

## Data Sheet

### RIPK1, GST-Th-Tag

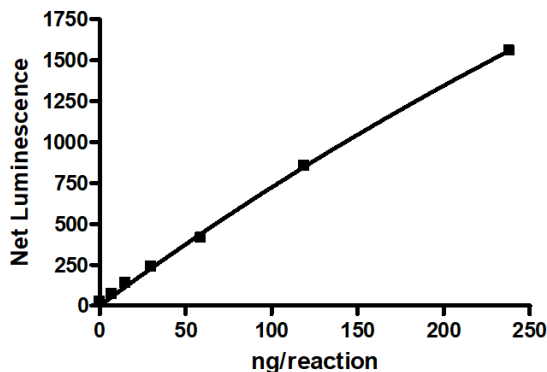
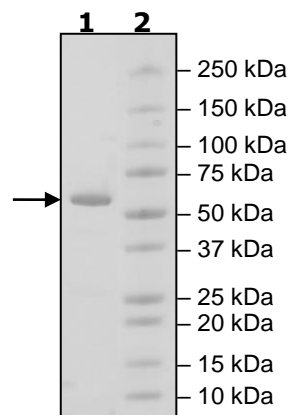
Human, Recombinant,  
N-terminal GST-Th-Tag**Catalog #:** 40371**Lot #:** 181106-G**Conc.:** 0.84 mg/ml**Formulated in:** 40 mM Tris-HCl, pH 8.0,  
110 mM NaCl, 2.2 mM KCl, and 20%  
Glycerol.**Stability:** At least 6 months at  $-80^{\circ}\text{C}$ .  
Avoid freeze/thaw cycles.**References:**

1. Hsu, H. *et al.*, *Immunity* **4**: 387-396, 1996.
2. Kim, J Y. *et al.*, *J Cell Sci.* 2011 Feb 15;**124**(Pt 4):647-56.

**Description:** Human RIPK1, also known as RCK and MAP3K19, GenBank Accession No. NM\_003804, a.a. 1-327, with N-terminal GST-tag and thrombin cleavage site, expressed in a Baculovirus-infected Sf9 cell expression system. MW=64 kDa.**Specific Activity:**  $\geq 3$  pmol/min/ $\mu\text{g}$ .Assay Conditions: RIPK1 reactions was done in 25  $\mu\text{l}$  containing 40 mM Tris, pH 7.4, 20 mM  $\text{MgCl}_2$ , 0.1 % BSA, 3 mM DTT, 0.2 mg/ml MBP and 20  $\mu\text{M}$  ATP at  $30^{\circ}\text{C}$  for 45 min. The amount of ATP transferred to the substrate was measured using ADP-Glo reagents (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:**  
2  $\mu\text{g}$  RIPK1  
**Lane 2:**  
Protein Marker**MW:** 64 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 **1.858.202.1401** 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)