

## Data Sheet

### TNK1, GST-Th-Tag

Human, Recombinant, N-terminal GST-Tag,  
 Thrombin Cleavage Site

**Catalog #:** 80019

**Lot #:** 170413-G2      **Conc.:** 0.08 mg/ml

**Formulated in:** 40 mM Tris-HCl, pH 8.0,  
 110 mM NaCl, 2.2 mM KCl, 0.04%  
 Tween20, 20% glycerol, 9mM Glutathione

**Stability:** At least 6 months at  $-80^{\circ}\text{C}$ . *Avoid freeze/thaw cycles. Storing diluted enzyme is not recommended, if necessary, use carrier protein (BSA 0.1 – 0.5%).*

**References:**

1. Felschow, D. (2000). *BBRC*. **273(1)**:294-301.
2. Azoitei, N. *et al.* (2010). *Oncogene*. **26**: 6536-6545.

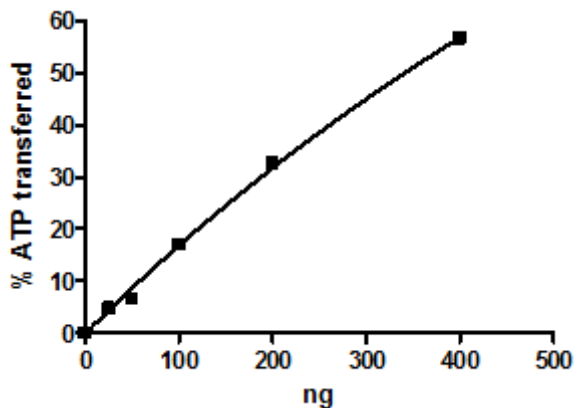
**Description:** TNK1 (Tyrosine Kinase Non Receptor 1), also known as CD38 Negative Kinase 1 (GenBank Accession No. NM\_003985), a.a. 106-390, with N-terminal GST-tag and Thrombin Cleavage Site, MW = 57 kDa, expressed in Sf9 cells via a baculovirus expression system.

**Specific Activity:**  $\geq 3.5$  pmol/min/ug  
 Assay Conditions: Assay was done in Kinase buffer containing 1 mM DTT using CSKtide substrate (0.2 mg/ml) and 10  $\mu\text{M}$  ATP. Reaction was done at  $30^{\circ}\text{C}$  for 50 min. Amount of ATP transferred was calculated using Kinase-Glo reagent (Promega)

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

## Quality Assurance

### Specific Activity

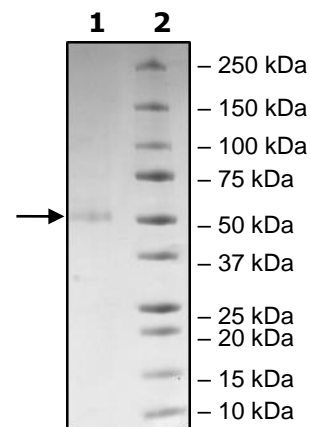


### 4-20% SDS-PAGE Coomassie staining

**Lane 1:**  
TNK1

**Lane 2:**  
Protein Marker

**MW:** 57 kDa  
**Purity:**  $\geq 90\%$



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

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



6042 Cornerstone Court W, Ste B  
San Diego, CA 92121  
**Tel:** 1.858.829.3082  
**Fax:** 1.858.481.8694  
**Email:** [info@bpsbioscience.com](mailto:info@bpsbioscience.com)

### Sequence

MSPILGYWKIKGLVQPTRLLEYLEEKYEEHLYERDEGDKWRNKKFELGLEFPNL  
PYYIDGDVKLTQSMAIIRYIADKHNMLGGCPKERAETSMLEGAVLDIRYGVSRIA  
YSKDFETLKVDLFLSKLPEMLKMFKDRLCHKTYLNGDHVTHPDFMLYDALDVVLYM  
DPMCLDAFPKLVCFKKRIEAI PQIDKYLKSSKYIAWPLQGWQATFGGGDHLVPRG  
SGLKCLIPGAVCRGELLGSGCFGVVHRGLWTLPSGKSVPVAVKSLRVGPEGPMG  
TELGDFLREVSVMNLEHHPVLRRLHGLVLGQPLQMVME LAPLGS LHARLTAPAPT  
PPLLVALLCFLRQLAGAMAYLGARGLVHRDLATRNLLLASPRTIKVADFGLVRP  
LGGARGRYVMGGPRPIPYAWCAPESLRHGAFSSASDVWMFGVTLWEMFSGGEEPW  
AGVPPYLILQRLEDRARLPRPPLCSRALYSLALRCWAPHPADRPSFSHLEGLLQE  
AGPSEACCVRD

**Red:** GST tag

**Blue:** Linker including Thrombin Cleavage Site

**Blue:** BamH1 site

**Black:** TNK1 (aa 106-390)

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