## **BOB-1 Monoclonal Mouse Antibody** (BOB1/2423)



## **Product Description**

BOB-1 is a transcriptional coactivator that specifically associates with either POU2F1/OCT1 or POU2F2/OCT2, boosting the promoter activity of those complexes. It is essential for the response of B-cells to antigens, required for the formation of germinal centers, and regilates IL-6 expression. BOB-1 expression in a variety of established B-cell lines, representing different stages of B-cell development, has suggested a constitutive, B-cell-specific expression pattern. LP cells in nodular lymphocyte predominant Hodgkin lymphoma, because they are germinal center-derived, are consistently immuno-positive for BOB-1. Conversely, only some cases of classical Hodgkin lymphoma show BOB-1 immuno-reactivity within the Hodgkin and Reed-Sternberg cells. Expression of BOB-1 has been reported in follicular center cell lymphoma, diffuse large B-cell lymphoma and some cases of acute myeloid leukemia. B-CLL, marginal zone lymphoma, and mantle cell lymphoma may show weak to moderate immunoreactivity.

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 2423, Anti-BOB.1 (BOB1/2423)

## Product attributes

| Product attributes                      |  |  |  |
|---|--|--|--|
| Antibody number                         | #2423  |  |  |
| Antibody reactivity (target)            | BOB.1  |  |  |
| Antibody type                           | Primary  |  |  |
| Host species                            | Mouse  |  |  |
| Clonality                               | Monoclonal   |  |  |
| Clone                                   | BOB1/2423  |  |  |
| sotype                                  | IgG2b, kappa   |  |  |
| Molecular weight                        | 35 kDa   |  |  |
| Synonyms                                | BOB-1; B-cell-specific coactivator OBF-1; OCT-binding factor 1; POU domain class 2-associating factor 1  |  |  |
| Human gene symbol                       | POU2AF1  |  |  |
| Entrez gene ID                          | 5450   |  |  |
| SwissProt                               | Q16633   |  |  |
| Unigene                                 | 654525   |  |  |
| mmunogen                                | Recombinant fragment (around aa 148-255) of human BOB1 (POU2AF1)protein (exact sequence is proprietary)  |  |  |
| Verified antibody<br>applications       | IHC (FFPE) (verified)  |  |  |
| Antibody target cellular<br>ocalization | Nucleus  |  |  |
| 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, immunohistology (formalin): 1-2 ug/mL for 30 minutes at RT, Staining of formalin-fixed tissues requires boiling tissue sections in 10 mM citrate buffer, pH 6.0, for 10-20 minutes followed by cooling at RT for 20 minutes, Optimal dilution for a specific application should be determined by user |  |  |
| Positive control                        | Ramos cells. Spleen 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<br>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                 | Cancer, 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                  | B-cells  |  |  |
| 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),<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 3, 2025 at 09:34:24 PM. Visit product page to check for updated information before use. Product link: <a href="https://biotium.com/product/bob-1-monoclonal-mouse-antibody-bob1-2423/">https://biotium.com/product/bob-1-monoclonal-mouse-antibody-bob1-2423/</a>