

Data Sheet

OX40 (CD134), Biotin-labeled, His-tag

Human, recombinant, with C-terminal Avi-tag™ and His-tag

Catalog #: 71310

Lot #: 150806-A **Conc.:** 0.44 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. Rogers, P.R. *et al.*, *Immunity*. 2001; **15(3)**: 445-455.
2. Ito, T. *et al.*, *J. Exp. Med.* 2005; **202(9)**: 1213-1223.

Description:

Human secreted OX40, also known as CD134, Tumor Necrosis Factor Receptor Superfamily Member 4, and TNFRSF4, GenBank Accession No. NM_003327, a.a. 29-216 with C-terminal Avitag™ and His-tag, expressed in a HEK293 cell expression system. This protein is enzymatically biotin-labeled using Avitag™ technology. MW = 23.6 kDa (monomer). This protein runs at a higher MW by SDS-PAGE due to glycosylation. Biotinylation was confirmed to be $>90\%$.

Application:

Useful for studying protein binding, screening small molecules, and *in vitro* cellular assays.

Quality Assurance

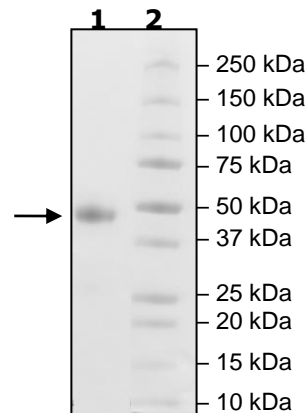
**4-20% SDS-PAGE
Coomassie staining**

Lane 1:
3 μg OX40

Lane 2:
Protein Marker

MW: 23.6 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