

Data Sheet

FCGR2B, Avi-His-Tag

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

Catalog #: 100089

Lot#: 180516

Conc.: 0.78 mg/ml

Formulated in: 8mM phosphate, pH 7.4, 110mM NaCl, 2.2mM KCl, and 20% glycerol.

Stability: At least 6 months at -80°C . Avoid freeze/thaw cycles. Protein may be diluted to $\geq 100 \mu\text{g/ml}$ in PBS + glycerol and stored at -80°C .

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.

Description:

Human FCGR2B, known as CD32B, and FCG2, GenBank Accession No. NM_004001.4, a.a. 46-217, with C-terminal AviTM-His-Tag, expressed in a HEK293 cell expression system. This protein runs at a higher MW by SDS-PAGE due to glycosylation. MW = 23 kDa.

Applications:

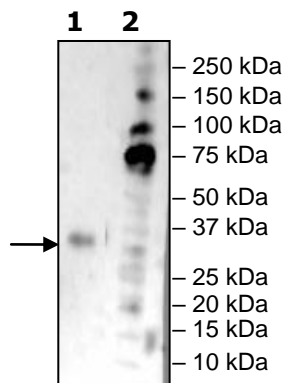
Useful as a positive control for western blot studies.

Quality Assurance

Western Blot

Lane 1:
FCGR2B
Lane 2:
Protein Marker

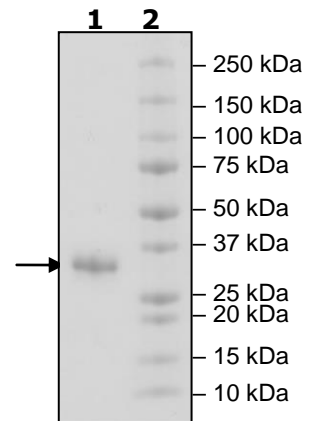
MW:
23 kDa+glycans



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

Lane 1:
5 μg 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 **1.858.829.3082** Fax **1.858.481.8694**
 Or you can Email us at: info@bpsbioscience.com
 Please visit our website at: www.bpsbioscience.com