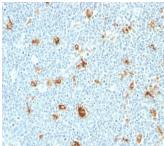


Calprotectin / Protein S100-A9 Monoclonal Mouse Antibody (CAGB/426)



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.

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