

Data Sheet

OX40L (CD252)

Human, recombinant, N-terminal His-tag

Catalog #: 71185

Lot #: 140708-D

Conc.: 1.93 mg/ml

Formulated in: 40 mM Tris-HCl, pH 8.0,
110 mM NaCl, 2.2 mM KCl, and 20%
glycerol.

Stability: >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 .*

Endotoxin: <1 EU/ μg

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:

Human OX40L, also known as tumor necrosis factor ligand superfamily member 4, TNFSF4, and CD252, GenBank Accession No. NM_003326, a.a. 51-183 with a N-terminal His-tag, expressed in a HEK293 cell expression system. MW = 18 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:

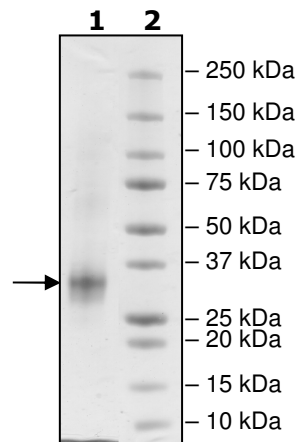
3 μg OX40L

Lane 2:

Protein Marker

MW: 18 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