

Data Sheet

Anti-CD47 Antagonist Antibody

Recombinant, Human Monoclonal
Anti-CD47

Catalog #: 79065-2

Lot #: 170619

Conc.: 1.16 mg/ml

Host Species | Isotype:

Human | Human IgG1

Formulated in: 8 mM 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.*

References:

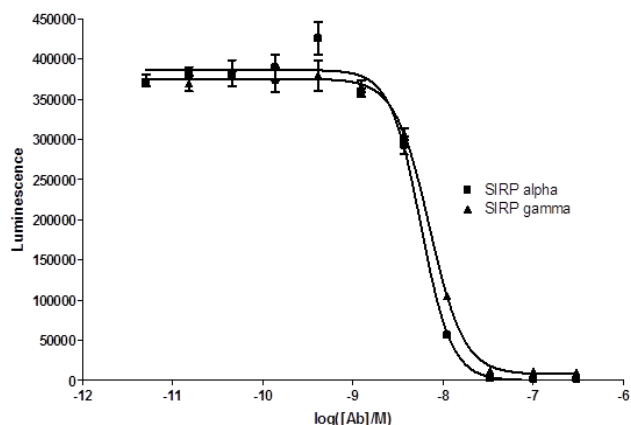
1. Vernon-Wilson, E.F. et al., Eur. J. Immunol. 2000; 30(8): 2130-2137.
2. Majeti, R. et al., Cell. 2009; 138(2): 286- 299.

Description: Anti-CD47 recombinant human antagonist antibody 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 as an antagonist to block CD47 signaling and binding to SIRP- α or SIRP- γ in biochemical assays.

CD47: SIRP α and SIRP γ



Coated wells in assay plate with either PBS (No Coat) or 100 ng/well of CD47-His in 50 μ l PBS (overnight at 4°C).

Anti CD47 was diluted to 1.5 μ M in Assay buffer. A series of (1:3) serial dilution was made in Assay buffer. To 30 μ l of Assay Buffer, 10 μ l of inhibitor was added to each well, except no coat and positive control wells.

10 μ l of SIRP α -biotin (600ng/reaction) or SIRP γ -biotin (40ng/reaction) in assay buffer was added to each well. Incubated for 2 hours at room temperature. Detected binding with Strep-HRP generated luminescence. Assay Buffer is PBS + 0.1% BSA.

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