Albumin Monoclonal Mouse Antibody (ALB/2355)



Product Description

This MAb is Monospecific to albumin and shows no significant cross-reaction with other human proteins. Albumin is a soluble, monomeric protein, which comprises about one half of the blood serum protein. Albumin functions primarily as a carrier protein for steroids, fatty acids, and thyroid hormones and plays a role in stabilizing extracellular fluid volume. Albumin is synthesized in the liver as preproalbumin, which has an N-terminal peptide that is removed before the nascent protein is released from the rough endoplasmic reticulum. The product, proalbumin, is in turn cleaved in the Golgi vesicles to produce the secreted form of albumin.

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 2355, Albumin Monoclonal Mouse Antibody (ALB/2355)

Product attributes

research-areas He	355 matology pumin		
	6,7		
Antibody reactivity (target) Alb	numin		
			
Antibody type Pri	mary		
Host species Mo	ouse		
Clonality Mo	pnoclonal		
Clone AL	B/2355		
Isotype IgG	G2b, kappa		
Molecular weight 66	66 kDa		
Synonyms AL	ALB; Albumin		
Human gene symbol AL	ALB		
Entrez gene ID 213	213		
SwissProt P0	P02768		
Unigene 418	418167		
Immunogen Re	Recombinant full-length human ALB protein		
Verified antibody IF (applications	(verified), IHC (FFPE) (verified)		
Antibody target cellular Second localization	creted (extracellular)		
Species reactivity Hu	man		
prir sec For	gher concentration may be required for direct detection using mary antibody conjugates than for indirect detection with condary antibody, Immunofluorescence: 2-4 ug/mL; ELISA: or coating, order Ab without BSA; Optimal dilution for a specific plication should be determined.		
Positive control Hu	trol Human serum. Liver or Hepatocellular Carcinoma		
Shipping condition Ro	om temperature		
	ore at 2 to 8 °C, Protect fluorescent conjugates from light, te: store BSA-free antibodies at -10 to -35 °C		
Regulatory status For	r research use only (RUO)		
formulation cor in F	njugates: 0.1 mg/mL in PBS/0.1% BSA/0.05% azide, HRP njugates: 0.1 mg/mL in PBS/0.05% BSA, Purified: 0.2 mg/mL PBS/0.05% BSA/0.05% azide, Purified, BSA-free: 1 mg/mL in IS without azide		
Validated in protein array Mo	nospecific		
	aranteed for at least 24 months from date of receipt when ared as recommended		
boy	oduct may contain either bovine serum albumin (BSA) from vine serum (Bos taurus), or recombinant BSA produced in inese hamster ovary cells. Inquire for the specific lot.		

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 03:19:30 PM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/albumin-monoclonal-mouse-antibody-alb-2355/