

## Data Sheet

### RIPK4, GST-Tag

Human, Recombinant, N-terminal GST-Tag

**Catalog #:** 79712

**Lot #:** 190214

**Conc.:** 0.50 mg/ml

**Formulated in:** 50 mM Tris-HCl, pH 7.5, 150 mM NaCl, 10mM glutathione, 0.1 mM EDTA, 0.25 mM DTT, 0.1 mM PMSF, 25% glycerol.

**Stability:** At least 6 months at  $-80^{\circ}\text{C}$ .  
Avoid freeze/thaw cycles.

#### References:

1. Mitchell, K. et al., *Am. J. Hum. Genet.* **90**: 69-75, 2012.
2. Rountree, R.B. et al., *J. Invest. Dem.* **130**: 102-112, 2010

#### Description:

Human RIPK4, also known as ANKK2, DIK, PKK, RIP4, and ANKRD3, GenBank Accession No. BC110618, a.a. 15-426, with N-terminal GST-tag, expressed in a Baculovirus-infected Sf9 cell expression system. MW = 57 kDa.

**Application:** Useful for protein binding studies, as an immunogen, or for Western blotting

## Quality Assurance

### 4-20% SDS-PAGE

#### Coomassie staining

#### Lane 1:

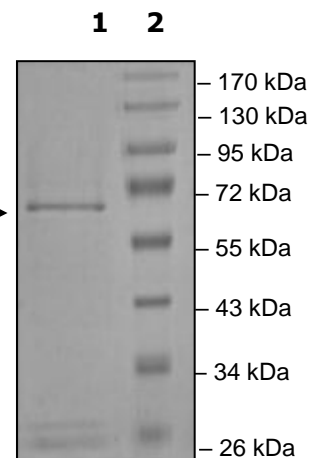
RIPK4 →

#### Lane 2:

Protein Marker

**MW:** 57 kDa

**Purity:**  $\geq 80\%$



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)