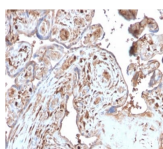


# S100A4 Recombinant Monoclonal Rabbit Antibody (S100A4/2750R)



## Product Description

S100A4 belongs to the S100 super-family of proteins containing 2 EF-hand calcium-binding domains. S100 genes include at least 25 members, including S100A1-S100A18, trichohyalin, filaggrin, repetin, S100P, and S100Z. S100A4 exerts its function via direct interaction with a number of proteins including P53, P63, non-muscle myosin IIA, alpha6beta4 integrin, and liprin b1. 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 2750, Anti-S100A4 (S100A4/2750R)**

Call us : [800-304-5357](tel:800-304-5357)

## Product attributes

Antibody number	#2750
Antibody reactivity (target)	S100A4
Antibody type	Primary
Host species	Rabbit
Clonality	Recombinant Monoclonal
Clone	S100A4/2750R
Isotype	IgG
Molecular weight	10-12 kDa
Synonyms	S100A4; 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 fragment of human S100A4 protein (around aa1-200) (exact sequence is proprietary)
Antibody target cellular localization	Nucleus & cytoplasm
Species reactivity	Human, Mouse
Verified antibody applications	Flow (intracellular) (verified), IF (verified), IHC (FFPE) (verified), WB (verified)
Antibody application notes	Higher concentration may be required for direct detection using primary antibody conjugates than for indirect detection with secondary antibody. Immunohistochemistry (formalin): 0.5-1.0 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. Optimal dilution for a specific application should be determined by user
Positive control	HeLa, A549 or A375 Cells. Placenta
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 research areas	Cancer
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
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	<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.

This datasheet was generated on February 1, 2026 at 02:23:42 AM. Visit product page to check for updated information before use.  
Product link: <https://biotium.com/product/s100a4-recombinant-monoclonal-rabbit-antibody-s100a4-2750r/>