

Data Sheet

RIPK5, GST-Tag

Human, Recombinant, N-terminal GST-Tag

Catalog #: 79713

Lot #: 190214

Conc.: 1.0 mg/ml

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

Stability: At least 6 months at -80°C .
Avoid freeze/thaw cycles.

References:

1. Peng, J. et al., *Biochem. Biophys. Acta.* **1759(11-12)**: 562-72., 2006.
2. Zha, J. et al., *Biochem., Biophys. Res. Commun.* **319**: 298-303, 2004.

Description:

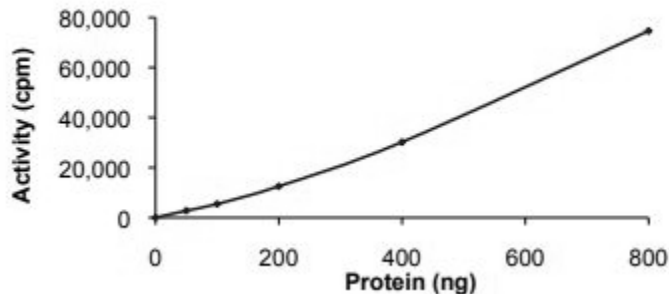
Human RIPK5, also known as DSTYK, DustyPK, HDCMD38P, KIAA0472, and RIP5, GenBank Accession No. NM_015375, a.a. full length, with N-terminal GST-tag, expressed in a Baculovirus-infected Sf9 cell expression system. MW = 140 kDa.

Specific Activity: 6 pmol/min/ μg .

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

Quality Assurance

Specific Activity

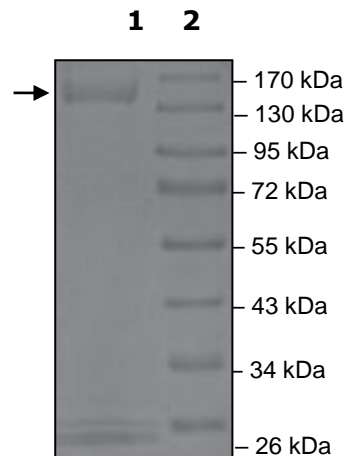


4-20% SDS-PAGE Coomassie staining

Lane 1:
RIPK5

Lane 2:
Protein Marker

MW: 140 kDa
Purity: $\geq 85\%$



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

Please visit our website at: www.bpsbioscience.com