

## Data Sheet

### Anti-CD47 Antagonist Antibody

Recombinant, Human Monoclonal  
Anti-CD47

**Catalog #:** 79065-1

**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.*

**Description:** Anti-CD47 recombinant human  
antagonist antibody recognizing the SIRP- $\alpha$  or  
SIRP- $\gamma$  binding region of human CD47. This  
antibody has not been tested for cross-reactivity  
with any other species.

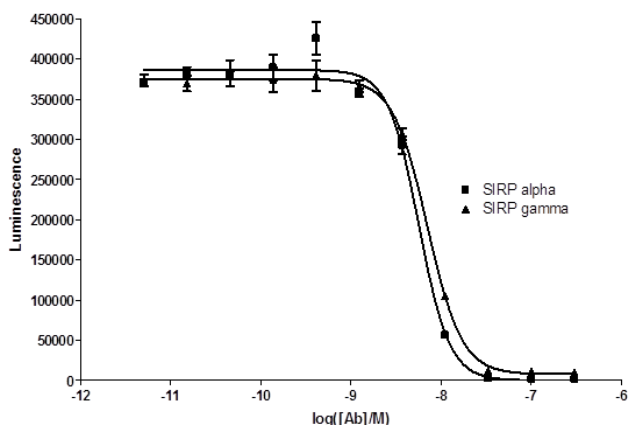
**Purification:** Protein A affinity chroma-  
tography from HEK293 supernatants.

**Application:** Useful as an antagonist to block  
CD47 signaling and binding to SIRP- $\alpha$  or  
SIRP- $\gamma$  in biochemical assays.

#### 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.

#### CD47: SIRP $\alpha$ and SIRP $\gamma$



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

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

10  $\mu$ l of SIRP $\alpha$  -biotin (600ng/reaction) or SIRP  $\gamma$ -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](mailto:info@bpsbioscience.com)

Please visit our website at: [www.bpsbioscience.com](http://www.bpsbioscience.com)