

## Data Sheet

### TAC1, Fc-Fusion (IgG1), Avi-Tag

Human, Recombinant, C-terminal Fc-Avi-Tag

**Catalog #:** 100283

**Lot#:** 190205

**Conc.:** 0.79 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  $-80^{\circ}\text{C}$ . Avoid freeze/thaw cycles.

**Description:** Human TAC1 also known as Tachykinin Precursor 1, Preprotachykinin Neuropeptide Gamma, Substance K, and Neurokinin 1, GenBank Accession No. NM\_012452, a.a. 23-166, fused at the C-terminus to the Fc portion of human IgG1 followed by a C-terminal Avi-tag™, expressed in a HEK293 cell expression system. MW=45 kDa.

#### 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.

---

### **Quality Assurance**

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

**Lane 1:**

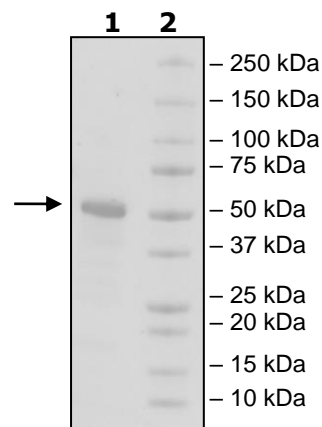
4 $\mu\text{g}$  TAC1

**Lane 2:**

Protein Marker

**MW:** 45 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](mailto:info@bpsbioscience.com)

Please visit our website at: [www.bpsbioscience.com](http://www.bpsbioscience.com)