

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 |
| 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 |
| 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, LI-COR Bioscience.