

## Data Sheet

### Frizzled-4, Fc Fusion

Human, recombinant, Fc fusion protein

Catalog #: 71126

Lot#: 140219                      Conc.: 0.97 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 (FZD4)-Fc fusion protein and WNT receptor, also known as FZD4, GenBank Accession No. BC\_114527, a.a. 37-180, expressed in a HEK293 cell expression system. MW = 42.9 kDa (monomer). This protein runs at a higher apparent M.W. by SDS-PAGE due to glycosylation.

#### Application:

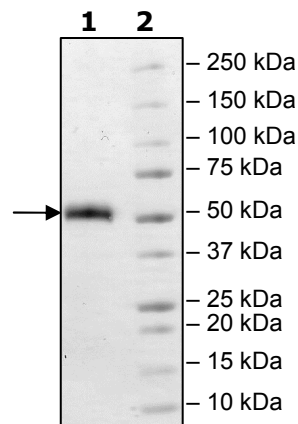
Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

## Quality Assurance

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

**Lane 1:**  
2 µg Fzd4  
**Lane 2:**  
Protein Marker

**MW:** 42.9 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](mailto:info@bpsbioscience.com)  
Please visit our website at: [www.bpsbioscience.com](http://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](mailto:info@bpsbioscience.com)

**Sequence:**

FGDEEERRCDPIRISM CQNLGYNVTKMPNLVGH ELQTD AELQLTTFTPLIQYGCSSQLQ  
FFLCSVYVPMCTEKINIPIGPCGGMCLSVKRRCEPVLKEFGFAWPESLNCSKFPPQND  
HNHMCMEGPGDEEVPLPHKTPIQPGEEIEGRMDPKSCDKTHTCPPCPAPELLGGPSV  
FLFPPKPKDTLMISRTPEVTCVVVDVSHEDPEVKFNWYVDGVEVHNAKTKPREEQYNS  
TYRVVSVLTVLHQDWLNGKEYKCKVSNKALPAPIEKTISKAKGQPREPQVYTLPPSRDE  
LTKNQVSLTCLVKGFYPSDIAVEWESNGQPENNYKTTTPVLDSDGSFFLYSKLTVDKSR  
WQQGNVFSCSVMHEALHNHYTQKSLSLSPGK

Black: FZD4 (aa 37-180)

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