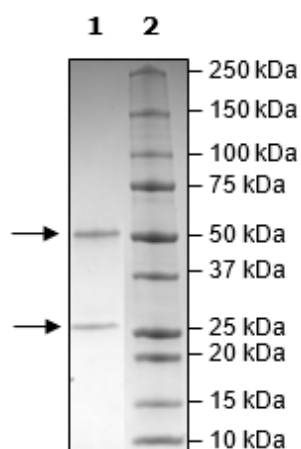


Product Information

Description:	Anti-CD137 recombinant human antibody recognizing the CD137 ligand binding region.
Concentration:	1.00 mg/ml
Species:	Human
Isotype:	IgG1
Clonality:	Monoclonal
Cross Reactivity:	This antibody not been tested for cross-reactivity with other species.
Formulated In:	8 mM phosphate, pH 7.4, 110 mM NaCl, 2.2 mM KCl, and 20% glycerol
Expression System:	HEK293
Format:	Aqueous buffer solution
Stability:	At least 12 months at -80°C. Avoid freeze/thaw cycles.
Storage:	-80°C
MW:	Light chain: 25 kDa; Heavy chain: 52 kDa
Purity:	≥90%
Purification:	Protein A affinity chromatography from HEK293 supernatants.
Applications:	This antibody is useful for activation of CD137 costimulatory pathway in T cells.

Quality Control Data

4-20% SDS-Page Coomassie Staining



CD137/NF-κB Reporter HEK293 Activity

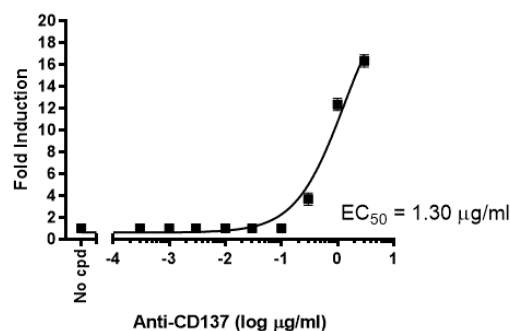


Fig. 1: Dose response of Anti-CD137 agonist antibody. CD137/NF-κB Reporter-HEK293 recombinant cells (BPS Bioscience #79289) were plated according to the cell line protocol and allowed to attach overnight. The next day a serial dilution of anti-CD137 agonist antibody was added to the cells. The plate was incubated for 24 hours before being assayed using the One-Step Luciferase Detection Reagent (BPS Bioscience #60690). Data shown is the fold induction compared to untreated wells.