

## Data Sheet

### NEK7, His-Tag

Human, recombinant, N-terminal His-tag

**Catalog #:** 100476

**Lot #:** 190916-1      **Conc.:** 1.70 mg/ml

**Description:** Human NEK7, also known as NIMA Related Kinase 7, GenBank Accession No. NM\_133494, a.a. 2-302 (end), with N-terminal His-tag, expressed in a HEK293 cell expression system. MW=37 kDa.

**Formulated in:** 20 mM Tris-HCl, pH 7.5, 200 mM NaCl, 0.5 mM TCEP, 10% glycerol, and 140 mM Imidazole.

**Stability:** At least 12 months at -80°C.  
Avoid freeze/thaw cycles.

#### References:

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

### Quality Assurance

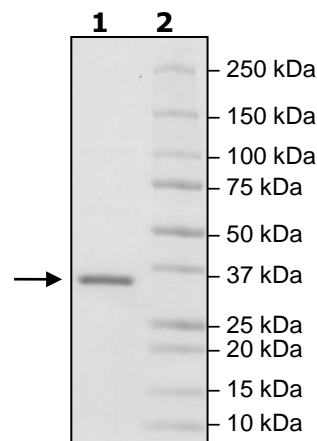
#### 4-20% SDS-PAGE Coomassie staining

**Lane 1:**  
NEK7

**Lane 2:**  
Protein Marker

**MW:** 37 kDa

**Purity:** ≥90%



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

To place your order, please contact us by Phone **1.858.202.1401** Fax **1.858.481.8694**

Or you can Email us at: [info@bpsbioscience.com](mailto:info@bpsbioscience.com)

Please visit our website at: [www.bpsbioscience.com](http://www.bpsbioscience.com)