MAML3 Monoclonal Mouse Antibody (MAML3/1303)

Product Description

MAML3 (mastermind-like protein 3) is a nuclear speckle protein that acts as a transcriptional coactivator for Notch receptors. The Notch signaling pathway influences cell fate by regulating the ability of precursor cells to properly respond to developmental signals. MAML3 is a member of the mastermind-like family of proteins that are human homologs of the Drosophila melanogaster mastermind protein. Through its N-terminal region, MAML3 interacts with the ankyrin repeats of the Notch proteins Notch 1, Notch 2, Notch 3 and Notch 4.This interaction leads to formation of a DNA-binding complex with the Notch proteins and RBP-J, a complex that can then induce HES1 gene expression. While the N-terminal domain of MAML3 is essential for proper Notch binding, the C-terminal domain of MAML3 is essential for transcriptional activation. Due to its involvement in cell signaling and transcriptional activation, upregulation of MAML3 is thought to be involved in oncogenesis.

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 1303, Anti-MAML3 (MAML3/1303)

Product attributes

Call us: 800-304-5357

Product attributes			
Antibody number	#1303		
Antibody reactivity (target)	MAML3		
Antibody type	Primary		
Host species	Mouse		
Clonality	Monoclonal		
Clone	MAML3/1303		
Isotype	IgG1		
Molecular weight	150-170 kDa		
Synonyms	CAGH3; ERDA3; GDN; MAML3; Mastermind-like protein 3; TNRC3		
Human gene symbol	MAML3		
Entrez gene ID	55534 (Human)		
SwissProt	Q96JK9 (Human)		
Unigene	586165 (Human)		
Immunogen	Recombinant human full-length MAML3 protein		
Antibody target cellular localization	Nucleus		
Verified antibody applications	IHC (FFPE) (verified)		
Species reactivity	Human		
Antibody application notes	Staining of formalin-fixed tissues is enhanced by boiling tissue sections in 10 mM Tris, 1 mM EDTA pH 9.0 for 10-20 min followed by cooling at RT for 20 min, Immunofluorescence 0.5-1 ug/mL, Flow Cytometry 0.5-1 ug/million cells/0.1 mL, Optimal dilution for a specific application should be determined by user		
Positive control	MCF-7 cells. Pancreas or Placenta. Cervix or Colon Carcinoma.		
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		
Shelf life	Guaranteed for at least 24 months from date of receipt when stored as recommended		
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 i PBS without azide		
Antibody research areas	Developmental biology, Signal transduction		
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.		

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 09:03:43 AM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/mouse-monoclonal-anti-maml3-maml31303/