

Data Sheet

Anti-CD137 Agonist Antibody

Recombinant, Human, Monoclonal

Catalog #: 79097-1

Lot #: 170718 **Conc.:** 0.08 mg/ml

Host Species | Isotype:

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. Avoid freeze/thaw cycles. Store diluted antibody with carrier protein (BSA 0.1 – 0.5%) at 4°C, stable for 2 weeks.

References:

1. Wilcox, R.A., et al., J. Clin. Invest. 2002; 109(5): 651-659.
2. Foell, J. et al., J. Clin. Invest. 2003; 111(10): 1505-1518.

Description: Anti-CD137 recombinant human antibody recognizing the CD137 ligand binding region. This antibody not been tested for cross-reactivity with other species.

Purification: Protein A affinity chromatography from HEK293 supernatants.

Application: CD137 agonist antibody prevents tumor recurrence and metastases in cancer patients. This antibody is useful for activation of CD137 costimulatory pathway in T cells.

Dose Responsive Curve of CD137/NFkB-HEK293 Cell Line

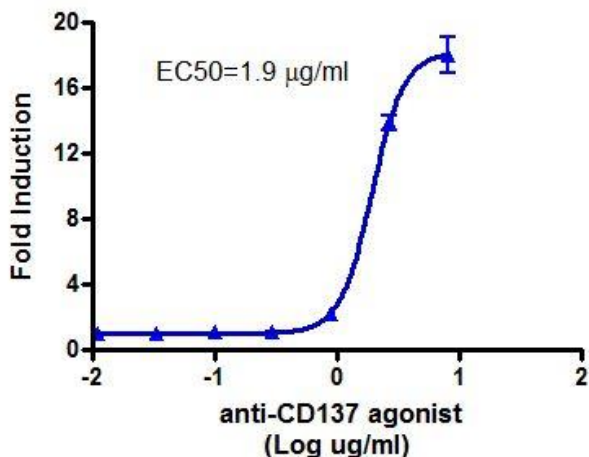


Fig.1 Anti-CD137 agonist (BPS #79097) was diluted and added to the cells (BPS #79289), then incubated at 37°C cell culture incubator for 6 and 24 hrs. respectively. After the treatment, Luciferase assay was performed using One-Step Luciferase Assay System (BPS #60690).

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