

Data Sheet

ROR1, Fc-Fusion (IgG1), Avi-Tag, Biotin-Labeled

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

Catalog #: 79482

Lot#: 180301

Conc.: 0.76 mg/ml

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.*

References:

1. Fukuda, T., *et al.*, *PNAS*. 2007 Dec;
105(8): 3047-3052.
2. Hudecek, M., *et al.*, *Clinical Cancer
Research*. 2013 Jun; **19(12)**: 3109.

Description:

Human ROR1, also known as receptor tyrosine kinase-like orphan receptor 1, neurotrophic tyrosine kinase, and receptor-related 1 (NTRKR1), GenBank Accession No. NM_005012, a.a 24-403, with C-terminal Avi-TagTM fused to the Fc portion of Human IgG1 and expressed in a HEK293 cell expression system. MW = 71 kDa. This protein runs at a higher apparent M.W. by SDS-PAGE due to glycosylation. This protein is enzymatically biotinylated using Avi-TagTM technology.

Quality Assurance

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

Lane 1:

4 μg ROR1

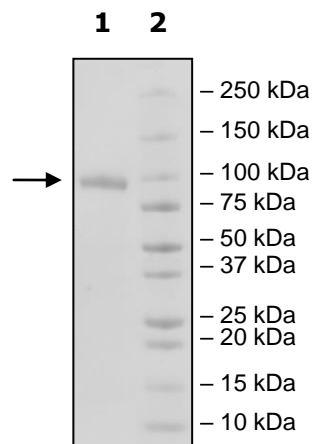
Lane 2:

Protein Marker

MW:

71kDa + glycans

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 **1.858.829.3082** Fax **1.858.481.8694**

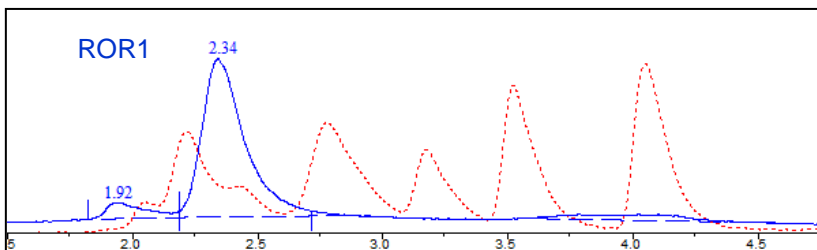
Or you can Email us at: info@bpsbioscience.com

Please visit our website at: www.bpsbioscience.com



6042 Cornerstone Court W, Ste 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