



Glowing products for science

CD19 Monoclonal Mouse Antibody (SJ25C1) - Biotium Choice

CD19 (clone SJ25C1) is a validated mouse monoclonal antibody that recognizes CD19, a transmembrane glycoprotein in B cells. The antibody belongs to the Biotium Choice list of select antibodies that have been validated and optimized in-house for optimal performance.



Product Description

CD19 (clone SJ25C1) is a mouse monoclonal antibody that recognizes the transmembrane glycoprotein CD19. This antibody belongs to the Biotium Choice list of select antibodies that have been validated and optimized in-house for optimal performance. The antibody is available conjugated to CF® Dyes, Astral™ tandem dyes, R-PE and PerCP. They are supplied in PBS, 0.1% BSA, 0.05% azide.

CD19 is a transmembrane glycoprotein that contains two extracellular immunoglobulin-like domains. CD19 is present in both benign and malignant B-cells and is considered to be the most reliable surface marker of this lineage over a wide range of maturational stages. In normal lymphoid tissue, CD19 is observed in germinal centers, in mantle zone cells, and in scattered cells of the inter-follicular areas. Anti-CD19 exhibits an overall immunoreactivity pattern similar to those of the antibodies against CD20 and CD22. However, in contrast to CD20, expression of CD19 is continuous throughout B-cell development and through terminal differentiation of B-cells into plasma cells. Anti-CD19 positivity is seen in the vast majority of B-cell neoplasms commonly at a lower intensity than normal B-cell counterparts. Plasma cell neoplasms are nearly always negative, as are T-cell neoplasms.

As an isotype control, we recommend [Isotype Control, Monoclonal Mouse IgG1 Kappa \(IGG1/1331\)](#) for this antibody.

Discover Sharper Signals and Unmatched Panel Flexibility with Biotium Choice Antibodies Powered by CF® Dyes and Astral Leap™ Tandem Dyes

Biotium Choice antibodies are carefully curated and validated in-house to offer exceptional signal-to-noise. Labeled with our advanced CF® Dyes and Astral Leap™ tandem dyes, they are our top-recommended antibodies for flow cytometry and other applications.

Biotium Choice Antibody Features

- Robust and validated clones against common targets
- Developed and optimized for flow cytometry and other applications
- Conjugated to bright, photostable [CF® Dyes](#) for superior signal and clarity
- Also available with [Astral Leap™ tandem dyes](#) for expanding multiplexing
- New antibody clones and dye conjugates continuously in development

[View our full catalog of Biotium Choice antibodies](#)

Call us : [800-304-5357](tel:800-304-5357)

Product attributes

| | |
|---------------------------------------|---|
| Antibody number | #P010 |
| Antibody reactivity (target) | CD19 |
| Biotium Choice Antibody | Primary |
| Antibody type | Biotium Choice Flow Antibodies, Primary |
| Host species | Mouse |
| Clonality | Monoclonal |
| Clone | SJ25C1 |
| Isotype | IgG1, kappa |
| Molecular weight | 95 kDa |
| Synonyms | B-Lymphocytes Surface Antigen B4; T-cell Surface Antigen Leu-12; CVID3; B4 |
| Human gene symbol | CD19 |
| Entrez gene ID | 930 |
| SwissProt | P15391 |
| Unigene | 652262 |
| Antibody target cellular localization | Plasma membrane |
| Verified antibody applications | Flow (surface) (verified) |
| Species reactivity | Human |
| Antibody application notes | Recommended concentration for flow cytometry: 5 uL per million cells/0.1 mL |
| Positive control | Human peripheral blood mononuclear cells (PBMCs) |
| Shipping condition | Room temperature |
| Storage Conditions | Store at 2 to 8 °C, Protect fluorescent conjugates from light |
| 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: PBS/0.1% BSA/0.05% azide |
| Antibody research areas | Immunology |
| Cell/tissue expression | B-cells, Lymphocytes |
| Volume per assay | 5 uL/test |
| 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. |

CD19 Monoclonal Mouse Antibody (SJ25C1) - Biotium Choice

| Conjugation | Ex/Em | Conc. | Size | Catalog No. | Dye Features |
|--|---|-----------|-------------------|-------------------------------|---|
| CF@405M 100 tests (500 uL) | 416/452 nm P010-405M-500 | 100 ug/mL | 25 tests (125 uL) | P010-405M-125 | CF@405M Features |
| CF@488A 100 tests (500 uL) | 490/516 nm P010-488A-500 | 100 ug/mL | 25 tests (125 uL) | P010-488A-125 | CF@488A Features |
| CF@575 100 tests (500 uL) | 575/601 nm P010-575-500 | 200 ug/mL | 25 tests (125 uL) | P010-575-125 | |
| RPE-Astral™616 100 tests (500 uL) | 496, 546, 566/617 nm P010-R616-500 | 50 ug/mL | 25 tests (125 uL) | P010-R616-125 | RPE-Astral™616 Features |
| CF@647Plus 100 tests (500 uL) | 652/668 nm P010-647P-500 | 25 ug/mL | 25 tests (125 uL) | P010-647P-125 | |
| CF@700 100 tests (500 uL) | 696/721 nm P010-700-500 | 100 ug/mL | 25 tests (125 uL) | P010-700-125 | CF@700 Features |
| APC-Astral™813 100 tests (500 uL) | 633, 638/813 nm P010-A813-500 | 50 ug/mL | 25 tests (125 uL) | P010-A813-125 | APC-Astral™813 Features |
| R-PE 100 tests (500 uL) | 496, 546, 565/576 nm P010-RPE-500 | 50 ug/mL | 25 tests (125 uL) | P010-RPE-125 | |
| PerCP 100 tests (500 uL) | 482/675 nm P010-PCP-500 | 100 ug/mL | 25 tests (125 uL) | P010-PCP-125 | |

This datasheet was generated on December 14, 2025 at 11:29:58 PM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/cd19-monoclonal-mouse-antibody-sj25c1/>