

Data Sheet

TREM2, Fc-fusion (IgG1), Avi-Tag (Mouse)

Mouse, Recombinant, C-terminal Fc-Avi-tag

Catalog #: 79458**Lot #:** 180207**Conc:** 2.06 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. Storing diluted enzyme is not recommended, if necessary, use carrier protein (BSA 0.1 – 0.5%).

References:

1. Guerreiro, R., *et al.*, *New England Journal of Medicine*. 2013; **368**: 117-127.
2. Jonsson, T., *et al.*, *New England Journal of Medicine*. 2013; **368**: 107-116.

Description:

Mouse TREM2, also known as Triggering receptor expressed on myeloid cells 2 and Trem2a, GenBank Accession No. NM_031254, a.a. 19-168, fused at the C-terminus to the Fc portion of Human IgG1 and with a C-terminal Avi-tag, expressed in a Hek293 expression system. MW=45 kDa.

Quality Assurance

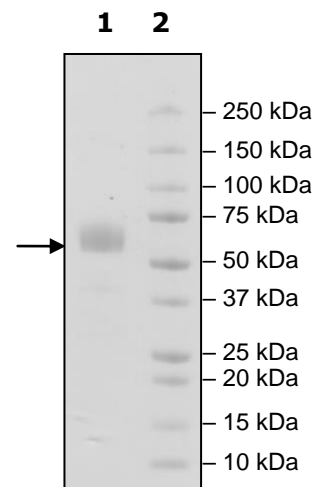
4-20% SDS-PAGE Coomassie staining

Lane 1:
3 μg TREM2(Mouse)

Lane 2:
Protein Marker

MW:
45 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 **1.858.202.1401** or **1.858.481.8694**, Fax **1.858.259.5734**

Or you can Email us at: info@bpsbioscience.com
Please visit our website at: www.bpsbioscience.com