

Angiotensin I Converting Enzyme / CD143 Monoclonal Mouse Antibody (9B9)

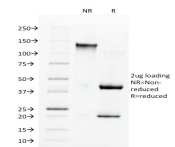
Product Description

Angiotensin-converting enzyme (ACE), also known as CD143, encodes an enzyme involved in catalyzing the conversion of angiotensin I into a physiologically active peptide angiotensin II. Angiotensin II is a potent vasopressor and aldosterone-stimulating peptide that controls blood pressure and fluid-electrolyte balance. This enzyme plays a key role in the renin-angiotensin system. Many studies have associated the presence or absence of a 287 bp Alu repeat element in this gene with the levels of circulating enzyme or cardiovascular pathophysiologies. Two most abundant alternatively spliced variants of this gene encode two isozymes – the somatic form and the testicular form that are equally active. Multiple additional alternatively spliced variants have been identified but their full length nature has not been determined.

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 1560, Angiotensin I Converting Enzyme / CD143 Monoclonal Mouse Antibody (9B9)



Call us : [800-304-5357](tel:800-304-5357) Email: techsupport@biotium.com

Product attributes

| | |
|---------------------------------------|---|
| Antibody number | #1560 |
| Antibody reactivity (target) | Angiotensin I Converting Enzyme, CD143 |
| Antibody type | Primary |
| Host species | Mouse |
| Clonality | Monoclonal |
| Clone | 9B9 |
| Isotype | IgG1, kappa |
| Molecular weight | 195 kDa |
| Synonyms | angiotensin 1 converting enzyme; angiotensin-converting enzyme; ACE; carboxycathepsin; CD143; dipeptidyl carboxypeptidase 1; Kininase II; peptidase P |
| Human gene symbol | ACE |
| Entrez gene ID | 1636 |
| SwissProt | P12821 |
| Unigene | 298469 |
| Immunogen | Full-length recombinant human ACE/CD143 protein |
| Verified antibody applications | IHC (FFPE) (verified) |
| Antibody target cellular localization | Plasma membrane |
| Species reactivity | Human |
| Positive control | Ubiquitously expressed, with highest levels in lung, kidney, heart, gastrointestinal system and prostate. |
| 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 |
| 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 August 26, 2025 at 05:26:49 AM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/angiotensin-i-converting-enzyme-cd143-monoclonal-mouse-antibody-9b9/>