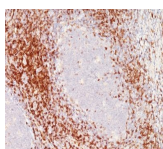


# CD6 Monoclonal Mouse Antibody (C6/372 + 3F7B5)



## 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 [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 1228, Anti-CD6 (C6/372 3F7B5)**

## 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% azide, Purified, 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.                                                                                                                                                                                                                                                                                                                                   |

| 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 December 23, 2024 at 03:55:30 PM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/monoclonal-anti-cd6-c6372-3f7b5/>