

Data Sheet

FCGR3B, Avi-His-Tag

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

Catalog #: 79016

Lot#: 180517

Conc.: 0.81 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. Fanciulli, M., et al., Nature Genetics. 2007; 39: 721-723.
2. Breunis, W. Human Mutation. 2009; 30(5):E640-E650.

Description:

Human FCGR3B, known as CD16b, and FCG3, GenBank Accession No. NM_000570.4, a.a. 20-208, 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 = 25 kDa.

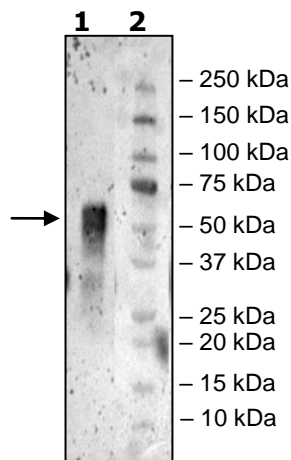
Applications:

Useful as a positive control for western blot studies.

Quality Assurance

Western Blot

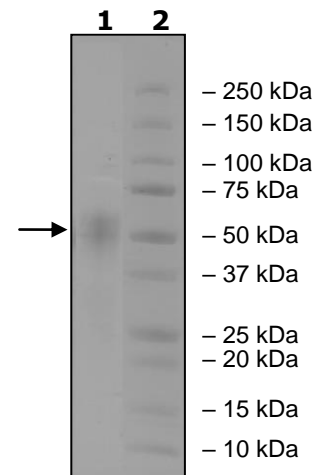
Lane 1:
FCGR3B
Lane 2:
Protein Marker



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

Lane 1:
4 μg FCGR3B
Lane 2:
Protein Marker

MW: 25 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