

## Data Sheet

### **TREM2, Fc-fusion (IgG1), Avi-Tag (Mouse), Biotin**

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

**Catalog #:** 79459

**Lot #:** 180208

**Conc:** 0.65 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  $-80^{\circ}\text{C}$ . *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. This protein was biotinylated using AviTag™ technology. MW=45 kDa.

## **Quality Assurance**

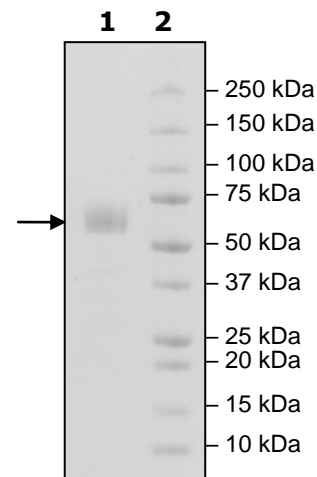
#### **4-20% SDS-PAGE Coomassie staining**

**Lane 1:**  
3µg TREM2(Mouse)

**Lane 2:**  
Protein Marker

**MW:**  
45kDa + glycans

**Purity:** ≥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](mailto:info@bpsbioscience.com)  
Please visit our website at: [www.bpsbioscience.com](http://www.bpsbioscience.com)