

## Data Sheet

### Frizzled-7, Fc Fusion

Human, recombinant, Fc fusion protein

**Catalog #:** 71139

**Lot#:** 140408      **Conc.:** 0.44 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 (FZD7)-Fc fusion protein and WNT receptor, also known as FZD7, GenBank Accession No. BC\_015915, a.a. 33-185, expressed in a HEK293 cell expression system. MW = 43.3 kDa (monomer). This protein runs at a higher apparent M.W. by SDS-PAGE due to glycosylation.

#### Application:

Useful for binding studies, screening inhibitors, and selectivity profiling.

## Quality Assurance

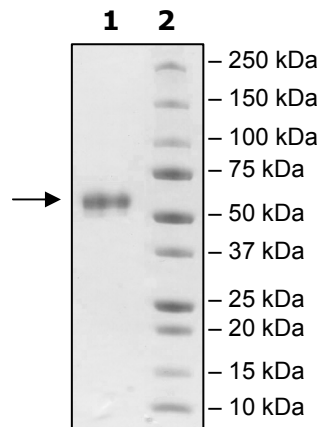
### 4-20% SDS-PAGE Coomassie staining

**Lane 1:**  
2 µg sFZD7

**Lane 2:**  
Protein Marker

**MW:** 43.3 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:**

QPYHGEGKGISVPDHGFCQPISIP LCTDIAYNQ TILPNLLGHTNQEDAGLEVHQFYPLVKV  
QCSPELRFFLC SMYAPVCTVLDQAIPPCRSLCERARQGCEALMNKFGFQWPERLRCE  
NFPVHGAGEICV GQNTSDGSGGGPGGGPTAYPTAPYL **IEGRMDPKSCDKTHTCPPCPA**  
**PELLGGPSVFLFPPKPKDTLMISRTPEVTCVVDVSHEDPEVKFNWYVDGVEVHNAKT**  
**KPREEQYNSTYRVVSVLTVLHQDWLNGKEYKCKVSNKALPAPIEKTISKAKGQPREPQ**  
**VYTLPPSRDELTKNQVSLTCLVKGFYPSDIAVEWESNGQPENNYKTTTPVLDSDGSFFL**  
**YSKLTVDKSRWQQGNV FSCSVMHEALHNHYTQKSLSLSPGK**

Black: FZD7 (a.a. 33-185)

Green: linker

Red: Fc of human IgG1 (a.a. 100-330) (#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)