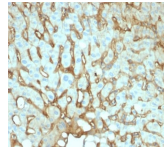


Albumin Monoclonal Mouse Antibody (ALB/2144)



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 prealbumin, 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 2144, Anti-Albumin (ALB/2144)**

Product attributes

research-areas	Hematology
Antibody number	#2144
Antibody reactivity (species)	Albumin
Antibody type	Primary
Host species	Mouse
Clonality	Monoclonal
Clone	ALB/2144
Isotype	IgG1, kappa
Molecular weight	66 kDa
Synonyms	ALB; Albumin
Human gene symbol	ALB
Entrez gene ID	213
SwissProt	P02768
Unigene	418167
Immunogen	Recombinant full-length human ALB protein
Antibody target cellular localization	Secreted (extracellular)
Verified antibody application	IHC (FFPE) (verified)
Species reactivity	Human
Antibody application notes	Immunohistology (formalin): 1-2 ug/mL for 30 minutes at RT. Staining of formalin-fixed tissues requires boiling tissue sections in 10 mM citrate buffer, pH 6.0, for 10-20 minutes followed by cooling at RT for 20 minutes, ELISA For coating, order Ab without BSA, Optimal dilution for a specific application should be determined by user
Positive control	Human serum. Liver or Hepatocellular 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 buffer 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
Antibody research areas	Hematology

Catalog Number Key for Primary Antibody Conjugates

Antibody # prefix	Conjugation	Ex/Em (nm)	Laser line	Detection channel	Dye Features
BNC04	CF@405S	404/431	405	DAPI (microscopy), AF405	CF@405S Features
BNC05	CF@405M	408/452	405	DAPI (microscopy), Pacific Blue™	CF@405M Features
BNC06	CF@405L	395/545	405	Pacific Orange™	CF@405L Features
BNC88	CF@488A	490/515	488	GFP, FITC	CF@488A Features
BNC14	CF@514	516/548	488	GFP, FITC	CF@514 Features
BNC43	CF@543	541/560	532	RFP, TRITC	CF@543 Features
BNC55	CF@555	555/565	532, 561	RFP, TRITC	CF@555 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
BNC60	CF@660C	667/685	633-640	Cy@5	CF@660C Features
BNC61	CF@660R	663/682	633-640	Cy@5	CF@660R Features
BNC80	CF@680	681/698	633-640	Cy@5.5, IR@700	CF@680 Features
BNC81	CF@680R	680/701	633-640	Cy@5.5, IR@700	CF@680R Features
BNC00	CF@700	695/720	633-640	AF700	CF@700 Features
BNC70	CF@770	770/797	785, 808	IR@800	CF@770 Features
BNCR	R-PE (PE)	496, 546, 565/578	488-561	PE	
BNCA	APC	650/660	594-633	APC	
BNCP	PerCP	482/677	488-561	PerCP, PE-Cy@5	
BNCB	Biotin	N/A	N/A	N/A	
BNCAP	Alkaline Phosphatase	N/A	N/A	N/A	
BNCH	HRP	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 GE Healthcare; IRDye, LI-COR, and Odyssey are registered trademarks of LI-COR Bioscience.