

Data Sheet

Anti-CD27 Agonist Antibody

Recombinant, Human Monoclonal
Anti-CD27

Catalog #: 100111

Lot #: 180619 **Conc.:** 1.6 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:

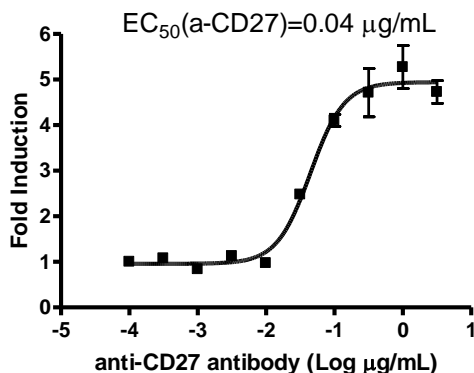
1. Vitale, L., *et al.*, *Clin Cancer Res.* 2012;
18(14): 3812-3821.

Description: Anti-CD27 agonist antibody binds to CD27 on T cells and promotes T cell response. It represents new opportunities for cancer immunotherapy.

Purification: Protein A affinity chromatography from HEK293 supernatants.

Application: This antibody crosslinked with FcGR2B is useful for activation of CD27 costimulatory pathway in T cells.

Agonist Activity of Anti-CD27 Antibody on CD27/NF-κB-Jurkat Cells (#79509)



Activation Assay: CD27/NF-κB-Jurkat cells (#79509) were cocultured with FcGR2B-CHO cells (#79511), and serial dilutions of anti-CD27 antibody were added to the cells, then incubated at 37°C incubator for 16 hrs. After the treatment, perform Luciferase assay using ONE-Step™ Luciferase assay system (#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