## CD6 Monoclonal Mouse Antibody (C6/372 + 3F7B5)



## **Product Description**

CD6 is a type I transmembrane glycoprotein that contains a 24-amino acid signal sequence, three extracellular scavenger receptor cysteine-rich (SRCR) domains, a membrane-spanning domain and a 44-amino acid cytoplasmic domain. The CD6 glycoprotein is tyrosine phosphorylated during TCR-mediated T cell activation. CD6 shows significant homology to CD5. CD6 is present on mature thymocytes, peripheral T cells and a subset of B cells. Antibodies to CD6 are used to deplete T cells from bone marrow transplants to prevent graft versus host disease.

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 <a href="mailto:order@biotium.com">order@biotium.com</a> to inquire about stock status and lead times before placing your order.

Catalog number key for antibody number 1228, Anti-CD6 (C6/372 3F7B5)

## Droduct attributes

| Product attributes                    |   |  |  |
|---------------------------------------|---|--|--|
| Antibody number                       | #1228   |  |  |
| Antibody reactivity (target)          | CD6   |  |  |
| Antibody type                         | Primary   |  |  |
| Host species                          | Mouse   |  |  |
| Clonality                             | Monoclonal  |  |  |
| Clone                                 | C6/372 + 3F7B5  |  |  |
| Isotype                               | IgG's   |  |  |
| Molecular weight                      | 90-130 kDa  |  |  |
| Synonyms                              | T12; TP120  |  |  |
| Human gene symbol                     | CD6   |  |  |
| Entrez gene ID                        | 923   |  |  |
| SwissProt                             | P30203  |  |  |
| Unigene                               | 744366  |  |  |
| Immunogen                             | Human recombinant CD6 protein (C6/372); Human rheumatoid synovial T cell line ST-1 (3F7B5)  |  |  |
| Antibody target cellular localization | Plasma membrane   |  |  |
| Verified antibody applications        | IF (verified), IHC (FFPE) (verified)  |  |  |
| 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, Immunohistology formalin-fixed 0.5-1 ug/mL, Staining of formalin-fixed tissues requires boiling tissue sections in 10 mM citrate buffer, pH 6.0, for 10-20 min followed by cooling at RT for 20 minutes, Flow Cytometry 0.5-1 ug/million cells/0.1 mL, Optimal dilution for a specific application should be determined by user |  |  |
| Positive control                      | CCRF-CEM, Jurkat cells, Tonsil  |  |  |
| 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% BSA-free: 1 mg/mL in PBS without azide  |  |  |
| Cell/tissue expression                | B-cells, T-cells  |  |  |
| Antibody research areas               | Immunology  |  |  |
| 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.   |  |  |

Call us: 800-304-5357 Email: btinfo@biotium.com

| Antibody # prefix | Conjugation           | Ex/Em (nm) | Laser line | Detection channel           | Dye Features     |
|-------------------|-----------------------|------------|------------|-----------------------------|------------------|
| BNC04             | CF®405S               | 404/431    | 405        | DAPI (microscopy),<br>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,<br>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 November 4, 2025 at 07:02:12 AM. Visit product page to check for updated information before use. Product link: <a href="https://biotium.com/product/monoclonal-anti-cd6-c6372-3f7b5/">https://biotium.com/product/monoclonal-anti-cd6-c6372-3f7b5/</a>