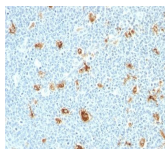


Calprotectin (MRP14)(S100A9) Monoclonal Mouse Antibody (CAGB/426)



Product Description

Calprotectin is a heteroduplex of Protein S100-A8/A9 (also called MRP-8/14, or Calgranulin A/B). This antibody recognizes the 14 kDa B component (Protein S100-A9) of the complex. Protein S100-A9 is a member of the S100 family of proteins containing 2 EF-hand calcium-binding motifs. Altered expression of this protein is associated with the disease cystic fibrosis. This antibody reacts with neutrophils, monocytes, macrophages, and squamous mucosal epithelia and has been shown as an important marker for identifying macrophages in tissue sections.

This antibody is available purified with BSA/azide at 200 ug/mL, or BSA/azide-free at 1 mg/mL. **Catalog number key for antibody number 0426, Anti-Calprotectin|S100-A9 (CAGB/426)**

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

Product attributes

Antibody number	#0426
Antibody reactivity (target)	Calprotectin, S100-A9
Antibody type	Primary
Host species	Mouse
Clonality	Monoclonal
Clone	CAGB/426
Isotype	IgM, kappa
Molecular weight	14 kDa
Synonyms	Calprotectin; Protein S100-A9; Myeloid-related protein 14; MRP-14; Calgranulin B; CAGB; Calprotectin L1H Subunit
Entrez gene ID	6280 (S100A9 / Calgranulin B / MRP-14)
SwissProt	P06702 (S100A9 / Calgranulin B / MRP-14)
Human gene symbol	S100A9
Unigene	112405 (S100A9 / Calgranulin B / MRP-14)
Immunogen	Recombinant human CAGB protein
Antibody target cellular localization	Cytoplasmic, Cytoskeleton
Cell/tissue expression	Granulocytes, Monocytes/macrophages
Species reactivity	Human
Verified antibody applications	IHC (FFPE) (verified)
Antibody application notes	Immunohistology formalin-fixed. Not suitable for frozen tissues. 0.5-1 ug/mL. Staining of formalin-fixed tissues requires boiling tissue sections in 10 mM citrate buffer, pH 6.0, for 10-20 min followed by cooling at RT for 20 minutes. Immunofluorescence 0.5-1 ug/mL. Flow Cytometry 0.5-1 ug/million cells/0.1 mL. Optimal dilution for a specific application should be determined by user
Positive control	Tonsil, lymph node, or spleen
Shipping condition	Room temperature
Storage Conditions	Store at 2 to 8 °C, 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	Immunology, Inflammation
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	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 December 7, 2025 at 06:41:16 PM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/monoclonal-anti-calgranulin-b-cagb426/>