Bcl-6 Monoclonal Mouse Antibody (BCL6/1718)



Product Description

Bcl-6 is a transcriptional repressor mainly required for germinal center (GC) formation and antibody affinity maturation which has different mechanisms of action specific to the lineage and biological functions. It forms complexes with different corepressors and histone deacetylases to repress the transcriptional expression of different subsets of target genes. Bcl-6 allows GC B-cells to proliferate very rapidly in response to T-cell dependent antigens and tolerate the physiological DNA breaks required for immunglobulin class switch recombination and somatic hypermutation without inducing a p53/TP53-dependent apoptotic response. Antibodies to Bcl-6 are helpful in a number of diagnostic settings, for example in the differential diagnosis of small B-cell lymphoma, and in distinguishing classical Hodgkin lymphoma from nodular lymphocyte predominant Hodgkin lymphoma (NLPHL). 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®-058 and CF®-05M 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 1718, Anti-Bcl-6 (BCL6/1718)

Product attributes

| Product attributes | | | |
|---------------------------------------|---|--|--|
| Antibody number | #1718 | | |
| Antibody reactivity (target) | Bcl-6 | | |
| Antibody type | Primary | | |
| Host species | Mouse | | |
| Clonality | Monoclonal | | |
| Clone | BCL6/1718 | | |
| Isotype | IgG1, kappa | | |
| Molecular weight | 95 kDa | | |
| Synonyms | Bcl-6; B-Cell Lymphoma 6 Protein; B-cell lymphoma 5 protein; Zinc Finger Protein 51; zinc finger transcription factor BCL6S; BCL6 Transcription Repressor | | |
| Human gene symbol | BCL6 | | |
| Entrez gene ID | 604 | | |
| SwissProt | P41182 | | |
| Unigene | 478588 | | |
| Immunogen | Recombinant human bcl-6 protein fragment (aa256-389) (exact sequence is proprietary) | | |
| Antibody target cellular localization | Nucleus | | |
| Verified antibody applications | IHC (FFPE) (verified) | | |
| Species reactivity | Human | | |
| Positive control | Raji or Ramos cells. Tonsil or Hodgkin's lymphoma. | | |
| 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 | Apoptosis, Cancer, Immunology, Transcription factor | | |
| 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. | | |
| Tumor expression | Leukemia/lymphoma | | |
| | | | |

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), 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 November 2, 2025 at 02:49:26 PM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/bcl-6-monoclonal-mouse-antibody-bcl6-1718/