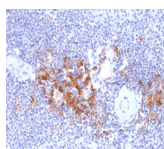


# TRAF1 Monoclonal Mouse Antibody (TRAF1/2770)



## 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](mailto:order@biotium.com) to inquire about stock status and lead times before placing your order.

**Catalog number key for antibody number 2770, Anti-TRAF1 (TRAF1/2770)**

## Product attributes

Antibody number	#2770
Antibody reactivity (target)	TRAF1
Antibody type	Primary
Host species	Mouse
Clonality	Monoclonal
Clone	TRAF1/2770
Isotype	IgG1, 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	<a href="#">CF®405S Features</a>
BNC88	CF®488A	490/515	488	GFP, FITC	<a href="#">CF®488A Features</a>
BNC68	CF®568	562/583	532, 561	RFP, TRITC	<a href="#">CF®568 Features</a>
BNC94	CF®594	593/614	561	Texas Red®	<a href="#">CF®594 Features</a>
BNC40	CF®640R	642/662	633-640	Cy®5	<a href="#">CF®640R Features</a>
BNC47	CF®647	650/665	633-640	Cy®5	<a href="#">CF®647 Features</a>
BNC74	CF®740	742/767	633-685	775/50	<a href="#">CF®740 Features</a>
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 July 18, 2025 at 12:43:09 PM. Visit product page to check for updated information before use. Product link: <https://biotium.com/product/traf1-monoclonal-mouse-antibody-traf1-2770/>