

Data Sheet

FCGR3A, Avi-His-Tag

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

Catalog #: 79013

Lot#: 180516

Conc.: 2.26 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 .

Description:

Human FCGR3A, known as Fc gamma RIIIA, and CD16a, GenBank Accession No. NM_001127593.1, a.a. 17-208, with N-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.

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.

Quality Assurance

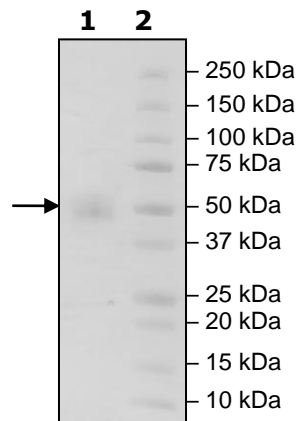
4-20% SDS-PAGE Coomassie staining

Lane 1:
3 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 **1.858.829.3082** Fax **1.858.481.8694**

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

Please visit our website at: www.bpsbioscience.com