

## ProQinase™ NEK3

NIMA (never in mitosis gene a)-related kinase 3

Recombinant Human Active Protein Kinase

HGNC Symbol: NEK3

Synonyms: HSPK36, HSPK 36, MGC29949

Product No.: 1624-0000-1

Lot: 003

**Description:** Human NEK3, full length, amino acids M<sub>1</sub>-R<sub>506</sub> (as in NCBI/Protein entry NP\_689933.1), N-terminal GST-HIS<sub>6</sub> fusion protein with a 3C cleavage site, expressed in Sf9 insect cells

**Product identity:** NEK3 Lot 003, was confirmed as NEK3 by mass spectroscopy LC-ESI-MS/MS

**Theoretical MW<sub>Fusion Protein</sub>:** 85,954 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

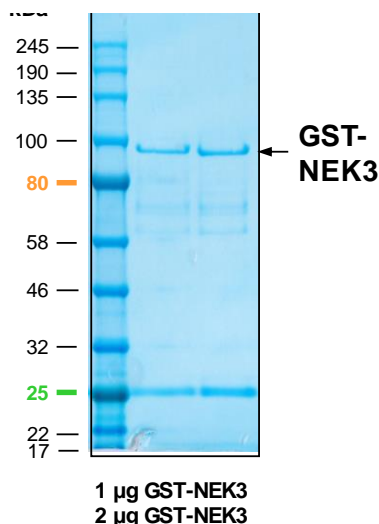
For complete recovery, mix well and spin before use. Product must not be stored in diluted solutions, aliquots below 10µl are not advisable. Avoid repeated freeze-thaw cycles!

**Protein concentration:** 0.263 µg/µl  
(Bradford method using BSA [Sigma, cat# A-7638, Lot 79H7641] as standard protein)

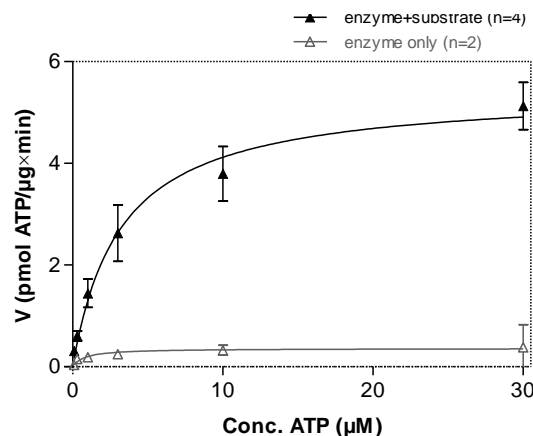
### Biochemical Parameters:

Specific kinase activity (P<sub>i</sub> transfer): 5.4 pmol/µg × min  
ATP-K<sub>M</sub>: 3.2 µM

NEK3 Lot 003:  
Coomassie stain



NEK3 Lot 003:  
Determination of V<sub>max</sub> and K<sub>M</sub> value for ATP



### Determination of K<sub>M</sub> value & Specific activity:

- Assay conditions:
  - 60 mM HEPES-NaOH, pH 7.5
  - 3 mM MgCl<sub>2</sub>
  - 3 mM MnCl<sub>2</sub>
  - 3 µM Na-orthovanadate
  - 1.2 mM DTT
  - 50 µg/ml PEG<sub>20,000</sub>
  - ATP (variable)
  - Substrate: CDC25C-derived peptide 80 µg/ml
  - Kinase: 2 µg/ml
- Filter binding assay
  - MSPH membrane (Millipore)

### Additional assay technology: NEK3 Lot 003

was also successfully tested by Reaction Biology for the use with the ADP-Glo™ Kinase assay from ADP-Glo assay conditions may vary from radiometric assay conditions, please inquire for assay details



## ProQinase™ NEK3

Product No.: 1624-0000-1

Recombinant Proteins

GST-NEK3 Recombinant Fusion Protein Amino Acid Sequence							
1	MSPILGYWKI	KGLVQPTRL	LEYLEEKYEE	HLYERDEGDK	WRNKKFELGL	EFPNLPYYID	60
61	GDVKLTQSMA	IIRYIADKHN	MLGGCPKERA	EISMLEGAVL	DIRYGVSRIA	YSKDFETLKV	120
121	DFLSKLP EML	KMFKDRLCHK	TYLNGDHVTH	PDFMLYDALD	VVLYMDPMCL	DAFPKLVCFK	180
181	KRIEAI PQID	KYLKSSKYIA	WPLQGWQATF	GGGDHPPKSD	PMGHHHHHG	RDSLEVL FQG	240
241	PLAMDDY MVL	RMIGEGSFGR	ALLVQL ESN	QMFAMKEIRL	PKSFSNTQNS	RKEAVLLAKM	300
301	KHPNIVAFKE	SFEAEGHLYI	VMEYCDGGDL	MQKIKQQK GK	LPEDMILNW	FTQMCLGVNH	360
361	IHKRVLHRD	IKSKNIFLTQ	NGKVKLGDFG	SARLLSNPMA	FACTYVGTPY	YVPPEIWENL	420
421	PYNNKSDIWS	LGCILYELCT	LKHPFQANSW	KNLILKVCQG	CISPLPSHYS	YELQFLVKQM	480
481	FKRNP SHRPS	ATLLSRGIV	ARLVQKCLPP	EIIMEYGEEV	LEEIKNSKHN	TPRKKNPSR	540
541	IRIALGNEAS	TVQEEEQDRK	GSHTDLESIN	ENLVESALRR	VNREEKGNKS	VHLRKASSPN	600
601	LHRRQWEK NV	PNTALTAL EN	ASILTSSLTA	EDDRGGSVIK	YSKNTTRKQW	LKETPD TLLN	660
661	ILKNADLSLA	FQTYTIYRPG	SEGFLKGPLS	EETEASDSVD	GGHDSVILDP	ERLEPGLDEE	720
721	DTDFEEEDDN	PDVSELKKR	AGWQGLCDR				780

1-218: GST   **Red**: HIS6-tag   **Green**: 3C cleavage site   **blue**:NEK3   **boxed**:variation from RefSeq

NEK3 wt <sup>1</sup> Amino Acid Sequence							
1	MDDYMLRMI	GE GSFGRALL	VQH ESNQMF	AMKEIRLPKS	FSNTQNSRKE	AVLLAKMKHP	60
61	NIVAFKESFE	AEGHLYIVME	YCDGGDLMQK	IKQQKGLF P	EDMILNWFTQ	MCLGVNHIHK	120
121	KRVLHRDIKS	KNIFLTQNGK	VKLGDFGSAR	LLSNPMAFAC	TYVGT PYYVP	PEIWENLPYN	180
181	NKSDIWSLGC	ILYELCTLKH	PFQANSWKNL	ILKVCQGCIS	PLPSHYSYEL	QFLVKQMFKR	240
241	NPSHRPSATT	LLSRGIVARL	VQKCLPPEII	MEYGEEVLEE	IKNSKHNTPR	KKTNP SRIRI	300
301	ALGNEASTVQ	EEEQDRK GSH	TDLESINENL	VESALRRVNR	EEKGNKSVHL	RKASSPNLHR	360
361	RQWEKNVPNT	ALTAL ENASI	LTSSLTAEDD	RGGSVIKYSK	NTTRKQWLKE	TPD TLLN ILK	420
421	NADLSLAFQT	YTIYRPGSEG	FLKGPLSEET	EASDSVDGGH	DSVILDPERL	EPGLDEEDTD	480
481	FEEEDDNPDW	VSELKKRAGW	QGLCDR				540

**blue**: NEK3 sequence expressed in recombinant protein   **Red**: variant in recombinant protein

<sup>1</sup>NCBI/Protein accession number NP\_689933.1  
H23L: SNP variation see NCBI/dbSNP ID: rs17482764