LMO2 Recombinant Monoclonal Rabbit Antibody (LMO2/3147R)



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 3147, Anti-LMO2 (LMO2/3147R)

Product attributes

Tumor expression

Product attributes					
Antibody number	#3147				
Antibody reactivity (target)	102				
Antibody type	Primary				
Host species	Rabbit				
Clonality	Recombinant Monoclonal				
Clone	LMO2/3147R				
Isotype	IgG				
Molecular weight	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 fragment (around aa 23-140) human LMO2 protein (exact sequence is proprietary)				
Antibody target cellular localization	Nucleus				
Verified antibody applications	Flow (intracellular) (verified), IF (verified)				
Species reactivity					
Species reactivity	Human				
Antibody application notes	Human Flow cytometry: 1-2 ug/million cells; Immunofluorescence: 1-2 ug/mL; Optimal dilution for a specific application should be determined., Higher concentration may be required for direct detection using primary antibody conjugates than for indirect detection with secondary antibody				
	Flow cytometry: 1-2 ug/million cells; Immunofluorescence: 1-2 ug/mL; Optimal dilution for a specific application should be determined., Higher concentration may be required for direct detection using primary antibody conjugates than for indirect				
Antibody application notes	Flow cytometry: 1-2 ug/million cells; Immunofluorescence: 1-2 ug/mL; Optimal dilution for a specific application should be determined, Higher concentration may be required for direct detection using primary antibody conjugates than for indirect detection with secondary antibody K562, Ramos or Raji cells. Placenta, pancreas or Hodgkin s				
Antibody application notes Positive control	Flow cytometry: 1-2 ug/million cells; Immunofluorescence: 1-2 ug/mL; Optimal dilution for a specific application should be determined, Higher concentration may be required for direct detection using primary antibody conjugates than for indirect detection with secondary antibody K562, Ramos or Raji cells. Placenta, pancreas or Hodgkin s lymphoma.				
Antibody application notes Positive control Shipping condition	Flow cytometry: 1-2 ug/million cells; Immunofluorescence: 1-2 ug/ml.; Optimal dilution for a specific application should be determined., Higher concentration may be required for direct detection using primary antibody conjugates than for indirect detection with secondary antibody K562, Ramos or Raji cells. Placenta, pancreas or Hodgkin s lymphoma. Room temperature Store at 2 to 8 °C, Protect fluorescent conjugates from light,				
Antibody application notes Positive control Shipping condition Storage Conditions	Flow cytometry: 1-2 ug/million cells; Immunofluorescence: 1-2 ug/mL; Optimal dilution for a specific application should be determined., Higher concentration may be required for direct detection using primary antibody conjugates than for indirect detection with secondary antibody K562, Ramos or Raji cells. Placenta, pancreas or Hodgkin s lymphoma. Room temperature Store at 2 to 8 °C, Protect fluorescent conjugates from light, Note: store BSA-free antibodies at -10 to -35 °C				
Antibody application notes Positive control Shipping condition Storage Conditions Regulatory status Antibody/conjugate	Flow cytometry: 1-2 ug/million cells; Immunofluorescence: 1-2 ug/mL; Optimal dilution for a specific application should be determined., Higher concentration may be required for direct detection using primary antibody conjugates than for indirect detection with secondary antibody K562, Ramos or Raji cells. Placenta, pancreas or Hodgkin s lymphoma. Room temperature Store at 2 to 8 °C, Protect fluorescent conjugates from light, Note: store BSA-free antibodies at -10 to -35 °C For research use only (RUO) 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/0.05% BSA/0.05% azide, Purifier 1 mg/mL in PBS/0.05%				
Antibody application notes Positive control Shipping condition Storage Conditions Regulatory status Antibody/conjugate formulation	Flow cytometry: 1-2 ug/million cells; Immunofluorescence: 1-2 ug/mL; Optimal dilution for a specific application should be determined., Higher concentration may be required for direct detection using primary antibody conjugates than for indirect detection with secondary antibody K562, Ramos or Raji cells. Placenta, pancreas or Hodgkin s lymphoma. Room temperature Store at 2 to 8 °C, Protect fluorescent conjugates from light, Note: store BSA-free antibodies at -10 to -35 °C For research use only (RUO) 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 Guaranteed for at least 24 months from date of receipt when				

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), 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 November 4, 2025 at 03:19:59 PM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/lmo2-recombinant-monoclonal-rabbit-antibody-lmo2-3147r/