Bovine Serum Albumin Monoclonal Mouse Antibody (ALB/398)

Product Description

This antibody reacts with a protein of ~66 kDa, identified as bovine serum albumin (BSA). It is a high affinity antibody and can be used for detection of traces of BSA. BSA is an abundant plasma protein in cows that is important for maintaining osmotic pressure in blood plasma for proper distribution of body fluids between intravascular compartments and body tissues. BSA is a common buffer component in immunoglobulin type assays due to good solubility characteristics for water, calcium, sodium, potassium, fatty acids, hormones and bilirubin. BSA makes up about half of the protein in plasma and represents the most stable and soluble protein in the plasma. It is a suitable reagent for laboratories developing immunoassays, mostly due to its availability, solubility and the numerous functional groups present for coupling. BSA contains several lysines that are capable of reacting with conjugation sites of linkers, making it applicable as a carrier protein for antigenic compounds.

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 0398, Anti-Bovine Serum Albumin (ALB/398)

Product attributes

Product attributes	
Antibody number	#0398
Antibody reactivity (target)	Bovine Serum Albumin
Antibody type	Primary
Host species	Mouse
Clonality	Monoclonal
Clone	ALB/398
Isotype	lgG1
Molecular weight	66 kDa
Synonyms	ALB; BSA; Growth inhibiting protein 20
Entrez gene ID	280717 (Cow)
SwissProt	P02769 (Cow)
Unigene	Bt.106669 (Cow)
Immunogen	Purified bovine serum albumin
Antibody target cellular localization	Secreted (extracellular), Vesicular
Species reactivity	Cow
Species reactivity Antibody application notes	For coating for ELISA, order Ab without BSA, Higher concentration may be required for direct detection using primary antibody conjugates than for indirect detection with secondary antibody, Optimal dilution and staining procedure for a specific application should be determined by user, Recommended starting concentrations for titration are 1-2 ug/mL for most applications, or 1 ug/million cells/100 uL for flow cytometry, Does not react with human, rabbit, dog, or guinea pig, others not tested
	For coating for ELISA, order Ab without BSA, Higher concentration may be required for direct detection using primary antibody conjugates than for indirect detection with secondary antibody, Optimal dilution and staining procedure for a specific application should be determined by user, Recommended starting concentrations for titration are 1-2 ug/mL for most applications, or 1 ug/million cells/100 uL for flow cytometry, Does not react with human, rabbit, dog, or guinea pig, others not
Antibody application notes	For coating for ELISA, order Ab without BSA, Higher concentration may be required for direct detection using primary antibody conjugates than for indirect detection with secondary antibody, Optimal dilution and staining procedure for a specific application should be determined by user, Recommended starting concentrations for titration are 1-2 ug/mL for most applications, or 1 ug/million cells/100 uL for flow cytometry, Does not react with human, rabbit, dog, or guinea pig, others not tested
Antibody application notes Positive control	For coating for ELISA, order Ab without BSA, Higher concentration may be required for direct detection using primary antibody conjugates than for indirect detection with secondary antibody. Optimal dilution and staining procedure for a specific application should be determined by user, Recommended starting concentrations for titration are 1-2 ug/mL for most applications, or 1 ug/million cells/100 L for flow cytometry, Does not react with human, rabbit, dog, or guinea pig, others not tested Bovine serum albumin. Cow liver
Antibody application notes Positive control Shipping condition	For coating for ELISA, order Ab without BSA, Higher concentration may be required for direct detection using primary antibody conjugates than for indirect detection with secondary antibody, Optimal dilution and staining procedure for a specific application should be determined by user, Recommended starting concentrations for titration are 1-2 ug/mL for most applications, or 1 ug/million cells/100 uL for flow cytometry, Does not react with human, rabbit, dog, or guinea pig, others not tested Bovine serum albumin. Cow liver Room temperature Store at 2 to 8 °C, Protect fluorescent conjugates from light,
Antibody application notes Positive control Shipping condition Storage Conditions	For coating for ELISA, order Ab without BSA, Higher concentration may be required for direct detection using primary antibody conjugates than for indirect detection with secondary antibody, Optimal dilution and staining procedure for a specific application should be determined by user, Recommended starting concentrations for titration are 1-2 ug/ml. for most applications, or 1 ug/million cells/100 ul. for flow cytometry, Does not react with human, rabbit, dog, or guinea pig, others not tested Bovine serum albumin. Cow liver Room temperature Store at 2 to 8 °C, Protect fluorescent conjugates from light, Note: store BSA-free antibodies at -10 to -35 °C Guaranteed for at least 24 months from date of receipt when

Call us: 800-304-5357 Email: btinfo@biotium.com

Antibody # prefix	Conjugation	Ex/Em (nm)	Laser line	Detection channel	Concentration	Supplied in	Dye Features
BNC04	CF®405S	404/431	405	DAPI (microscopy), AF405	Place custom order	PBS/0.05% azide/no BSA	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
BNCB	Biotin	N/A	N/A	N/A			
BNUM	Purified, BSA-free	N/A	N/A	N/A	1 mg/mL	PBS, no BSA, no azide	

The prefix indicates conjugation, followed by the antibody number and size suffix.

Alexa Fluor, Pacific Blue 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 6, 2025 at 05:47:51 AM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/monoclonal-anti-bovine-serum-albumin-bsa-alb398/