

Data Sheet

VSIG8, Fc Fusion, Avi-Tag

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

Catalog #: 40743

Lot #: 170512-G **Conc.:** 0.42 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.*

Reference(s):

1. Rice, R. *et al. Journal of Invest. Dermatology* (2011) **131**: 9. 1936-1938.

Description: Human VSIG8, also known as V-set and immunoglobulin domain containing 8, Genbank accession number NM_001013661, a.a. 22-263, fused at the C-terminus with the Fc portion of human IgG1, with C-terminal Avi-tag™, expressed in a HEK293 expression system. MW=56 kDa.

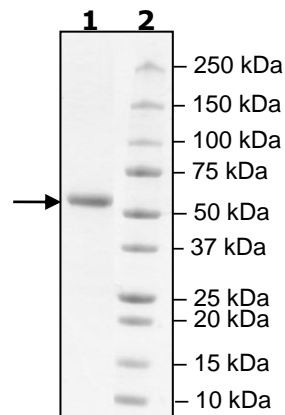
Quality Assurance

4-20% SDS-PAGE Coomassie staining

Lane 1:
2 μg VSIG8

Lane 2:
Protein Marker

MW: 56 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