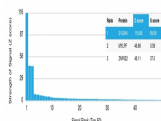


# S100A4 Monoclonal Mouse Antibody (CPTC-S100A4-3)



## 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](mailto: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

<b>Antibody number</b>	#2224
<b>Antibody reactivity (target)</b>	Calvasculin, Metastasin, S100A4
<b>Antibody type</b>	Primary
<b>Host species</b>	Mouse
<b>Clonality</b>	Monoclonal
<b>Clone</b>	CPTC-S100A4-3
<b>Isotype</b>	IgG2c, kappa
<b>Molecular weight</b>	12 kDa
<b>Synonyms</b>	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
<b>Human gene symbol</b>	S100A4
<b>Entrez gene ID</b>	6275
<b>SwissProt</b>	P26447
<b>Unigene</b>	654444
<b>Immunogen</b>	Recombinant human S100A4 full length protein
<b>Antibody target cellular localization</b>	Nucleus & cytoplasm
<b>Verified antibody applications</b>	Flow (intracellular) (verified), IF (verified)
<b>Species reactivity</b>	Human
<b>Expected antibody applications</b>	IHC (FFPE) (published for clone)
<b>Antibody application notes</b>	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
<b>Positive control</b>	T98G, HeLa, A549 or A375 cells. Placenta, small intestine.
<b>Shipping condition</b>	Room temperature
<b>Storage Conditions</b>	Store at 2 to 8 °C, Protect fluorescent conjugates from light, Note: store BSA-free antibodies at -10 to -35 °C
<b>Regulatory status</b>	For research use only (RUO)
<b>Antibody/conjugate formulation</b>	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
<b>Validated in protein array</b>	Monospecific
<b>Shelf life</b>	Guaranteed for at least 24 months from date of receipt when stored as recommended
<b>Product origin</b>	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.
<b>Tumor expression</b>	Breast cancer

Antibody # prefix	Conjugation	Ex/Em (nm)	Laser line	Detection channel	Dye Features
BNC04	CF®405S	404/431	405	DAPI (microscopy), AF405	<a href="#">CF®405S Features</a>
BNC88	CF®488A	490/515	488	GFP, FITC	<a href="#">CF®488A Features</a>
BNC68	CF®568	562/583	532, 561	RFP, TRITC	<a href="#">CF®568 Features</a>
BNC94	CF®594	593/614	561	Texas Red®	<a href="#">CF®594 Features</a>
BNC40	CF®640R	642/662	633-640	Cy®5	<a href="#">CF®640R Features</a>
BNC47	CF®647	650/665	633-640	Cy®5	<a href="#">CF®647 Features</a>
BNC74	CF®740	742/767	633-685	775/50	<a href="#">CF®740 Features</a>
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 November 27, 2024 at 08:32:55 AM. Visit product page to check for updated information before use.

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