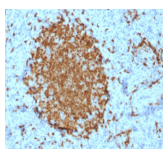


BOB-1 Monoclonal Mouse Antibody (BOB1/2421)



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 to inquire about stock status and lead times before placing your order.

Catalog number key for antibody number 2421, BOB.1 Monoclonal Mouse Antibody (BOB1/2421)

Product attributes

Antibody number	#2421
Antibody reactivity (target)	BOB.1
Antibody type	Primary
Host species	Mouse
Clonality	Monoclonal
Clone	BOB1/2421
Isotype	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
Immunogen	Recombinant fragment (around aa 148-255) of human BOB1 (POU2AF1) protein (exact sequence is proprietary)
Antibody target cellular localization	Nucleus
Verified antibody applications	Flow (intracellular) (verified), IHC (FFPE) (verified), WB (verified)
Species reactivity	Human
Positive control	Raji, 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
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
Shelf life	Guaranteed for at least 24 months from date of receipt when stored as recommended
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

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 July 15, 2025 at 01:56:07 PM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/bob-1-monoclonal-mouse-antibody-bob1-2421/>