

Data Sheet

FCGR3A (F176V), Avi-His-Tag

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

Catalog #: 100079

Lot#: 180517

Conc.: 1.44 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°C . Avoid freeze/thaw cycles.

References:

1. Zhang, W., *et al.*, *Journal of Clinical Oncology*. 2007; **25(24)**: 3712-3718.
2. Dall'Ozzo, S., *et al.*, *Cancer Research*. 2004; **64(13)**: 4664-4669.

Description: Human FCGR3A, known as Fc gamma RIIIA, and CD16a with F176V mutation, GenBank Accession No. NM_001127593.1, a.a. 17-208, with a C-terminal Avi-His-Tag™, expressed in a HEK293 cell expression system. MW = 25 kDa. This protein runs at a higher MW by SDS-PAGE due to glycosylation.

Quality Assurance

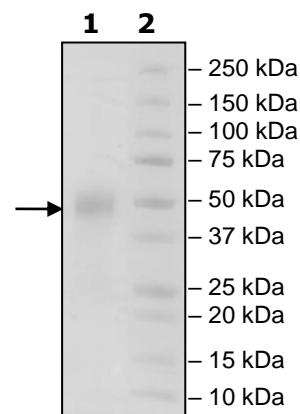
4-20% SDS-PAGE Coomassie staining

Lane 1:
4 μg FCGR3A

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 **1.858.202.1401** Fax **1.858.481.8694**

Or you can Email us at: info@bpsbioscience.com

Please visit our website at: www.bpsbioscience.com