

CD257 / BAFF / TNFSF13B Monoclonal Mouse Antibody (C257/1638)

Product Description

BAFF/CD257 is a type II transmembrane protein, and a member of the tumor necrosis factor ligand superfamily. It is proteolytically cleaved to form a soluble protein. BAFF, like other TNF ligand family members, forms homo-trimers. The predicted molecular weight of a BAFF monomer is approximately 31 kDa. It is expressed on peripheral blood B and T lymphocytes, monocytes, macrophages, and dendritic cells and is upregulated by IFN-gamma and down-regulated by PMA/ionomycin treatment. BAFF stimulates B and T cell immunity and regulates humoral immunity by binding TACI and BCMA receptors; BAFF prevents apoptosis and promotes B cell survival by binding the BAFF-specific receptor (BAFFR/BR3).

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 1638, Anti-BAFF (C257/1638)

Product attributes

Antibody number	#1638
Antibody reactivity (target)	BAFF, CD257, TNFSF13B
Antibody type	Primary
Host species	Mouse
Clonality	Monoclonal
Clone	C257/1638
Isotype	IgG1
Molecular weight	31 kDa (Monomer)
Synonyms	ApoL related ligand TALL 1; B lymphocyte stimulator; B-cell-activating factor; BAFF; BLYS; CD257; Dendritic cell-derived TNF-like molecule; DTL; PRO738; TALL-1; THANK; TNFSF13B; TNFSF20; Tumor necrosis factor ligand superfamily member 13b; Tumor necrosis factor ligand superfamily member 20; ZTNF4
Human gene symbol	TNFSF13B
Entrez gene ID	10673
SwissProt	Q9Y275
Unigene	525157
Immunogen	Purified recombinant human full-length CD257 protein
Antibody target cellular localization	Plasma membrane
Species reactivity	Human
Antibody application notes	For coating for ELISA, order Ab without BSA. Higher concentration may be required for direct detection using primary antibody conjugates than for indirect detection with secondary antibody. Optimal dilution and staining procedure for a specific application should be determined by user. Recommended starting concentrations for titration are 1-2 ug/mL for most applications, or 1 ug/million cells/100 uL for flow cytometry
Positive control	Jurkat cells. Tonsil and lymph node
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
Shelf life	Guaranteed for at least 24 months from date of receipt when stored as recommended
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
Antibody research areas	Immunology
Product origin	Product may contain either bovine serum albumin (BSA) from bovine serum (Bos taurus), or recombinant BSA produced in Chinese hamster ovary cells. Inquire for the specific lot.
Cell/tissue expression	B-cells, Dendritic cells, Monocytes/macrophages, 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 April 25, 2025 at 04:59:59 AM. Visit product page to check for updated information before use. Product link: <https://biotium.com/product/mouse-monoclonal-anti-cd257-baff-tnfsf13b-c2571638/>