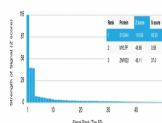


S100A4 Monoclonal Mouse Antibody (CPTC-S100A4-3)



Product Description

S100A4 belongs to the S100 super-family of proteins containing 2 EF-hand calcium-binding domains. S100A4 has been implicated in the progression and prognosis of several forms of human cancer, e. g. breast, colorectal, gastric, pancreatic and bladder cancer, SCLC and oesophageal squamous cell carcinoma, among others. Poor prognosis associated with high S100A4 expression is accompanied by clear signs of disease progression, e. g. high histological and clinical grades and involvement of lymph nodes. Also indicative of poor prognosis is high S100A4 expression coupled with reduced E-cadherin expression in pancreatic, oral squamous cell carcinoma and in melanoma. S100A4 expression is inversely related with expression of metastasis suppressor nm23 and with prognosis of breast cancer. S100A4 is overexpressed in highly metastatic cancers, which makes it useful as a marker of tumor progression.

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 2224, Anti-Calvasculin|Metastasin|S100A4 (CPTC-S100A4-3)

Product attributes

Antibody number	#2224
Antibody reactivity (target)	Calvasculin, Metastasin, S100A4
Antibody type	Primary
Host species	Mouse
Clonality	Monoclonal
Clone	CPTC-S100A4-3
Isotype	IgG2c, kappa
Molecular weight	12 kDa
Synonyms	S100 calcium-binding protein A4; Calvasculin; CAPL; Fibroblast specific protein 1 (FSP1); Leukemia multidrug resistance associated protein; Malignant transformation suppression 1 (MTS1); Metastasin; Placental calcium-binding protein
Human gene symbol	S100A4
Entrez gene ID	6275
SwissProt	P26447
Unigene	654444
Immunogen	Recombinant human S100A4 full length protein
Antibody target cellular localization	Nucleus & cytoplasm
Verified antibody applications	Flow (intracellular) (verified), IF (verified)
Species reactivity	Human
Expected antibody applications	IHC (FFPE) (published for clone)
Antibody application notes	Flow cytometry: 1-2 ug/million cells; Immunofluorescence: 1-2 ug/mL; Western blot 1-2 ug/mL; Optimal dilution for a specific application should be determined., Higher concentration may be required for direct detection using primary antibody conjugates than for indirect detection with secondary antibody
Positive control	T98G, HeLa, A549 or A375 cells. Placenta, small intestine.
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 in 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.
Tumor expression	Breast cancer

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.

References

Note: References for this clone sold by other suppliers may be listed for expected applications.

Stroke (2013) 44: 1456-1458. (IHC, FFPE)

This datasheet was generated on July 12, 2025 at 02:38:16 AM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/s100a4-monoclonal-mouse-antibody-cptc-s100a4-3/>