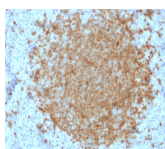


# BOB-1 Monoclonal Mouse Antibody (BOB1/2422)



## 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 regulates 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 [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 2422, BOB.1 Monoclonal Mouse Antibody (BOB1/2422)**

## Product attributes

<b>Antibody number</b>	#2422
<b>Antibody reactivity (target)</b>	BOB.1
<b>Antibody type</b>	Primary
<b>Host species</b>	Mouse
<b>Clonality</b>	Monoclonal
<b>Clone</b>	BOB1/2422
<b>Isotype</b>	IgG2b, kappa
<b>Molecular weight of antigen</b>	35 kDa
<b>Synonyms</b>	BOB-1; B-cell-specific coactivator OBF-1; OCT-binding factor 1; POU domain class 2-associating factor 1
<b>Human gene symbol</b>	POU2AF1
<b>Entrez gene ID</b>	5450
<b>SwissProt</b>	Q16633
<b>Unigene</b>	654525
<b>Immunogen</b>	Recombinant fragment (around aa 148-255) of human BOB1 (POU2AF1) protein (exact sequence is proprietary)
<b>Antibody target cellular localization</b>	Nucleus
<b>Verified antibody applications</b>	IHC (FFPE) (verified), WB (verified)
<b>Species reactivity</b>	Human
<b>Antibody application notes</b>	Staining of formalin-fixed tissues requires heating tissue sections in 10mM Tris with 1mM EDTA, pH 9.0, for 45 min at 95 °C followed by cooling at RT for 20 minutes. Immunohistology (formalin): 0.1-0.2 ug/mL for 30 minutes at RT, Western Blot 0.5-1 ug/mL. Optimal dilution for a specific application should be determined by user
<b>Positive control</b>	Ramos cells. Spleen or Tonsil
<b>Shipping condition</b>	Room temperature
<b>Storage Conditions</b>	Store at 2 to 8 °C. Protect fluorescent conjugates from light. Note: store BSA-free antibodies at -10 to -35 °C
<b>Regulatory status</b>	For research use only (RUO)
<b>Antibody/conjugate formulation</b>	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
<b>Validated in protein array</b>	Monospecific
<b>Shelf life</b>	Guaranteed for at least 24 months from date of receipt when stored as recommended
<b>Antibody research areas</b>	Cancer, Immunology
<b>Product origin</b>	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.
<b>Cell/tissue expression</b>	B-cells
<b>Tumor expression</b>	Leukemia/lymphoma

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>
BNC74	CF®740	742/767	633-685	775/50	<a href="#">CF®740 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, and Odyssey are registered trademarks of LI-COR Bioscience.

This datasheet was generated on June 4, 2026 at 01:33:31 AM. Visit product page to check for updated information before use. Product link: <https://biotium.com/product/bob-1-monoclonal-mouse-antibody-bob1-2422/>