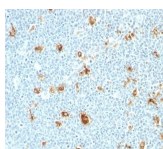


# Calprotectin (MRP14)(S100A9) Monoclonal Mouse Antibody (CAGB/426)



## 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)**

## Product attributes

|  |   |
|--|---|
| <b>Antibody number</b>                       | #0426   |
| <b>Antibody reactivity (target)</b>          | Calprotectin, S100-A9   |
| <b>Antibody type</b>                         | Primary   |
| <b>Host species</b>                          | Mouse   |
| <b>Clonality</b>                             | Monoclonal  |
| <b>Clone</b>                                 | CAGB/426  |
| <b>Isotype</b>                               | IgM, kappa  |
| <b>Molecular weight</b>                      | 14 kDa  |
| <b>Synonyms</b>                              | Calprotectin; Protein S100-A9; Myeloid-related protein 14; MRP-14; Calgranulin B; CAGB; Calprotectin L1H Subunit  |
| <b>Entrez gene ID</b>                        | 6280 (S100A9 / Calgranulin B / MRP-14)  |
| <b>SwissProt</b>                             | P06702 (S100A9 / Calgranulin B / MRP-14)  |
| <b>Human gene symbol</b>                     | S100A9  |
| <b>Unigene</b>                               | 112405 (S100A9 / Calgranulin B / MRP-14)  |
| <b>Immunogen</b>                             | Recombinant human CAGB protein  |
| <b>Antibody target cellular localization</b> | Cytoplasmic, Cytoskeleton   |
| <b>Cell/tissue expression</b>                | Granulocytes, Monocytes/macrophages   |
| <b>Species reactivity</b>                    | Human   |
| <b>Verified antibody applications</b>        | IHC (FFPE) (verified)   |
| <b>Antibody application notes</b>            | 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 |
| <b>Positive control</b>                      | Tonsil, lymph node, or spleen   |
| <b>Shipping condition</b>                    | Room temperature  |
| <b>Storage Conditions</b>                    | Store at 2 to 8 °C, Note: store BSA-free antibodies at -10 to -35 °C  |
| <b>Shelf life</b>                            | Guaranteed for at least 24 months from date of receipt when stored as recommended   |
| <b>Regulatory status</b>                     | For research use only (RUO)   |
| <b>Antibody research areas</b>               | Immunology, Inflammation  |
| <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>Product origin</b>                        | Product may contain either bovine serum albumin (BSA) from bovine serum ( <i>Bos taurus</i> ), 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 November 26, 2024 at 05:45:13 PM. Visit product page to check for updated information before use.  
Product link: <https://biotium.com/product/monoclonal-anti-calgranulin-b-cagb426/>