

Data Sheet

RIPK3, GST-Tag

Human, Recombinant, N-terminal GST-Tag

Catalog #: 79710**Lot #:** 190213**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°C .
Avoid freeze/thaw cycles.

References:

1. Kasof, G. M. et al., *FEBS Lett.* **47**: 285-291, 2000.
2. Oberst, A. et al., *Nature* **471**: 3636-367. 2011.

Description:

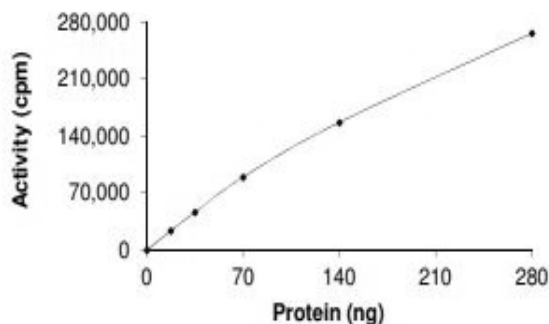
Human RIPK3, also known as Receptor Interacting Serine/Threonine Kinase 3 and RIP3, GenBank Accession No. NM_006871, a.a. full length, with N-terminal GST-tag, expressed in a Baculovirus-infected Sf9 cell expression system. MW = 96 kDa.

Specific Activity: ≥ 44 pmol/min/ μg .

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

Quality Assurance

Specific Activity



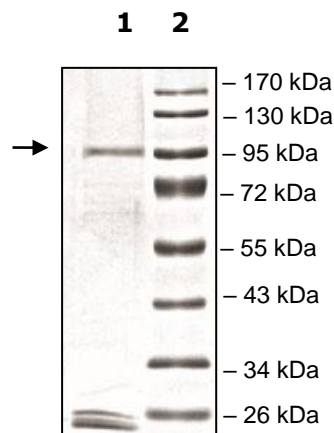
4-20% SDS-PAGE

Coomassie staining

Lane 1:
RIPK3

Lane 2:
Protein Marker

MW: 96 kDa
Purity: $\geq 70\%$



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