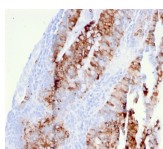


CD262 / DR5 Monoclonal Mouse Antibody (DR5/3381)



Product Description

Tumor necrosis factor (TNF) is a pleiotropic cytokine whose function is mediated by two distinct cell surface receptors, designated TNF-R1 and TNF-R2, which are expressed on most cell types. TNF function is primarily mediated through TNF-R1 signaling. Both receptors belong to the growing TNF receptor superfamily which includes Fas antigen and CD40. TNF-R1 contains a cytoplasmic motif, termed the death domain, that has been found to be necessary for the transduction of the apoptotic signal. The death domain is also found in several other receptors, including Fas, DR2 (or TRUNDD), DR3 (death receptor 3), DR4 and DR5. TRUNDD, DR4 and DR5 are receptors for the apoptosis-inducing cytokine TRAIL. A non-death domain-containing receptor, designated decoy receptor (DcR1 or TRID), also specifically associates with TRAIL and may play a role in cellular resistance to apoptotic stimuli.

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 3381, Anti-CD262|DR5 (DR5/3381)

Call us : [800-304-5357](tel:800-304-5357)

Product attributes

Antibody number	#3381
Antibody reactivity (target)	CD262, DR5
Antibody type	Primary
Host species	Mouse
Clonality	Monoclonal
Clone	DR5/3381
Isotype	IgG1, kappa
Molecular weight of antigen	48 kDa
Synonyms	Fas like protein; Apoptosis inducing protein TRICK2A/2B; Apoptosis inducing receptor TRAIL R2; Cytotoxic TRAIL receptor 2; Death domain containing receptor for TRAIL/Apo 2L; Death receptor 5 (DR5); p53 regulated DNA damage inducible cell death receptor (killer); TNF related apoptosis inducing ligand receptor 2 (TRICK2; Tumor necrosis factor receptor like protein ZTNFR9; Tumor necrosis factor receptor superfamily member 10B; ZTNFR9
Human gene symbol	TNFRSF10B
Entrez gene ID	8795
SwissProt	521456
Unigene	661668
Immunogen	Recombinant human DR5 protein fragment (around aa266-393) (exact sequence is proprietary)
Verified antibody applications	IHC (FFPE) (verified)
Antibody target cellular localization	Plasma membrane
Species reactivity	Human
Positive control	Human colon cancer or endometrial tissue (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 (<i>Bos taurus</i>), or recombinant BSA produced in Chinese hamster ovary cells. Inquire for the specific lot.

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 June 21, 2026 at 10:48:00 AM. Visit product page to check for updated information before use.
 Product link: <https://biotium.com/product/cd262-dr5-monoclonal-mouse-antibody-dr5-3381/>