

Data Sheet

NEK9

Human, recombinant, N-terminal GST tag

Catalog #: 40142

Lot #: 111129-5 **Conc.:** 0.19 mg/ml

Formulated in: 40 mM Tris-HCl, pH 8, 110 mM NaCl, 2.2 mM KCl, 4 mM Glutathione, 20% glycerol, and 3 mM DTT.

Stability: >12 months at -80°C

References:

1. Roig, J., *et al. Genes Dev.* **16**: 1640-1658 (2002).
2. Belham, C., *et al. Curr. Biol.* **15**: 1155-1167 (2001).

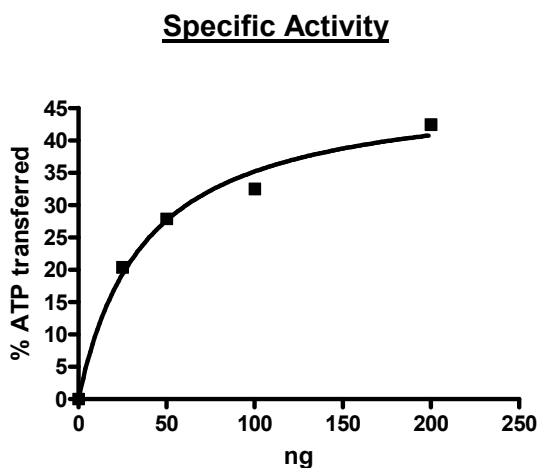
Description: Human NEK9, also known as NERCC (GenBank Accession No. NM_033116), a.a 1-979 (end), truncation of a.a. 346-731, with N-terminal GST tag, MW =92 kDa, expressed in Sf9 insect cells via a baculovirus expression system.

Specific Activity: 147 pmol/min/μg

Assay Buffer: Enzyme reaction is conducted in a Kinase Buffer (50 mM Tris pH 7.4, 20 mM MgCl₂, 0.1 mg/ml BSA and 1 mM DTT) containing 20 μM ATP and 0.2 mg/ml MBP for 30 minutes at 30°C. Enzyme activity is measured using Kinase-Glo plus reagents (Promega).

Application: Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

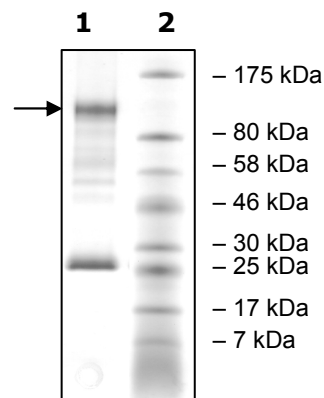
Quality Assurance



10% SDS-PAGE **Coomassie staining**

Lane 1:
1.1 μg NEK9
Lane 2:
Protein Marker
BioLabs (#P7708L)

MW: 92 kDa
Purity: ≥38%



OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.829.3082** Fax **1.858.481.8694**

Or you can Email us at: info@bpsbioscience.com

Please visit our website at: www.bpsbioscience.com



6044 Cornerstone Court W, Ste E
San Diego, CA 92121
Tel: 1.858.829.3082
Fax: 1.858.481.8694
Email: info@bpsbioscience.com

Amino Acid Sequence:

MSPILGYWKIKGLVQPTRLLLEYLEEKYEEHLYERDEGDKWRNKKFELGLEFPNLPYYID
GDVCLTQSMARIYIADKHNMLGGCPKERAEISMLEGAVLDIRYGVSRIAYSKDFETLKVDF
LSKLPPEMLKMFEDRLCHKTYLNGDHVTHPDFMLYDALDVVLYMDPMCLDAFPKLVCFKK
RIEAIPQIDKYLKSSKYIAWPLQGQWQATFGGGDHPPKSDPLVPRGSPGFMMSVLGEYER
HCDSINSDFGSESGGCGDSSPGPSASQGPRAGGGAAEQEELHYIPIRVLGRGAFGEAT
LYRRTEDDSLWVWKEVDLTRLSEKERRDALNEIVILALLQHDNIIAYYNHFMDNTLLIEL
EYCNGGNLYDKILRQKDKLFEEMVVWYLFQIVSAVSCIHKAGILHRDIKTLNIFLTKANL
IKLGDYGLAKKLNSEYSMAETLVGTPYYMSPELCQGVKYNFKSDIWAIVGCVIFELLTLK
RTFDATNPLNLCVKIVQGIRAMEVDSSQYSLELIQMVHSCLDQDPEQRPTADELLDRPL
LRKRRREMEEKVTLNAPTKRPRSSTVTEAPIAVVTSRRTKTIRSNSSGLSIGTVFQSSSP
GGGGGGGGGEEEDSQESETPDPSGGFRGTMEADRGMEGLISPTTEAMGNSNGASSS
CPGWLKELNAEFIPMPDPSPLSAFSESEKDTLPYEELQGLKVASEAPLEHKPQV
EASSPRLNPAVTCAGKGTPLTPACACSSLQVEVERLQGLVLKCLAEQQKLQENLQI
FTQLQKLNKKLEGGQQVGMHSGTQTAKKEEMEMDPKPDLDSDSWCLLGTDSRPSL

OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.829.3082** Fax **1.858.481.8694**
Or you can Email us at: info@bpsbioscience.com
Please visit our website at: www.bpsbioscience.com