

ProQinase™ NEK8

NIMA related kinase 8

Recombinant Human Active Protein Kinase

HGNC Symbol: NEK8

Synonyms: NPHP9

Product No.: 2015-0000-1

Lot: 001

Description: Human NEK8, N-terminal fragment, amino acids M₁-L₂₆₅ (as in [NCBI/Protein](#) entry NP_835464.1), N-terminal GST-HIS₆ fusion protein with a 3C cleavage site, expressed in Sf9 insect cells

Product identity: NEK8 Lot 001, has been verified by mass spectrometry LC-ESI-MS/MS

Theoretical MW_{Fusion Protein}: 58198 Da

Expression host: Sf9 insect cells

Purification: GST-Affinity Chromatography

Activation: in vitro auto activation

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: 0.57 µg/µl

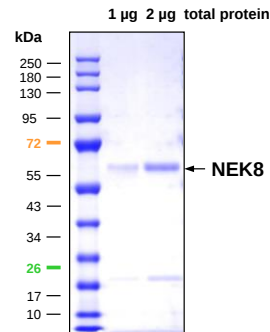
(Bradford method using BSA [Sigma, cat# A-7638, Lot 79H7641] as standard protein)

Biochemical Parameters:

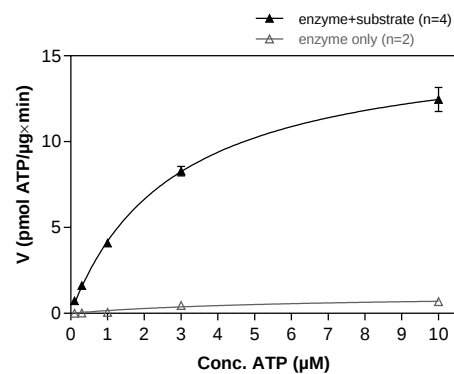
Specific kinase activity (P_i transfer): 16 pmol/µg*min

ATP-K_M: 2.8 µM

NEK8 Lot 001:
Coomassie stain



NEK8 Lot 001:
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: Casein 40 µg/ml
- Kinase: 4 µg/ml

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

Recombinant Proteins

Sequence information

GST-NEK8 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	VVLYMDPMC	DAFPKLVCFK	180
181	KRIEAIPIQID	KYLKSSKYIA	WPLQGWQATF	GGGDHPPKSD	PMGHHHHHHG	RDSLEVLFGG	240
241	MEKYERIRV	VGRGAFGIVH	LCLRKADQKL	VIIKQIPVEQ	MTKEERQAAQ	NECQVLKLLN	300
301	HPNVIEWEYEN	FLEDKALMIA	MEYAPGGTLL	EFIQKRCNSL	LEEETILHFF	VQILLALHHV	360
361	HTHLILHRDL	KTQNILLDKH	RMVVKIGDFG	ISKILSSKSK	AYTVVGTPCY	ISPCEGCKP	420
421	YNQKSDIWAL	GCVLVELASL	KRAFEAANLP	ALVLKIMSGT	FAPISDRYSP	ELRQLVLSLL	480
481	SLEPAQRPL	SHIMAQPLCI	RALLNL				540

1-218: GST Red: HIS6-tag Green: 3C cleavage site blue: NEK8 fragment

NEK8 wt ¹ Amino Acid Sequence							
1	MEKYERIRV	GRGAFGIVHL	CLRKADQKL	VIIKQIPVEQ	MTKEERQAAQ	NECQVLKLLN	60
61	PNVIEWEYEN	FLEDKALMIA	MEYAPGGTLL	EFIQKRCNSL	LEEETILHFF	VQILLALHHV	120
121	THLILHRDL	KTQNILLDKH	RMVVKIGDFG	ISKILSSKSK	AYTVVGTPCY	ISPCEGCKP	180
181	NQKSDIWAL	GCVLVELASL	KRAFEAANLP	ALVLKIMSGT	FAPISDRYSP	ELRQLVLSLL	240
241	LEPAQRPL	SHIMAQPLCI	RALLNL				300
301	GPVPAIPPP	LSSVYAWGGG	LGTPLRLPML	NTEVVQVAAG	RTQKAGVTRS	GRLILWEAPP	360
361	LGAGGGSLLP	GAVEQPQPQF	ISRFLEGQSG	VTIKHVACGD	FFTACLTDG	IIMTFGSGSN	420
421	GCLGHGSLTD	ISQPTIVEAL	LGYEMVQVAC	GASHVLALST	ERELFAWGRG	DSGRLGLGTR	480
481	ESHSCPQQVP	MPPGQEAQRV	VCGIDSSMIL	TVPGQALACG	SNRFNKLGLD	HLSLGEEPVP	540
541	HQQVEEALSF	TLLGSAPLDQ	EPLLSIDLGT	AHSAAVTASG	DCYTFGSNQH	GQLGTNTRRG	600
601	SRAPCKVQGL	EGIKMAMVAC	GDAFTVAIGA	ESEVYSWGKG	ARGRLGRRDE	DAGLPRPVQL	660
661	DETHPYTVTS	VSCCHGNTLL	AVRSVTDEPV	PP			720

blue: NEK8 sequence expressed in recombinant protein

¹[NCBI/Protein](https://www.ncbi.nlm.nih.gov/Protein/835464.1) accession number NP_835464.1