

Data Sheet

CD112, His-tag, Biotin-labeled

Human, recombinant, C-terminal His-tag,
biotin labeled

Catalog #: 71234

Lot #: 150625-1 **Conc.:** 0.58 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. Protein may be diluted to
 $\geq 100 \mu\text{g/ml}$ in PBS + glycerol and stored at
 -80°C .*

References:

1. Bottino, C., *et al.*, *J. Exp. Med.* 2003;
198(4): 557-567.
2. Pende, D., *et al.*, *Blood.* 2005; 105(5):
2066-2073.

Description:

Human secreted CD112, also known as Nectin-2, Herpes virus entry mediator B, HVEB, PVRL2, and PRR2, GenBank Accession No. NM_002856, a.a. 32-360 expressed in a HEK293 cell expression system, with C-terminal His and AviTM tags, this protein is enzymatically biotinylated using Avi-tagTM technology. MW = 38.3 kDa. This protein runs at a higher MW by SDS-PAGE due to glycosylation.

Application:

Useful for studying protein binding and for screening small molecules and antibodies.

Quality Assurance

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

Lane 1:

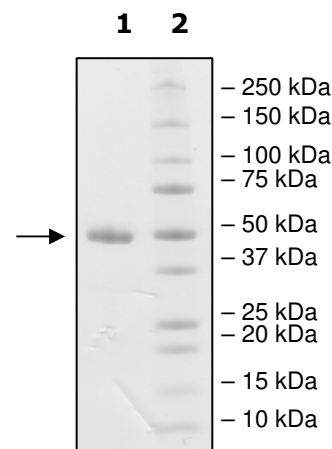
3 μg CD112

Lane 2:

Protein Markers

MW: 38.3 kDa

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**

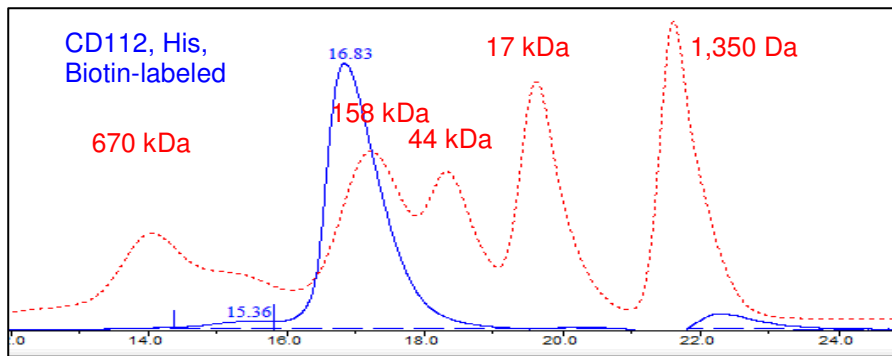
Or you can Email us at: info@bpsbioscience.com

Please visit our website at: www.bpsbioscience.com



6044 Cornerstone Court W, Ste E
San Diego, CA 92121
Tel: 1.858.829.3082
Fax: 1.858.481.8694
Email: info@bpsbioscience.com

Gel Filtration Curve



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**](mailto:info@bpsbioscience.com)
Please visit our website at: [**www.bpsbioscience.com**](http://www.bpsbioscience.com)