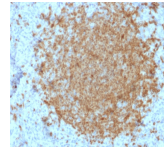


BOB-1 Monoclonal Mouse Antibody (BOB1/2422)



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 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

Antibody number	#2422
Antibody reactivity (target)	BOB.1
Antibody type	Primary
Host species	Mouse
Clonality	Monoclonal
Clone	BOB1/2422
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	IHC (FFPE) (verified), WB (verified)
Species reactivity	Human
Antibody application notes	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
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
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
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
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.