

Data Sheet

CD28H, Fc fusion, Biotin-labeled

Human, recombinant, C-terminal Fc fusion (Human IgG1), C-terminal AviTag™

Catalog#: 71239

Lot#: 150901-G **Conc.:** 0.31 mg/ml

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

Stability: At least 6 months at -80°C. Avoid freeze/thaw cycles. Protein may be diluted to ≥ 100 µg/ml in PBS + glycerol and stored at -80°C.

References:

1. Zhu, Y., *et al.*, *Nat. Commun.* 2013; **4**: 2043.
2. Zhu, Y., *et al.*, *J. Immunol.* 2015; **194**: 142.2.

Description:

Human CD28 homolog (CD28H), also known as Transmembrane and immunoglobulin domain-containing protein 2 (TMIGD2) and immunoglobulin and proline-rich receptor 1 (IGPR-1). GenBank Accession No. NM_144615, a.a. 23-128 fused at the C-terminus to the Fc portion of human IgG1 followed by a C-terminal AviTag™ expressed in a HEK293 expression system. MW = 40 kDa. This protein is enzymatically biotinylated using Avi-tag™ technology. Biotinylation is confirmed to be greater than 90%. 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.

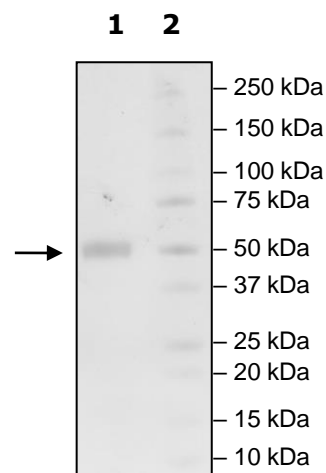
Quality Assurance

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

Lane 1:
 3 µg CD28H

Lane 2:
 Protein Marker

MW: 40 kDa+glycans
Purity: ≥90%



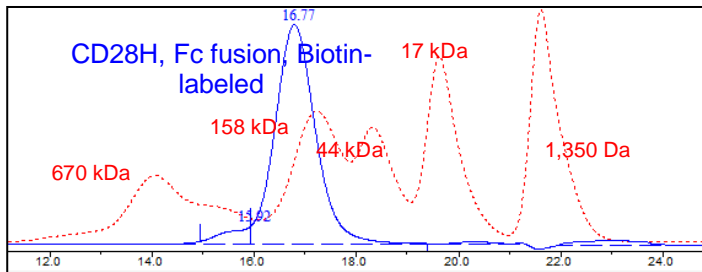
OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.202.1401** Fax **1.858.481.8694**
 Or you can Email us at: support@bpsbioscience.com
 Please visit our website at: www.bpsbioscience.com



6405 Mira Mesa Blvd Ste 100
San Diego, CA 92121
Tel: 1.858.202.1401
Fax: 1.858.481.8694
Email: support@bpsbioscience.com

Gel Filtration Curve



<5% aggregation

OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.202.1401** Fax **1.858.481.8694**
Or you can Email us at: support@bpsbioscience.com
Please visit our website at: www.bpsbioscience.com