

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

Catalog: 79458

Lot: 180420-G

## Product Information

<b>Construct:</b>	TREM2 (19-168-Fc(IgG1)-Avi)
<b>Concentration:</b>	0.32 mg/ml
<b>Species:</b>	Mouse
<b>Formulated In:</b>	8 mM Phosphate, pH 7.4, 110 mM NaCl, 2.2 mM KCl, and 20% Glycerol
<b>Expression System:</b>	HEK293
<b>Format:</b>	Aqueous buffer solution
<b>Stability:</b>	At least 6 months at -80°C. Avoid freeze/thaw cycles.
<b>Storage:</b>	-80°C
<b>Genbank Accession:</b>	NM_031254
<b>MW:</b>	45 kDa + glycans
<b>Glycosylation:</b>	This protein runs at a higher MW by SDS-PAGE due to glycosylation.
<b>Purity:</b>	≥90%

## Quality Control Data

### 4-20% SDS-Page Coomassie Staining

