

Data Sheet

OX40, Fc-Fusion, Avi-Tag, Biotin-Labeled (Mouse)

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

Catalog #: 79357

Lot #: 171130-G **Conc.:** 0.44 mg/ml

Formulated in: 8mM Phosphate, 110mM NaCl, 2.2mM KCl, pH7.4, 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. Linton, P.J., *et al.*, *J. Exp. Med.* 2003; **197(7)**: 875-883.
2. Arestides, R.S.S., *et al.*, *Eur. J. Immunol.* 2002; **32(10)**: 2874-2880.

Description: Mouse OX40, also known as Tumor Necrosis Factor Receptor Superfamily Member 4, TNFRSF4, and CD134, GenBank Accession No. NM_011659, a.a. 20-211 with C-terminal Avi-Tag fused to the Fc portion of Human IgG1 and expressed in a HEK293 cell expression system. This protein is enzymatically biotinylated using Avi-Tag™ technology. This protein runs at a higher MW by SDS-PAGE due to glycosylation. MW = 52 kDa.

Quality Assurance

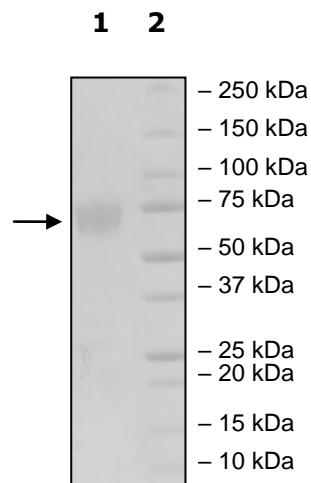
4-20% SDS-PAGE Coomassie staining

Lane 1:
5 μg CD40(Mouse)

Lane 2:
Protein Marker

MW:
52 kDa + 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**

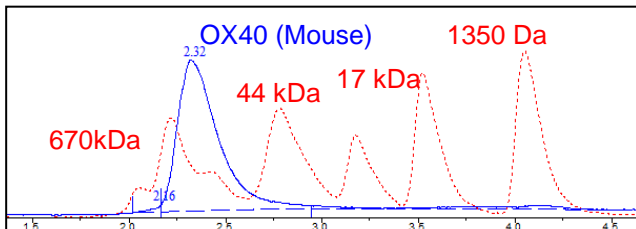
Or you can Email us at: info@bpsbioscience.com

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