

CD11a Monoclonal Mouse Antibody (DF1524)

Product Description

Recognizes a protein of 180 kDa, identified as CD11a (Leucocyte Workshop IV; Code 1524). CD11a complex with the 2 subunit of the integrin family, CD18, to form the cell surface heterodimer, LFA-1 or CD11a /C18 (aLbL). LFA-1 is expressed on all leukocytes including lymphocytes, monocytes, and granulocytes. It is involved in leukocyte adhesion to its ligands including intercellular adhesion molecule-1 (ICAM-1 or CD54), ICAM-2 (CD102), ICAM-3 (CD50) and Telencephalin (TLN) and play a role in most immune/inflammatory responses. This MAb potently blocks LFA-1 dependent homotypic cell aggregation. 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 0312, Anti-CD11a (DF1524)**

References

Note: References for this clone sold by other suppliers may be listed for expected applications.

1. J Neuropathol Exp Neurol (2000) 59(1): 18-28. (functional studies)
2. FASEB J (2003) 17(11), 1-21. (Flow)
3. Blood (2006) 107(8): 3212-3220. (IF)

Product attributes

| | |
|---------------------------------------|--|
| Antibody number | #0312 |
| Antibody reactivity (target) | CD11a |
| Antibody type | Primary |
| Host species | Mouse |
| Clonality | Monoclonal |
| Clone | DF1524 |
| Isotype | IgG2b, kappa |
| Molecular weight | 180 kDa |
| Synonyms | Antigen CD11a (p180), lymphocyte function associated antigen 1, alpha polypeptide; Integrin alpha-L; ITGAL; Leukocyte Adhesion Glycoprotein LFA1 Alpha Chain; LFA 1 alpha (LFA1A); Ly15; Ly21; p180 |
| Human gene symbol | ITGAL |
| Entrez gene ID | 3683 |
| SwissProt | P20701 |
| Unigene | 174103 |
| Immunogen | Human CD11a protein |
| Antibody target cellular localization | Plasma membrane |
| Expected antibody applications | Flow, surface (published for clone), Functional studies (published for clone), IF (published for clone) |
| Species reactivity | Human |
| Antibody application notes | Higher concentration may be required for direct detection using primary antibody conjugates than for indirect detection with secondary antibody, Immunofluorescence: 0.5-1 ug/mL, Flow Cytometry 0.5-1 ug/million cells/0.1 mL |
| Positive control | JY, HUT-78, MOLT-4, HL60, U937, and human leukocytes. Human lymph nodes and tonsils. |
| 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 | Hematopoietic cells |

| 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 |
| 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, LI-COR Bioscience.