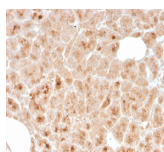


LMO2 Recombinant Monoclonal Mouse Antibody (rLMO2/1971)



Product Description

The LMO2 protein has a central and crucial role in hematopoietic development and is highly conserved. It has a particular function in normal and lymphatic endothelial cells involving the regulation of angiogenesis and lymph-angiogenesis. Immunohistochemical studies have also demonstrated expression of LMO2 in both normal germinal center B-cells and germinal center-derived B-cell lymphomas, including follicular lymphoma and diffuse large B-cell lymphoma. The use of anti-LMO2 is valuable as a tool in the identification of lymphomas of B-cell origin. LMO2 is useful in differentiating follicular lymphoma (LMO2+) from nodal marginal zone lymphoma (LMO2-). It also is positive in Hodgkin's and Burkitt's lymphomas.

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 3806, Anti-LMO2 (rLMO2/1971)

Call us : [800-304-5357](tel:800-304-5357)

Product attributes

Antibody number	#3806
Antibody reactivity (target)	LMO2
Antibody type	Primary
Host species	Mouse
Clonality	Monoclonal, Recombinant Monoclonal
Clone	rLMO2/1971
Isotype	IgG1, kappa
Molecular weight of antigen	24 kDa
Synonyms	Cysteine-rich protein TTG-2; LIM domain only protein 2; LMO-2; RBTN L1; RBTN2; Rhombotin like 1; Rhombotin-2 (RHOM2); T-cell translocation protein 2; TTG2
Human gene symbol	LMO2
Entrez gene ID	4005
SwissProt	P25791
Unigene	34560
Immunogen	Recombinant human LMO2 protein fragment (around aa 23-140) (exact sequence is proprietary)
Verified antibody applications	IHC (FFPE) (verified)
Antibody target cellular localization	Nucleus
Species reactivity	Human
Positive control	K562, Ramos or Raji cells. Placenta, pancreas or Hodgkin s lymphoma.
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
Shelf life	Guaranteed for at least 24 months from date of receipt when stored as recommended
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, Endothelial cells, Lymphatics
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 June 24, 2026 at 11:34:27 AM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/lmo2-recombinant-monoclonal-mouse-antibody-rlmo2-1971/>