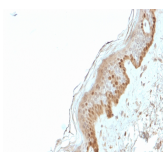


S100A2 / S100 Calcium Binding Protein A2 Monoclonal Mouse Antibody (CPTC-S100A2-2)



Product Description

S100A2 belongs to the S100 family of proteins containing 2 EF-hand calcium-binding motifs. S100A2 may function as a modulator against excess calcium accumulation in normal human mammary epithelial cells and also have a role in suppressing tumor cell growth. This protein may have a tumor suppressor function. Chromosomal rearrangements and altered expression of this gene have been implicated in breast cancer. Cytoplasmic overexpression may also be of prognostic significance when observed in oral cancer patients, and the S100A2 has also been identified as significantly down-regulated in gastric cancer.

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 to inquire about stock status and lead times before placing your order.

Catalog number key for antibody number 2591, Anti-S100 Calcium Binding Protein A2|S100A2 (CPTC-S100A2-2)

Product attributes

| | |
|---------------------------------------|---|
| Antibody number | #2591 |
| Antibody reactivity (target) | S100 Calcium Binding Protein A2, S100A2 |
| Antibody type | Primary |
| Host species | Mouse |
| Clonality | Monoclonal |
| Clone | CPTC-S100A2-2 |
| Isotype | IgG2b, kappa |
| Molecular weight | 11.12 kDa |
| Synonyms | CAN19; MGC111539; S100L |
| Human gene symbol | S100A2 |
| Entrez gene ID | 6273 |
| SwissProt | P29034 |
| Unigene | 516484 |
| Immunogen | Recombinant full-length human S100A2 protein |
| Verified antibody applications | Flow (intracellular) (verified), IF (verified), IHC (FFPE) (verified), WB (verified) |
| Antibody target cellular localization | Nucleus & cytoplasm |
| Species reactivity | Human |
| Positive control | HeLa, A549, HACAT, BxPC3 and A431 cell lysates; Skin and squamous cell cervical carcinoma. |
| 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 |
| Regulatory status | For research use only (RUO) |
| 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 |
| Validated in protein array | Monospecific |
| Shelf life | Guaranteed for at least 24 months from date of receipt when stored as recommended |
| 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 April 23, 2025 at 11:45:46 PM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/s100a2-s100-calcium-binding-protein-a2-monoclonal-mouse-antibody-cptc-s100a2-2/>