

Data Sheet

Frizzled-1, Fc Fusion

Human, recombinant, Fc fusion protein

Catalog #: 71123

Lot#: 140226 **Conc.:** 0.29 mg/ml

Formulated in: Sterile-filtered (0.2 µm) solution of 40 mM Tris, pH 8.0, 110 mM NaCl, 2.2 mM KCl, and 20% glycerol.

Stability: >6 months at -80°C. Avoid freeze/thaw cycles.

References:

1. Moffat LL., *et al.*, *Development*. 2014 Feb; **141(3)**:617-28.
2. Surana, R., *et al.*, *Biochim Biophys Acta*. 2014 Jan; **1845(1)**:53-65.

Description:

Secreted frizzled receptor (FZD1)-Fc fusion protein and WNT receptor, also known as FZD1, GenBank Accession No.BC_051271, a.a. 73-253, expressed in a HEK293 cell expression system. MW = 46.4 kDa (monomer). This protein runs at a higher apparent M.W. by SDS-PAGE due to glycosylation.

Application:

Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

Quality Assurance

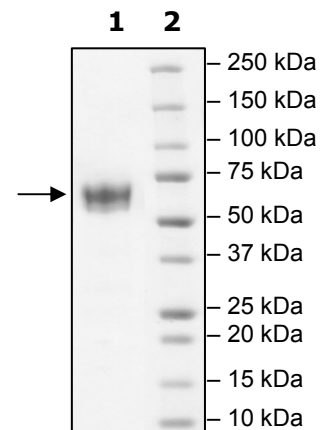
**4-20% SDS-PAGE
Coomassie staining**

Lane 1:
2 µg sFZD1-Fc

Lane 2:
Protein Marker

MW: 46.4 kDa

Purity: ≥90%



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

To place your order, please contact us by Phone **1.858.829.3082** Fax **1.858.481.8694**

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

Please visit our website at: www.bpsbioscience.com



6044 Cornerstone Court W, Ste E
San Diego, CA 92121
Tel: 1.858.829.3082
Fax: 1.858.481.8694
Email: info@bpsbioscience.com

Sequence:

QAAGQGPGQGPGPGQQPPPPPPQQQQSGQQYNGERGISVDPDHGYCQPISIPLCTDIAY
NQTIMPNULLGHTNQEDAGLEVHQFYPLVKVQCSAELKFFLCSMYAPVCTVLEQALPPC
RSLCERARQGCEALMNKFGFQWPDTLKCEKFPVHGAGELCVGQNTSDKGTPTPSLLP
EFWTSNPQHIEGRMDPKSCDKTHTCPPCPAPELLGGPSVFLFPPKPKDTLMISRTPEV
TCVVVDVSHEDPEVKFNWYVDGVEVHNAKTKPREEQYNSTYRVVSVLTVLHQDWLNG
KEYKCKVSNKALPAPIEKTISKAKGQPREPQVYTLPPSRDELTKNQVSLTCLVKGFYPS
DIAVEWESNGQPENNYKTTPPVLDSDGSFFLYSKLTVDKSRWQQGNVFCFSVMHEAL
HNHYTQKSLSLSPGK

Black: FZD1 (aa 73-253)

Green: linker

Red: Fc of human IgG1 (aa100-330) (#P01857)

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

To place your order, please contact us by Phone **1.858.829.3082** Fax **1.858.481.8694**
Or you can Email us at: info@bpsbioscience.com
Please visit our website at: www.bpsbioscience.com