

Data Sheet

ROR2, Fc-Fusion (IgG1), Avi-Tag, Biotin-Labeled

Human, Recombinant, C-terminal Fc-Avi-tag,
Biotin-Labeled

Catalog #: 100046

Lot #: 180426 **Conc.:** 1.23 mg/ml

Description: Human ROR2, also known as tyrosine-protein kinase transmembrane receptor, neurotrophic tyrosine kinase, and receptor related 2, GenBank Accession No. NM_004560, a.a. 28-403, with C-terminal Avi-Tag fused to the Fc region of Human IgG1, and expressed in a Hek293 expression system. MW=71 kDa. This protein is enzymatically biotinylated using Avi-Tag™ technology.

Formulated in: 8 mM Phosphate, pH 7.4, 110 mM NaCl, 2.2 mM KCl, and 20% glycerol

Stability: At least 6 months at -80°C. Avoid freeze/thaw cycles. Storing diluted enzyme is not recommended, if necessary, use carrier protein (BSA 0.1 – 0.5%).

References:

1. DeChiara, T.M. *et al.*, *Nature Genet.* 2000; **24**: 271-274.
2. Nomi, M. *et al.*, *Molec. Cell Biol.* 2001; **21**: 8239-8335.

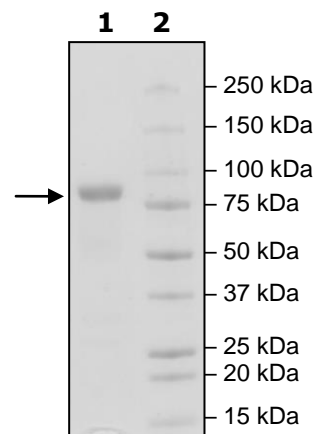
Quality Assurance

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

Lane 1:
4µg ROR2

Lane 2:
Protein Marker

MW:
71 kDa +glycans
Purity: ≥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**

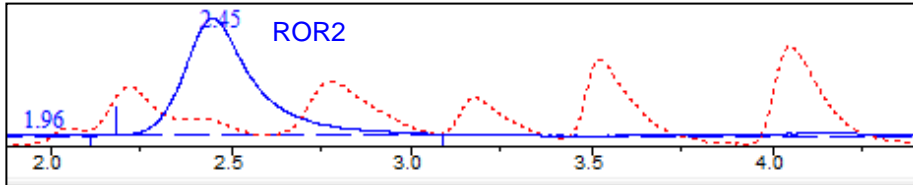
Or you can Email us at: info@bpsbioscience.com

Please visit our website at: www.bpsbioscience.com



6042 Cornerstone Court West, Suite B
San Diego, CA 92121
Tel: 1.858.829.3082
Fax: 1.858.481.8694
Email: info@bpsbioscience.com

Gel Filtration Curve



<10% aggregation

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
Please visit our website at: www.bpsbioscience.com