

Data Sheet

VSIG3, Avi-His-Tag

Human, Recombinant protein, C-terminal
Avi-His-Tag

Catalog #: 100349

Lot #: 190424-2

Conc.: 1.07 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. Katoh, L., *et al.*, *International Journal of Oncology*. 2003; **23**: 525-531.
2. Watanbe, T., *et al.*, *Cancer Science*. 2005; **10**: 17858-17863.

Description:

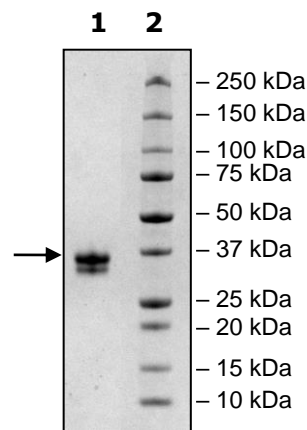
Human VSIG3, also known as V-set, immunoglobulin domain containing 3, and IgSF11, GenBank Accession No. NM_152538.3, a.a. 22-245, with C-terminal Avi-His-Tag™, expressed in a HEK293 cell expression system. MW = 26 kDa.

Quality Assurance

4-20% SDS-PAGE Coomassie staining

Lane 1:
3 μg VSIG3
Lane 2:
Protein Marker

MW: 26 kDa
Purity: $\geq 70\%$



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