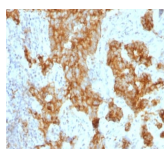


# CD11c Monoclonal Mouse Antibody (ITGAX/1243)



## Product Description

This antibody recognizes a protein of 145 kDa, identified as CD11c. CD11c (ITGAX), a member of the leukointegrin family, shares the same beta subunit with other members of the leukocyte adhesion molecule family, which includes CD11a (LFA-1), CD11b (MAC-1) and CD11d (ITGAD), but has a unique alpha chain. CD11c has been shown to play a role in phagocytosis, cell migration, and cytokine production by monocytes/macrophages as well as induction of T-cell proliferation by Langerhans cells. CD11c is expressed prominently on the plasma membranes of monocytes, tissue macrophages, NK cells, and most dendritic cells (DCs). A lower level of expression is also observed on neutrophils as a result of its high level of expression on most DCs. An antibody to CD11c may aid in identification of lesions with histiocytic origin. It may also been used as a marker for hairy cell leukemia in paraffin-embedded tissues.

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 1243, Anti-CD11c (ITGAX/1243)**

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

## Product attributes

|                                       |   |
|---------------------------------------|---|
| Antibody number                       | #1243   |
| Antibody reactivity (target)          | CD11c   |
| Antibody type                         | Primary   |
| Host species                          | Mouse   |
| Clonality                             | Monoclonal  |
| Clone                                 | ITGAX/1243  |
| Isotype                               | IgG2b, kappa  |
| Molecular weight of antigen           | 145 kDa   |
| Synonyms                              | CD11 antigen-like family member C; Complement component 3 receptor 4 subunit; Integrin alpha-X; integrin, alpha X (antigen CD11C (p150), alpha polypeptide); Leu M5 alpha subunit; Leukocyte adhesion glycoprotein p150 95 alpha chain; Myeloid membrane antigen alpha subunit; p150/95   |
| Human gene symbol                     | ITGAX   |
| Entrez gene ID                        | 3687  |
| SwissProt                             | P20702  |
| Unigene                               | 248472  |
| Immunogen                             | Recombinant human ITGAX protein fragment (aa 637-827) (exact sequence is proprietary)   |
| 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                      | THP-1 cells. Tonsil or 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   |
| 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   |
| Validated in protein array            | Monospecific  |
| Antibody research areas               | Cell adhesion, Immunology   |
| 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.  |
| 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 June 13, 2026 at 08:04:00 AM. Visit product page to check for updated information before use. Product link: <https://biotium.com/product/cd11c-monoclonal-mouse-antibody-itgax-1243/>