

Data Sheet

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

Mouse, Recombinant, C-terminal Fc-Avi-Tag
Catalog#: 72549
Lot#: 170830 **Conc.:** 0.15 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 cell expression system. This protein is fused at the C-terminus to the Fc region of human IgG1. 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

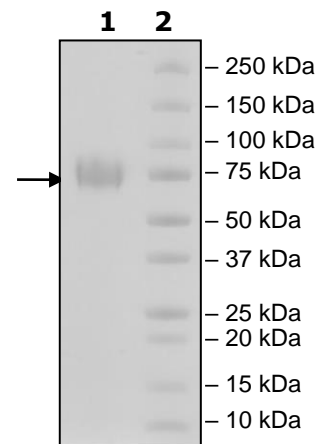
**4-20% SDS-PAGE
Coomassie staining**

Lane 1:
4 μg Tim3(Mouse)

Lane 2:
Protein Marker

MW: 47 kDa+glycans

Purity: $\geq 90\%$



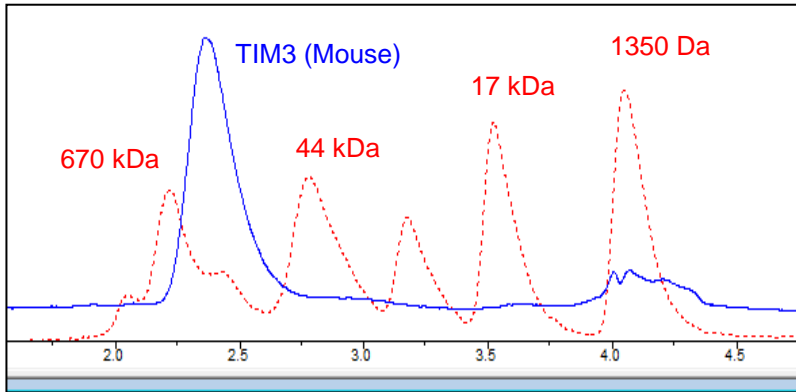
Gel Filtration Curve

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



6042 Cornerstone Court W, Ste B
San Diego, CA 92121
Tel: 1.858.829.3082
Fax: 1.858.481.8694
Email: info@bpsbioscience.com



<10% Aggregation

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