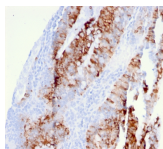


CD262 / DR5 Monoclonal Mouse Antibody (DR5/3381)



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 | 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 (Bos taurus), 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 December 11, 2025 at 10:24:46 PM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/cd262-dr5-monoclonal-mouse-antibody-dr5-3381/>