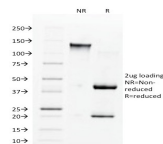


Angiotensin I Converting Enzyme / CD143 Monoclonal Mouse Antibody (9B9)



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 pathophysiology. 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)**

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; carboxypeptidase; 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

Antibody # prefix	Conjugation	Ex/Em (nm)	Laser line	Detection channel
BNC04	CF®405S	404/431	405	DAPI (microscopy), AF405
BNC88	CF®488A	490/515	488	GFP, FITC
BNC68	CF®568	562/583	532, 561	RFP, TRITC
BNC94	CF®594	593/614	561	Texas Red®
BNC40	CF®640R	642/662	633-640	Cy®5
BNC47	CF®647	650/665	633-640	Cy®5
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

Dye Features

[CF®405S Features](#)

[CF®488A Features](#)

[CF®568 Features](#)

[CF®594 Features](#)

[CF®640R Features](#)

[CF®647 Features](#)

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, LI-COR Bioscience.

This datasheet was generated on May 14, 2024 at 02:42:41 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/>