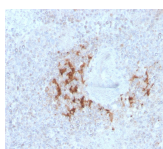


TRAF1 Monoclonal Mouse Antibody (TRAF1/3298)



Product Description

This MAbs recognizes a protein of 52 kDa, which is identified as TNFR1 (TNFR-associated factor 1). CD30-positive lymphoproliferations of the skin comprise 30% of all primary cutaneous T-cell lymphomas (CTCLs). Besides borderline cases this group includes lymphomatoid papulosis (LyP) and primary cutaneous anaplastic large T-cell lymphoma (cALCL). Although the two entities overlap clinically, histopathologically, immunopathologically and genetically, they differ considerably in their prognosis. In particular, common feature of both cases is histologically the presence of atypical lymphoid CD30-positive T blasts and genetically a clonal T-cell-receptor rearrangement. However, both cases differ considerably in their clinical course: Lesions of LyP regress spontaneously, whereas those of cALCL persist and may progress and spread. Moreover, LyP patients do not benefit from an aggressive radio- and/or chemotherapeutic approach, in contrast to patients with cALCL. Besides, LyP and cALCL differ strongly in the expression of TRAF1 (tumor necrosis factor receptor (TNFR)-associated factor 1), a component of TNFR signaling: Whereas tumor cells of most LyP cases (ca. 84%) show a strong TRAF1 expression, tumor cells of cALCL reveal TRAF1 expression in only a few cases (ca. 7%). Antibody to TRAF1 is highly useful for the differentiation of LyP and cALCL in patients with cutaneous CD30-positive lymphoproliferations.

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 3298, Anti-TRAF1 (TRAF1/3298)

Product attributes

| | |
|---------------------------------------|---|
| Antibody number | #3298 |
| Antibody reactivity (target) | TRAF1 |
| Antibody type | Primary |
| Host species | Mouse |
| Clonality | Monoclonal |
| Clone | TRAF1/3298 |
| Isotype | IgG2b, kappa |
| Molecular weight | 52 kDa |
| Synonyms | EBI6; EBV induced protein 6; Epstein-Barr virus-induced protein 6; TNF receptor-associated factor 1; TRAF1 |
| Human gene symbol | TRAF1 |
| Entrez gene ID | 7185 |
| SwissProt | Q13077 |
| Unigene | 531251 |
| Immunogen | Recombinant fragment of human TRAF1 protein (around aa 73-219) (exact sequence is proprietary) |
| Verified antibody applications | IHC (FFPE) (verified) |
| Antibody target cellular localization | Cytoplasm, Nucleus |
| Species reactivity | Human |
| Positive control | HeLa or 293T cells. Hodgkin s Lymphoma (IHC). |
| 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 (Bos taurus), or recombinant BSA produced in Chinese hamster ovary cells. Inquire for the specific lot. |
| Tumor expression | Leukemia/lymphoma |

| 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 August 7, 2025 at 07:26:14 PM. Visit product page to check for updated information before use. Product link: <https://biotium.com/product/traf1-monoclonal-mouse-antibody-traf1-3298/>