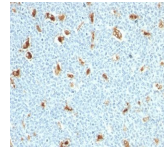


S100A9 Monoclonal Mouse Antibody (S100A9/1011)

Recognizes a 14 kDa protein, identified as S100A9 (also known as Calgranulin B or MRP-14); expressed by granulocytes, monocytes and by tissue macrophages.



Product Description

Recognizes a 14 kDa protein, identified as S100A9 (also known as Calgranulin B or MRP-14); expressed by granulocytes, monocytes and by tissue macrophages. The protein encoded by this gene is a member of the S100 family of proteins containing 2 EF-hand calcium-binding motifs. S100 proteins are localized in the cytoplasm and/or nucleus of a wide range of cells, and involved in the regulation of a number of cellular processes such as cell cycle progression and differentiation. Altered expression of this protein is associated with the disease cystic fibrosis. This MAb reacts with neutrophils, monocytes and macrophages, and has been shown as an important marker for identifying macrophages in tissue sections. Among cells that are now recognized as macrophages are histiocytes, Kupffer cells, osteoclasts, microglial cells, synovial type A cells, interdigitating cells, and Langerhans cells (in normal tissues) and epithelioid cells and Langerhans-type and foreign-body-type multinucleated giant cells (in inflamed tissues). This antibody is available purified with BSA/azide at 200 ug/mL, or BSA/azide-free at 1 mg/mL.

References

Cross, S.S., et al. 2005. Expression of S100 proteins in normal human tissues and common cancers using tissue microarrays: S100A6, S100A8, S100A9 and S100A11 are all overexpressed in common cancers. *Histopathology* 46: 256-269

Product attributes	
Antibody number	1011
Reactivity (target)	S100A9
Antibody type	Primary
Host species	Mouse
Clonality	Monoclonal
Clone	S100A9/1011
Isotype	IgM, kappa
Molecular weight	14 kDa
Synonyms	S100-A9; MRP14; calgranulin-B; calprotectin L1H subunit; leukocyte L1 complex heavy chain; migration inhibitory factor-related protein 14; S100 calcium-binding protein A9 (calgranulin B)
Human gene symbol	S100A9
Entrez gene ID	6280
SwissProt	P06702
Unigene	112405
Immunogen	Recombinant human S100A9 protein
Cellular localization	Cytoplasmic, Membrane/cell surface, Secreted (extracellular)
Species reactivity	Human
Applications	Immunofluorescence, Immunohistology (formalin), Flow cytometry
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	Liver or histiocytoma
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)
Supplied As	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, BSA-free: 1 mg/mL in PBS without azide, Purified: 0.2 mg/mL in PBS/0.05% BSA/0.05% azide
Antibody research areas	Inflammation
Cell/tissue expression	Monocytes/macrophages