

Data Sheet

TIM-3, (Mouse), Fc-Fusion (Human) Avi-Tag, Biotin

Mouse, Recombinant, C-terminal Fc-Avi-Tag,
 Biotin-Labeled
Catalog#: 77000
Lot#: 170817 **Conc.:** 0.72 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. Protein may be diluted to
 $\geq 100 \mu\text{g/ml}$ in PBS + glycerol and stored at
 -80°C .*

References:

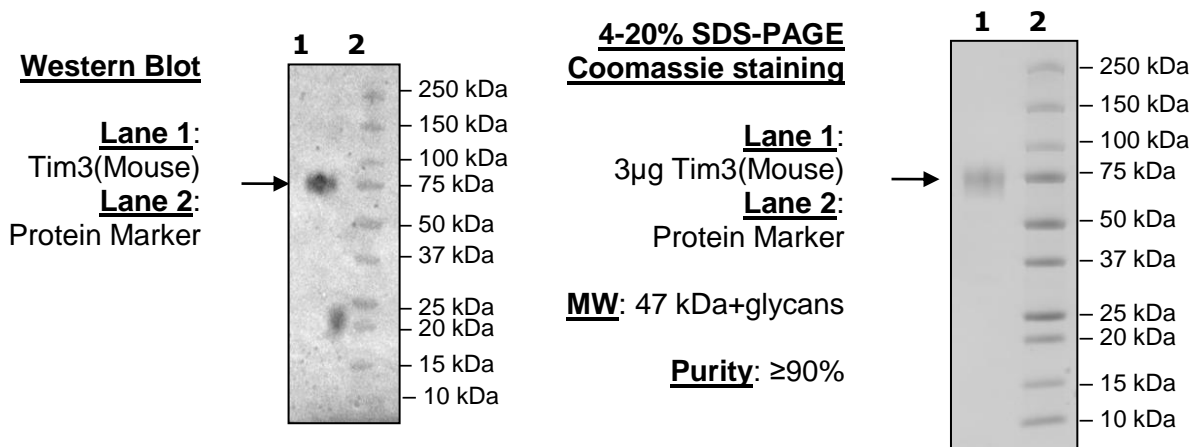
1. Zhu, C. *et al.*, *Nat. Immunol.* 2005;
6:1245-1252.
2. Anderson, A.C. *et al.*, *Science.* 2007;
318(5853):1141-1143.

Description:

Mouse TIM-3, also known as T-cell immunoglobulin mucin receptor 3, T-cell membrane protein 3, TIMD-3, Hepatitis A virus cellular receptor 2, HAVCR-2, and KIM-3. GenBank Accession No. NM_032782, a.a. 22-191, expressed in a HEK293 cells expression system. This protein is fused at the C-terminus to the Fc region of human IgG1 and is enzymatically biotinylated using Avi-tag™ technology. MW = 47 kDa. This protein runs at a higher MW by SDS-PAGE due to glycosylation.

Applications: Useful as a positive control for western blot studies.

Quality Assurance



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