

Data Sheet

VSIG3, Fc-fusion (IgG1), Avi-Tag, Biotin-Labeled

Human, Recombinant, C-terminal
Fc fusion protein, Avi-Tag, Biotin-Labeled

Catalog #: 100044

Lot #: 180412

Conc.: 0.91 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. Katoh, L., *et al.*, *International Journal of Oncology*. 2003; **23**: 525-531.
2. Watanbe, T., *et al.*, *Cancer Science*. 2005; **10**: 17858-17863.

Description:

Human VSIG3, also known as V-set and immunoglobulin domain containing 3, IgSF11, GenBank Accession No. NM_152538.3, a.a. 22-244, with C-terminal Avi-TagTM fused to Fc region of Human IgG1, expressed in a HEK293 cell expression system. MW = 52 kDa. This protein is enzymatically biotinylated using Avi-TagTM technology.

Quality Assurance

4-20% SDS-PAGE Coomassie staining

Lane 1:

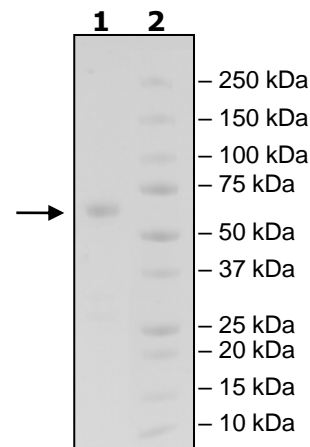
4 μg VSIG8

Lane 2:

Protein Marker

MW: 52 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: support@bpsbioscience.com

Please visit our website at: www.bpsbioscience.com