

Data Sheet

CTLA4 Neutralizing Antibody

Recombinant, Human Monoclonal Antibody
 Anti-CTLA4

Catalog#: 71212

Lot#: 160803

Conc.: 1.67 mg/ml

Host Species | Isotype:

Human | IgG1k

Formulated in: 8 mM phosphate, 110 mM NaCl, 2.2 mM KCl, pH 7.4, and 20% glycerol

Endotoxin Level: <1 EU/μg

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. Linsley, P.S., *et al.*, *J. Exp. Med.* 1991; **174(3)**: 561-569.
2. Hurwitz, A.A., *et al.*, *Canc. Res.* 2000; **60**: 2444-2448.

Description: Neutralizing recombinant human monoclonal antibody recognizing the extracellular domain of CTLA4 and blocking binding of CTLA4 to B7.1 and B7.2. This antibody recognizes human CTLA4 but does not bind to murine CTLA4. This antibody has not been tested for cross-reactivity with other species.

IC50: 5 nM using CTLA4:B7-1[Biotinylated] Inhibitor Screening Assay Kit (BPS Bioscience, Cat. #72009).

Assay Conditions: Antibody was added to CTLA4:B7.1 binding assay at various concentrations. Assay was done according to BPS Cat. #72009 protocol.

Purification: Protein A affinity chromatography from HEK293 supernatants.

Applications: Useful for inhibition of CTLA4:B7 interactions.

Image(s):

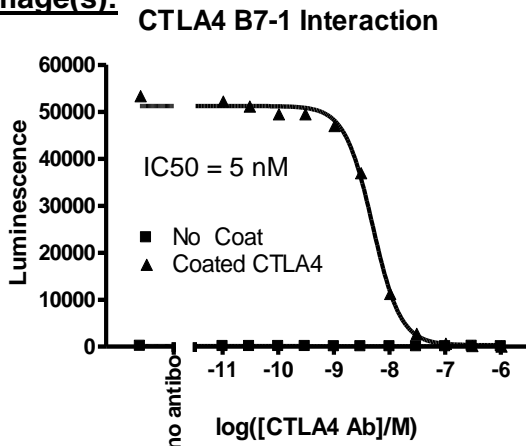


Figure 1. Wells were coated with 100 ng of CTLA4 (BPS Cat. #71149). 10 nM B7-1, Biotin-labeled (BPS Cat. #71114) and varying concentrations of CTLA4 Neutralizing Antibody (BPS Cat. #70041) were added and plate was incubated at room temperature for 2 hours. Straptavidin-HRP was added, followed by HRP substrate and chemiluminescence was read.

OUR PRODUCTS ARE FOR IN VITRO 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