

Data Sheet

Anti-CD28 Agonist Antibody

Recombinant, Monoclonal Anti-CD28
Catalog #: 100182
Lot #: 180821 **Conc.:** 1.05 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.*

References:

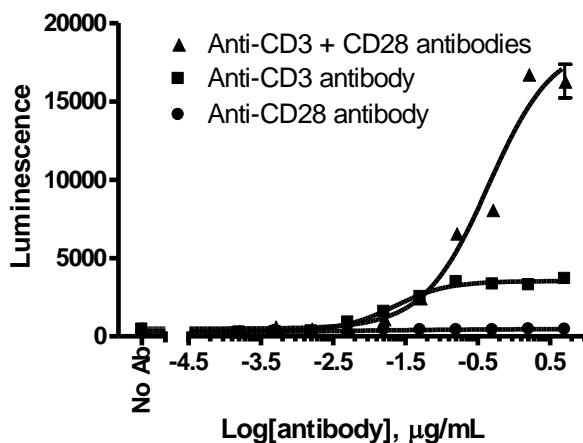
1. Takemura, S., et al., *FEBS Letters*, 2000; **476**: 266-271.

Description: The T cell-specific receptor CD28 is an important regulator of the immune system. The ligation of CD28 and its ligands CD80/CD86 on the APC provides a critical costimulatory signal alongside T cell receptor (TCR) for T cell activation. The costimulatory signal can also be delivered by the interaction between CD28 and its agonist antibody.

Purification: Protein A affinity chromatography from HEK293 supernatants.

Application: This antibody is useful for activation of CD28 costimulatory pathway in T cells

Costimulatory Activity of Anti-CD28 Antibody on IL-2 reporter/Jurkat Cells



Activation Assay: Serial dilutions of anti-CD3 (BPS Bioscience, #71274) and CD28 antibodies were added to the IL-2 reporter-Jurkat cells, then incubated at 37°C incubator for 16 hrs. After the treatment, perform Luciferase assay using One-Step Luciferase assay system (BPS Bioscience, #60690)

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