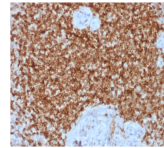


## Bcl-2 Recombinant Monoclonal Rabbit Antibody (BCL2/1878R)



### Product Description

This antibody recognizes a protein of 25-26 kDa, identified as Bcl-2. This antibody shows no cross-reaction with Bcl-x or Bax proteins. Bcl-2 suppresses apoptosis in a variety of cell systems including factor-dependent lymphohematopoietic and neural cells. It regulates cell death by controlling the mitochondrial membrane permeability. Bcl-2 appears to function in a feedback loop system with caspases. It inhibits caspase activity either by preventing the release of cytochrome c from the mitochondria and/or by binding to the apoptosis-activating factor (APAF-1). Expression of Bcl-2 inhibits programmed cell death (apoptosis). In most follicular lymphomas, neoplastic germinal centers express high levels of Bcl-2 protein, whereas the normal or hyperplastic germinal centers are negative. Consequently, this antibody is valuable when distinguishing between reactive and neoplastic follicular proliferation in lymph node biopsies. It may also be used in distinguishing between those follicular lymphomas that express Bcl-2 protein and the small number in which the neoplastic cells are Bcl-2 negative.

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 1878, Anti-Bcl-2 (BCL2/1878R)**

### Product attributes

|                                       |                                                                                                                                                                                                   |
|---------------------------------------|---------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------|
| Antibody number                       | #1878                                                                                                                                                                                             |
| Antibody reactivity (target)          | Bcl-2                                                                                                                                                                                             |
| Antibody type                         | Primary                                                                                                                                                                                           |
| Host species                          | Rabbit                                                                                                                                                                                            |
| Clonality                             | Recombinant Monoclonal                                                                                                                                                                            |
| Clone                                 | BCL2/1878R                                                                                                                                                                                        |
| Isotype                               | IgG                                                                                                                                                                                               |
| Molecular weight                      | 25-26 kDa                                                                                                                                                                                         |
| Synonyms                              | Apoptosis regulator Bcl-2; B-cell CLL/lymphoma-2                                                                                                                                                  |
| Human gene symbol                     | BCL2                                                                                                                                                                                              |
| Entrez gene ID                        | 596                                                                                                                                                                                               |
| SwissProt                             | P10415                                                                                                                                                                                            |
| Unigene                               | 150749                                                                                                                                                                                            |
| Immunogen                             | Recombinant full-length human bcl-2 protein                                                                                                                                                       |
| Verified antibody applications        | IHC (FFPE) (verified), WB (verified)                                                                                                                                                              |
| Antibody target cellular localization | Endoplasmic reticulum, Mitochondria, Nuclear membrane                                                                                                                                             |
| Species reactivity                    | Human                                                                                                                                                                                             |
| Positive control                      | Jurkat, K562, HL-60, or HeLa Cells. Tonsil or follicular lymphomas.                                                                                                                               |
| 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 |
| Antibody research areas               | Apoptosis, Cancer                                                                                                                                                                                 |

| 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>  |
| 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, LI-COR Bioscience.