

Data Sheet

CD47, Fc Fusion, Avi-Tag, Biotin-Labeled (Mouse)

Mouse, Recombinant, Fc fusion, C-Terminal
Avi-tag, Biotin-Labeled

Catalog #: 72514

Lot #: 170518-G

Conc.: 0.53 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. Protein may be diluted to $\geq 100 \mu\text{g/ml}$ in PBS + glycerol and stored at -80°C .*

References:

1. Vernon-Wilson, E.F., *et al.*, *Eur. J. Immunol.* 2000; **30(8)**: 2130-2137.
2. Majeti, R. *et al.*, *Cell.* 2009; **138(2)**: 286-299.

Description:

Mouse CD47, also known as leukocyte surface antigen CD47, Rh-related antigen, integrin-associated protein (IAP), antigenic surface determinant protein OA3, and MER6. GenBank Accession No. NM_010581.3, a.a. 19-158 fused at the C-terminus to the Fc portion of human IgG1 and C-terminal Avi-tag expressed in a HEK293 cell expression system. The protein is enzymatically biotinylated using AviTag™ technology. MW = 44 kDa. This protein runs at a higher MW by SDS-PAGE due to glycosylation.

Applications: Useful for the study of protein-protein interaction and in vitro and in vivo protein function studies.

Quality Assurance

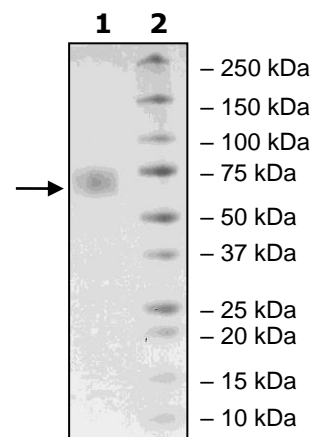
**4-20% SDS-PAGE
Coomassie staining**

Lane 1:
3 μg CD47

Lane 2:
Protein Marker

MW: 44 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