

Data Sheet

LTBR, Mouse, Fc Fusion

Mouse, recombinant, Fc fusion protein

Catalog #: 71122

Lot #: 140318-2 **Conc.:** 2.76 mg/ml

Formulated in: 10 mM phosphate, 138 mM NaCl, 2.7 mM KCl, pH 7.4.

Stability: At least 6 months at -80°C . *Avoid freeze/thaw cycles.*

References:

1. Crowe, P.D., *et al.*, *Science*. 1994 Apr 29;**264**(5159):707-10.
2. Wu, M.Y., *et al.*, *J Biol Chem*. 1999 Apr 23;**274**(17):11868-73.

Description:

Mouse secreted lymphotoxin beta receptor (mLTBR)-Fc fusion protein, also known as, Tumor Necrosis Factor C Receptor (TNFCR), and CD18, GenBank Accession No. NM_010736, a.a 28-221, expressed in a HEK293 cell expression system. MW = 48.8 kDa (monomer). Endotoxin level ~ 0.0007 EU/ μg . This protein runs at a higher apparent M.W. by SDS-PAGE due to glycosylation.

Application:

Useful for the study of protein-protein interaction and in vitro and in vivo protein function studies.

Quality Assurance

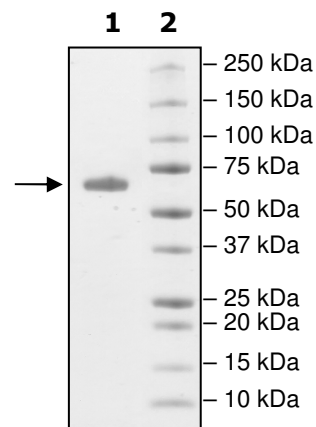
4-20% SDS-PAGE Coomassie staining

Lane 1:
2 μg TBR-Fc

Lane 2:
Protein Marker

MW: 48.8 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

Sequence:

KPQAPELRGSQPQLVPPYRIENQTCWDQDKEYYEPMHDVCCSRCPPGEFVFAVCSR
SQDTVCKTCPHNSYNEHWNHLSTCQLCRPCDIVLGFEEVAPCTSDRKAECRCQPGMS
CVYLDNECVHCEEERLVLCQPGTEAEVTDEIMDTDVNCVPCKPGHFQNTSSPRARCQ
PHTRCEIQGLVEAAPGTSYSDTICKNPPEPGAMPKSCDKTHTCPPCPAPPELLGGPSVFL
FPPKPKDTLMISRTPEVTCVVVDVSHEDPEVKFNWYVDGVEVHNAKTKPREEQYNSTY
RVVSVLTVLHQDWLNGKEYKCKVSNKALPAPIEKTISKAKGQPREPQVYTLPPSRDEL
KNQVSLTCLVKGFYPSDIAVEWESNGQPENNYKTTTPVLDSDGSFFLYSKLTVDKSRW
QQGNVFSCSVMHEALHNHYTQKSLSLSPGK

Green: mouse CD8a

Black: mouse LTBR (a.a. 28-221)

Red: Fc of human IgG1 (a.a.100-330) (#P01857)

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