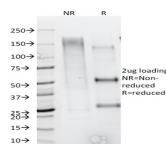


CD79b Monoclonal Mouse Antibody (IGB/1844)



Product Description

CD79 (also designated Ig α /Ig β) is a heterodimer composed of α chains, designated CD79A or MB-1, and β chains, designated CD79B or B29. The B cell antigen receptor complex (BCR) is formed by the association of CD79 with a membrane immunoglobulin, such as IgM or IgD. The membrane immunoglobulins IgM and IgD achieve surface expression and antigen presentation function in response to CD79 association. The cytoplasmic tails of both CD79A and CD79B contain an ITAM (immuno-receptor tyrosine-based activation) motif, which acts to initiate the BCR signaling reactions by binding to and activating tyrosine kinases.

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 1844, Anti-CD79b (IGB/1844)

Product attributes

| | |
|--|---|
| Antibody number | #1844 |
| Antibody reactivity (target) | CD79b |
| Antibody type | Primary |
| Host species | Mouse |
| Clonality | Monoclonal |
| Clone | IGB/1844 |
| Isotype | IgG1, kappa |
| Molecular weight of antigen | 39 kDa |
| Synonyms | AGM6; B-cell antigen receptor complex-associated protein beta chain; B-cell-specific glycoprotein B29; B29; B29/Ig-beta/CD79b; CD79b; Ig-beta; IGB; Immunoglobulin-associated B29 protein |
| Human gene symbol | CD79B |
| Entrez gene ID | 974 |
| SwissProt | P40259 |
| Unigene | 89575 |
| Immunogen | Recombinant human CD79b protein fragment (around aa 29-159) (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 | Daudi or Ramos cells. Germinal center B-cells in a lymph node or 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 |
| Validated in protein array | Monospecific |
| Antibody research areas | 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. |
| Cell/tissue expression | NK cells, T-cells |
| 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 | 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, 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 22, 2026 at 03:48:53 PM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/cd79b-monoclonal-mouse-antibody-igb-1844/>