

## Data Sheet

### FCGR2B, Avi-His-Tag, Biotin-Labeled

Human, Recombinant, C-terminal Avi-His-Tag  
Biotin labeled

**Catalog #:** 100474

**Lot#:** 190801

**Conc.:** 1.02 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  $-80^{\circ}\text{C}$ . *Avoid  
freeze/thaw cycles.*

**Description:** Human FCGR2B, known as CD32B, and FCG2, GenBank Accession No. NM\_004001.4, a.a. 46-217, with C-terminal Avi-His-Tag, expressed in a HEK293 cell expression system and enzymatically biotin-labeled using Avi-tag™ technology. Biotinylation is confirmed to be  $\geq 90\%$ . MW = 23 kDa. This protein runs at a higher MW by SDS-PAGE due to glycosylation.

#### References:

1. Kyogoku, C., *et al.*, *Arthritis and Rheumatology*. 2002; **46(5)**: 1242-1254.
2. Su, K. *Journal of Immunology*. 2004; **172(11)**:4385-4389.

### Quality Assurance

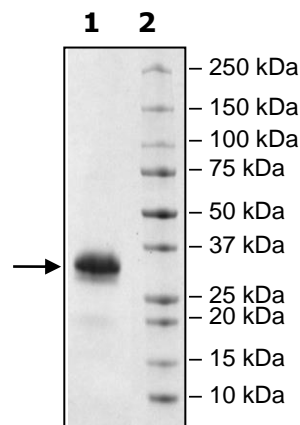
#### 4-20% SDS-PAGE Coomassie staining

**Lane 1:**  
4  $\mu\text{g}$  FCGR2B

**Lane 2:**  
Protein Marker

**MW:** 23 kDa+glycans

**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](mailto:info@bpsbioscience.com)

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