

Data Sheet

TAC1, Fc Fusion (IgG1) Avi-Tag, Biotin Labeled

Human, Recombinant, Fc-Avi-Tag,
Biotin Labeled

Catalog #: 100284

Lot #: 190214

Conc.: 0.44 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.

References:

1. Coste, A. T. *et al.*, *Eukaryotic Cell*.
2004; **3(6)**: 1639-1652.
2. Smith, S. T. *et al.*, *Nature Cell Bio*.
2004; **6**: 162-167.

Description: Human TAC1, also known as Hs.2563, NK2, NKNA, NPK, TAC2, and tachykinin precursor 1. GenBank Accession No. NM_0012452, a.a. 23-166, fused at the C-terminus of the Fc portion of Human IgG1, with C-terminal Avi-tag™, expressed in a HEK293 cell expression system. MW = 48 kDa.

Quality Assurance

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

Lane 1:

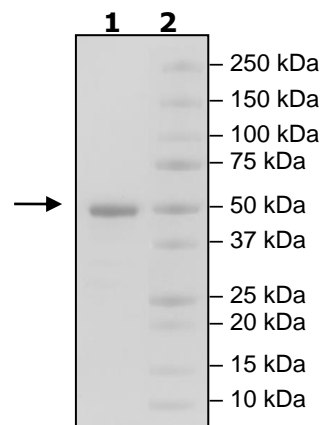
4 μg TAC1

Lane 2:

Protein Marker

MW: 48 kDa

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 **1.858.202.1401** Fax **1.858.481.8694**

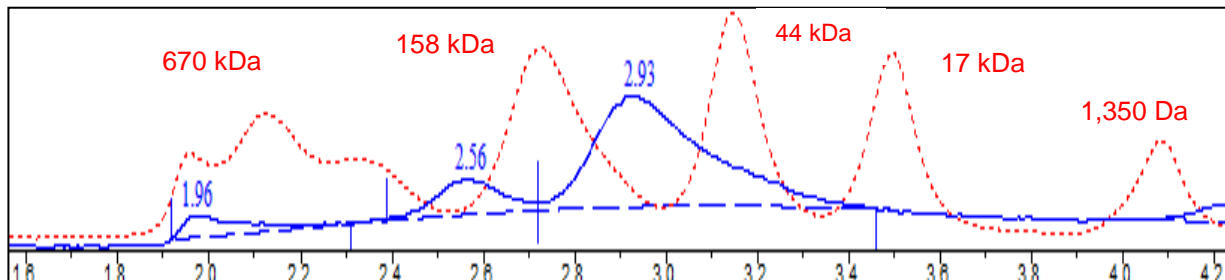
Or you can Email us at: info@bpsbioscience.com

Please visit our website at: www.bpsbioscience.com



6042 Cornerstone Court W, Ste B
San Diego, CA 92121
Tel: 1.858.202.1401
Fax: 1.858.481.8694
Email: info@bpsbioscience.com

Gel Filtration:



< 10% Aggregation

OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.202.1401** Fax **1.858.481.8694**
Or you can Email us at: info@bpsbioscience.com
Please visit our website at: www.bpsbioscience.com