

Data Sheet

Anti-CD47 Antibody, PE-Labeled

Recombinant, Human Monoclonal
Anti-CD47

Catalog #: 79105-2

Lot #: 170707

Conc.: 2.10 mg/ml

Host Species | Isotype:

Human | Human IgG1

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

Stability: Stable for at least 12 months at
-80°C. *Protect from light. Avoid freeze/thaw
cycles. Store diluted antibody with carrier
protein (BSA 0.1 – 0.5%) at 4°C, stable for 2
weeks.*

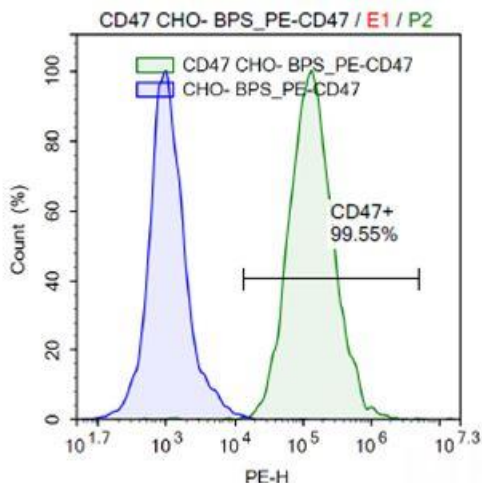
Description: R-Phycoerythrin-labeled anti-
CD47 recombinant human monoclonal
recognizing the SIRP- α or SIRP- γ binding
region of human CD47. This antibody has not
been tested for cross-reactivity with any other
species.

Purification: Protein A affinity
chromatography from HEK293 supernatants.

Application: Useful for labeling cells
expressing CD47 for flow cytometry and
immunofluorescence microscopy.

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.



Control CHO cells and CD47 expressing stable CHO cells were treated with human PE anti-CD47 antibody.

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