S100A2 / S100 Calcium Binding Protein A2 Monoclonal Mouse Antibody (CPTC-S100A2-2)



Product Description

S100A2 belongs to the S100 family of proteins containing 2 EF-hand calcium-binding motifs. S100A2 may function as a modulator against excess calcium accumulation in normal human mammary epithelial cells and also have a role in suppressing tumor cell growth. This protein may have a tumor suppressor function. Chromosomal rearrangements and altered expression of this gene have been implicated in breast cancer. Cytoplasmic overexpression may also be of prognostic significance when observed in oral cancer patients, and the S100A2 has also been identified as significantly down-regulated in gastric cancer.

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 2591, Anti-S100 Calcium Binding Protein A2|S100A2 (CPTC-S100A2-2)

Product attributes

Call us: 800-304-5357

Product attributes				
Antibody number	#2591			
Antibody reactivity (target)	S100 Calcium Binding Protein A2, S100A2			
Antibody type	Primary			
Host species	Mouse			
Clonality	Monoclonal			
Clone	CPTC-S100A2-2			
Isotype	IgG2b, kappa			
Molecular weight	11.12 kDa			
Synonyms	CAN19; MGC111539; S100L			
Human gene symbol	S100A2			
Entrez gene ID	6273			
SwissProt	P29034			
Unigene	516484			
Immunogen	Recombinant full-length human S100A2 protein			
Verified antibody applications	Flow (intracellular) (verified), IF (verified), IHC (FFPE) (verified), WB (verified)			
Antibody target cellular localization	Nucleus & cytoplasm			
Species reactivity	Human			
Positive control	HeLa. A549, HACAT, BxPC3 and A431 cell lysates; Skin and squamous cell cervical 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			
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 p PBS/0.05% BSA/0.05% azide, Purified, BSA-free: 1 mg/mL in PBS without azide			
Validated in protein array	Monospecific			
Shelf life	Guaranteed for at least 24 months from date of receipt when stored as recommended			
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.			

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 30, 2025 at 04:25:18 AM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/s100a2-s100-calcium-binding-protein-a2-monoclonal-mouse-antibody-cptc-s100a2-2/