE	960
961	TTGRAPKLQP	PLPPEPPEPN	KSPPLTSLKE	ESGMWEPLPL	SSLEPAPARN	PSSPERKATV	1020
1021	PEQELQQLEI	ELFLNSLSQP	FSLEEQEQL	SCLSIDSLSL	SDDSEKNPSK	ASQSSRDTLS	1080
1081	SGVHSWSSQA	EARSSSWNMV	LARGRPTDTP	SYFNGVKVQI	QSLNGEHLHI	REFHRVKVGD	1140
1141	IATGISSQIP	AAAFSLVTKD	GQPVRYDMEV	PDSGIDLQCT	LAPDGSFAWS	WRVKHGQLEN	1200
1201	RP						1260

1-218: GST Red: HIS6-tag Pink: Thrombin cleavage site blue: NIK

NIK wt ¹ Amino Acid Sequence							
1	MAVMEMACPG	APGSAGVQQK	ELPKAKEKTP	PLGKKQSSVY	KLEAVEKSPV	FCGKWEILND	60
61	VITKGTAKG	SEAGPAAISI	IAQAECENSQ	EFSPTFSERI	FIAGSKQYSQ	SESLDQIPNN	120
121	VAHATEGKMA	RVCWKGKRRS	KARKKRKKKS	SKSLAHAGVA	LAKPLPRTPE	QESCTIPVQE	180
181	DESPLGAPYV	RNTPQFTKPL	KEPGLGQLCF	KQLGEGLRPA	LPRSELHCLI	SPLQCLNHVW	240
241	KLHHPQDGGP	LPLPPTHFPY	SRLPHFPFPH	PLQPWKPHPL	ESFLGKLACV	DSQKPLPDPH	300
301	LSKLACVDS	KPLPGPHLEP	SCLSRGAHEK	FSVEEYLVA	LQGSVSSGQA	HSLTSLAKTW	360
361	AARGSRREP	SPKTEDNEGV	LLTEKLKPD	YEYREEVHWA	THQLRLGRGS	FGEVHRMEDK	420
421	QTGFQCAVKK	VRLEVFRAGE	LMACAGLTSP	RIVPLYGAVR	EGPWVNIEME	LLEGSLGQL	480
481	VKEQGCLPED	RALYYLGQAL	EGLEYLHRR	ILHGDVKADN	VLLSSDGS	ALCDFGHAVC	540
541	LQPDGLGKSL	LTGDYIPGTE	THMAPEVLG	RSCDAKVDVW	SSCCMMLHML	NGCHPWTQFF	600
601	RGPLCLKIAS	EPPPVRREIP	SCAPLTAQAI	QEGLRKEPIH	RVSAAELGK	VNRLQQVVG	660
661	LKSPWRGEYK	EPRHPPPNQA	NYHOTLHAQP	RELSRAPAGP	RPAEETTGRA	PKLQPPLPPE	720
721	PPEPNKSPPL	TLSKEESGMW	EPLPLSSLEP	APARNPSSPE	RKATVPEQEL	QQLEIELFLN	780
781	SLSQPFSL	EQILSCLSI	DSLSDSDSE	KNPSKASQSS	RTLSSGVHS	WSSQAEARSS	840
841	SWNMVLARGR	PTDTPSYFNG	VKVQIQSLNG	EHLHIREFHR	VKVGDIATGI	SSQIPAAAFS	900
901	LVTKDGPVR	YDMEVPDGI	DLQCTLAPDG	SFAWSWRVKH	GQLENRP		960

blue: NIK sequence expressed in recombinant protein

¹NCBI/Protein accession number NP_003945.2