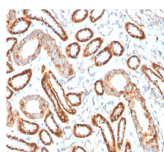


CD137 / 4-1BB / TNFRSF9 Monoclonal Mouse Antibody (4-1BB/3201)



Product Description

CD137 belongs to the tumor necrosis factor receptor family and delivers a costimulatory signal to T lymphocytes. It is expressed on activated T cells and binds an inducible ligand that is found on B cells, macrophages and dendritic cells. Interactions between CD137 and its ligand are involved in antigen presentation and the generation of cytotoxic T cells. CD137 antibody may improve cancer treatment, and has been implicated in breast cancer, melanoma and lymphoma.

Primary antibodies are available purified, or with a selection of fluorescent CF® dyes and other labels. CF® dyes offer exceptional brightness and photostability. See the [CF® Dye Brochure](#) for more information. Note: Conjugates of blue fluorescent dyes like CF®405S and CF®405M are not recommended for detecting low abundance targets, because blue dyes have lower fluorescence and can give higher non-specific background than other dye colors.

Stock status: Because Biotium offers a large number of antibody and conjugation options, primary antibody conjugates may be made to order. Typical lead times are up to one week for CF® dye and biotin conjugates, and up to 2-3 weeks for fluorescent protein and enzyme conjugates. Please email order@biotium.com to inquire about stock status and lead times before placing your order.

Catalog number key for antibody number 3201, Anti-4-1BB|CD137,|TNFRSF9 (4-1BB/3201)

Product attributes

Antibody number	#3201
Antibody reactivity (target)	4-1BB, CD137,, TNFRSF9
Antibody type	Primary
Host species	Mouse
Clonality	Monoclonal
Clone	4-1BB/3201
Isotype	IgG1, kappa
Molecular weight	32 kDa (monomer); 85 kDa (dimer)
Synonyms	4-1BB Ligand Receptor T Cell; Antigen 4-1BB Homolog; CDw137; HLDA VI; Homolog of Mouse 4 1BB; induced by lymphocyte activation (ILA); Interleukin activated receptor homolog of Mouse Ly63; T-cell antigen 4-1BB / ILA homolog; Tumor necrosis factor receptor superfamily member 9
Human gene symbol	TNFRSF9
Entrez gene ID	3604
SwissProt	Q07011
UniGene	654459; 86447
Immunogen	A recombinant fragment (around aa 19-188) of human CD137 / 4-1BB / TNFRSF9 protein (exact sequence is proprietary)
Verified antibody applications	IF (verified), IHC (FFPE) (verified), WB (verified)
Antibody target cellular localization	Plasma membrane
Species reactivity	Human
Positive control	HEK293 cells, Jurkat whole cell lysate or recombinant human CD137 protein (WB). Human kidney cancer, liver, stomach or thyroid cancer tissue (IHC-P). Expressed on the surface of activated T- cells.
Shipping condition	Room temperature
Storage Conditions	Store at 2 to 8 °C, Protect fluorescent conjugates from light, Note: store BSA-free antibodies at -10 to -35 °C
Regulatory status	For research use only (RUO)
Antibody/conjugate formulation	Conjugates: 0.1 mg/mL in PBS/0.1% BSA/0.05% azide, HRP conjugates: 0.1 mg/mL in PBS/0.05% BSA, Purified: 0.2 mg/mL in PBS/0.05% BSA/0.05% azide, Purified, BSA-free: 1 mg/mL in PBS without azide
Validated in protein array	Monospecific
Shelf life	Guaranteed for at least 24 months from date of receipt when stored as recommended
Product origin	Product may contain either bovine serum albumin (BSA) from bovine serum (<i>Bos taurus</i>), or recombinant BSA produced in Chinese hamster ovary cells. Inquire for the specific lot.
Cell/tissue expression	T-cells

Antibody # prefix	Conjugation	Ex/Em (nm)	Laser line	Detection channel	Dye Features
BNC04	CF®405S	404/431	405	DAPI (microscopy), AF405	CF®405S Features
BNC88	CF®488A	490/515	488	GFP, FITC	CF®488A Features
BNC68	CF®568	562/583	532, 561	RFP, TRITC	CF®568 Features
BNC94	CF®594	593/614	561	Texas Red®	CF®594 Features
BNC40	CF®640R	642/662	633-640	Cy®5	CF®640R Features
BNC47	CF®647	650/665	633-640	Cy®5	CF®647 Features
BNC74	CF®740	742/767	633-685	775/50	CF®740 Features
BNCB	Biotin	N/A	N/A	N/A	
BNUB	Purified	N/A	N/A	N/A	
BNUM	Purified, BSA-free	N/A	N/A	N/A	

Alexa Fluor, Pacific Blue, Pacific Orange, and Texas Red are trademarks or registered trademarks of Thermo Fisher Scientific; Cy is a registered trademark of Cytiva; IRDye, LI-COR, and Odyssey are registered trademarks of LI-COR Bioscience.

This datasheet was generated on February 12, 2025 at 01:16:08 AM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/cd137-4-1bb-tnfrsf9-monoclonal-mouse-antibody-4-1bb-3201/>