



Glowing products for science

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

CDw78 Monoclonal Mouse Antibody (DF1588)

Product Description

This antibody reacts with a carbohydrate determinant, identified as CDw78 (Workshop V; Code CDw78.2). Epitope recognized by this MAb and Leu21 is the same. It is present on some immature and some peripheral blood B cells, and some B cell lines. Its expression is increased after B-cell activation in vitro. CDw78 is also found on tissue macrophages and on epithelial cells, but not on T cells, NK cells, monocytes, granulocytes and myeloid tissues. It appears that CDw78 is a determinant on a major histocompatibility complex (MHC) class II subpopulation. **Catalog number key for antibody number 0353, Anti-CDw78 (DF1588)**

Product attributes

| | |
|---------------------------------------|---|
| Antibody number | #0353 |
| Antibody reactivity (target) | CDw78 |
| Antibody type | Primary |
| Host species | Mouse |
| Clonality | Monoclonal |
| Clone | DF1588 |
| Isotype | IgG1, kappa |
| Molecular weight | Not Known |
| Synonyms | Leu21; Ba antigen; LO panBa |
| Human gene symbol | Not Known |
| Entrez gene ID | Not Known |
| SwissProt | Not Known |
| Unigene | Not Known |
| Immunogen | Human leucocytes |
| Antibody target cellular localization | Plasma membrane |
| Species reactivity | Human |
| Antibody application notes | For coating for ELISA, order Ab without BSA, Higher concentration may be required for direct detection using primary antibody conjugates than for indirect detection with secondary antibody, Optimal dilution and staining procedure for a specific application should be determined by user, Recommended starting concentrations for titration are 1-2 ug/mL for most applications, or 1 ug/million cells/100 uL for flow cytometry |
| Positive control | Raji and human peripheral blood lymphocytes. Human tonsil and lymph node. |
| 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 |
| Shelf life | Guaranteed for at least 24 months from date of receipt when stored as recommended |
| Cell/tissue expression | Macrophage/monocyte, B-cells, Epithelial cells |
| 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 research areas | Immunology |

| 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 12, 2025 at 01:49:28 AM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/cdw78-monoclonal-mouse-antibody-df1588/>