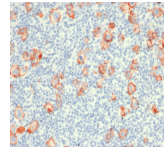


CD30 / TNFRSF8 Recombinant Monoclonal Mouse Antibody (rCD30/412)



Product Description

This antibody recognizes a single chain glycoprotein of 105/120 kDa, identified as CD30/Ki-1. CD30 is synthesized as a 90 kDa precursor, which is processed in the Golgi complex into a membrane-bound phosphorylated mature 105/120 kDa glycoprotein. In Hodgkin's disease, CD30/Ki-1 antigen is expressed by mononuclear-Hodgkin and multinucleated Reed-Sternberg cells. It is also expressed by the tumor cells of a majority of anaplastic large cell lymphomas as well as by a varying proportion of activated T and B cells. This MAb distinguishes large cell lymphomas derived from activated lymphoid cells from histiocytic malignancies and lymphomas derived from resting and precursor lymphoid cells or from anaplastic carcinomas. About one third of the Ki-1 positive lymphomas lack the leukocyte common antigen (CD45). 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 1919, Anti-CD30 / TNFRSF8 (rCD30/412)**

Product attributes

Antibody number	#1919
Antibody reactivity (target)	CD30, TNFRSF8
Antibody type	Primary
Host species	Mouse
Clonality	Recombinant Monoclonal
Clone	rCD30/412
Isotype	IgG2a, kappa
Molecular weight	105-120 kDa
Synonyms	CD30L receptor; Cytokine receptor CD30; Ki-1 antigen; Lymphocyte activation antigen CD30; Tumor necrosis factor receptor superfamily member 8 (TNFRSF8)
Human gene symbol	TNFRSF8
Entrez gene ID	943
SwissProt	P28908
Unigene	1314
Immunogen	Recombinant full-length human TNFRSF8 protein
Verified antibody applications	IHC (FFPE) (verified)
Antibody target cellular localization	Plasma membrane
Species reactivity	Human
Positive control	Hodgkin's lymphoma
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	Lymphocytes
Tumor expression	Leukemia/lymphoma

Antibody # prefix	Conjugation	Ex/Em (nm)	Laser line	Detection channel
BNC04	CF@405S	404/431	405	DAPI (microscopy), AF405
BNC88	CF@488A	490/515	488	GFP, FITC
BNC68	CF@568	562/583	532, 561	RFP, TRITC
BNC94	CF@594	593/614	561	Texas Red [®]
BNC40	CF@640R	642/662	633-640	Cy@5
BNC47	CF@647	650/665	633-640	Cy@5
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

Dye Features

- [CF@405S Features](#)
- [CF@488A Features](#)
- [CF@568 Features](#)
- [CF@594 Features](#)
- [CF@640R Features](#)
- [CF@647 Features](#)

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.