

Data Sheet

TIM-3, Fc Fusion

Human, Recombinant, Fc fusion protein

Catalog #: 71151

Lot #: 140414

Conc.: 0.42 mg/ml

Formulated in: 40 mM Tris, pH 8.0, 110 mM NaCl, 2.4 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 .

Endotoxin Level: $<0.1 \text{ EU}/\mu\text{g}$

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:

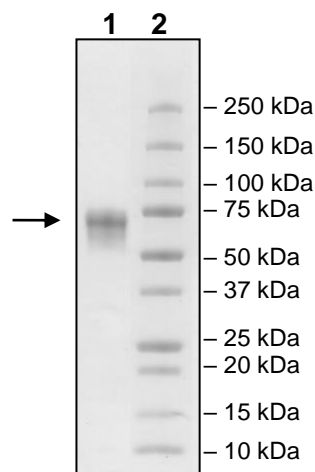
Human secreted TIM-3, Fc fusion protein, also known as T-cell immunoglobulin mucin receptor 3, T-cell membrane protein 3, T-cell and immunoglobulin and mucin domain-containing protein 3, TIMD-3, Hepatitis A virus cellular receptor 2, and HAVCR-2. GenBank Accession No. NM_032782.4, a.a. 22-200 expressed in a HEK293 cell expression system. MW = 47 kDa. This protein runs at a higher MW by SDS-PAGE due to glycosylation.

Quality Assurance

4-20% SDS-PAGE Coomassie staining

Lane 1:
3 μg TIM-3
Lane 2:
Protein Marker

MW: 47 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 **1.858.829.3082** 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

Sequence

SEVEYRAEVGQNAYLPCFYTPAAPGNLVPVCWGKGACPVFECGNVLRDTERDVNY
WTSRYWLNDFRKGDVSLTIENVTLADSGIYCCRIQIPGIMNDEKFNKLVKPAKVTP
APTRQRDFTAAPRMLTTRGHGPAETQTLGSLPDINLTQISTLANELRDSRLANDLRD
SGATIRIEGRMDPKSCDKTHTCPPCPAPELLGGPSVFLFPPKPKDTLMISRTPEVTCVV
VDVSHEDPEVKFNWYVDGVEVHNAKTKPREEQYNSTYRVVSVLTVLHQDWLNGKEYK
CKVSNKALPAPIEKTISKAKGQPREPQVYTLPPSRDELTKNQVSLTCLVKGFYPSDIAVE
WESNGQPENNYKTTTPVLDSGDSFFLYSKLTVDKSRWQQGNVVFSCVMHEALHNHY
TQKSLSLSPGK

Black: Tim3 (a.a. 22-200)

Green: linker

Red: Fc of human IgG1 (a.a.100-330)

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