

## **CD27 Monoclonal Mouse Antibody** (LPFS2/1611)

## **Product Description**

This antibody recognizes a protein of a disulfide-linked 120 kDa dimer, identified as CD27. It is expressed on the majority of peripheral T cells, medullary thymocytes, memory-type B cells, and natural killer cells. It is a transmembrane phosphoglycoprotein that belongs to the tumor necrosis factor receptor (TNFR) superfamily. CD27 binds to its ligand CD70, a member of the TNF family, and induces T-cell co-stimulation and B-cell activation. It also interacts with TRAFs and is involved in activation of NFkB and SAPK/JNK and induces apoptosis.

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 1611, Anti-CD27 (LPFS2/1611)

## Product attributes

Call us: 800-304-5357

| Product attributes                    |  |  |  |
|---------------------------------------|--|--|--|
| Antibody number                       | #1611  |  |  |
| Antibody reactivity (target)          | CD27   |  |  |
| Antibody type                         | Primary  |  |  |
| Host species                          | Mouse  |  |  |
| Clonality                             | Monoclonal   |  |  |
| Clone                                 | LPFS2/1611   |  |  |
| Isotype                               | IgG1   |  |  |
| Molecular weight                      | 120 kDa  |  |  |
| Synonyms                              | LPFS2; S152; T cell activation antigen S152; T-cell activation antigen CD27; T14; TNFRSF7; TNFSF7; Tp55; Tumor necrosis factor receptor superfamily member 7   |  |  |
| Human gene symbol                     | CD27   |  |  |
| Entrez gene ID                        | 939  |  |  |
| SwissProt                             | P26842   |  |  |
| Unigene                               | 355307   |  |  |
| Immunogen                             | Recombinant human CD22 protein (exact sequence is proprietary)   |  |  |
| Antibody target cellular localization | Plasma membrane  |  |  |
| Verified antibody applications        | Flow (verified), 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): 0.5-1 ug/mL, Staining of formalin-fixed tissues requires boiling tissue sections in 10 mM Tris buffer with 1 mM EDTA pH 9.0 for 10-20 min followed by cooling at RT for 20 min, Flow Cytometry 0.5-1 ug/million cells/0.1 mL, Optimal dilution for a specific application should be determined by user |  |  |
| Positive control                      | Human peripheral blood lymphocytes. Human tonsil and lymph node.   |  |  |
| Shipping condition                    | Room temperature   |  |  |
| Storage Conditions                    | itions Store at 2 to 8 °C, Protect fluorescent conjugates from light,<br>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   |  |  |
| 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.  |  |  |
| Cell/tissue expression                | Lymphocytes  |  |  |
| •                                     | Product may contain either bovine serum albumin (BSA) from bovine serum (Bos taurus), or recombinant BSA produced in   |  |  |
| Cell/tissue expression                | Lymphocytes  |  |  |

| 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 30, 2025 at 03:28:10 AM. Visit product page to check for updated information before use. Product link: <a href="https://biotium.com/product/mouse-monoclonal-anti-cd27-lpfs21611/">https://biotium.com/product/mouse-monoclonal-anti-cd27-lpfs21611/</a>