

Data Sheet

NEK7, GST-Tag

Human, recombinant, N-terminal GST tag

Catalog #: 40141

Lot #: 130319-1 **Conc.:** 0.22 mg/ml

Formulated in: 40 mM Tris-HCl, pH 8.0,
 110 mM NaCl, 2.2 mM KCl, 16 mM
 Glutathione, 3 mM DTT and 20% glycerol.

Stability: At least 12 months at -80°C

References:

1. Kandli, M., *et al. Genomics* **68**: 187-196 (2000).
2. Belham, C., *et al. Curr. Biol.* **15**: 1155-1167 (2001).

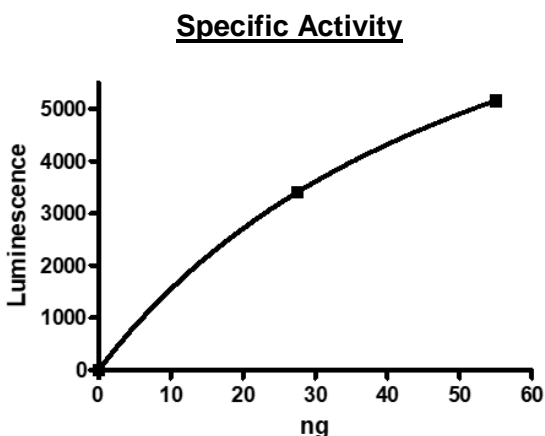
Description: Human NEK7, also known as NIMA Related Kinase 7 (GenBank Accession No. NM_133494, a.a. 1-302 (end)), full length, with N-terminal GST tag, MW= 61 kDa, expressed in Sf9 insect cells via a baculovirus expression system.

Specific Activity: 47 pmol/min/μg

Assay conditions: Assay was done in a reaction buffer (25 μl) containing 40 mM Tris, pH 7.4, 20 mM MgCl₂, 0.1 mg/ml BSA, 100 μM ATP, 0.2 mg/ml Casein and NEK7 for 50 min at 30°C. NEK7 activity was measured using ADP-Glo reagents (Promega).

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

Quality Assurance

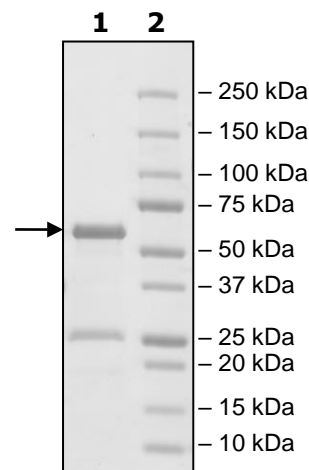


**4-20% SDS-PAGE
Coomassie staining**

Lane 1:
2.98 μg NEK7

Lane 2:
Protein Marker

MW: 61 kDa
Purity: ≥72%



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