

PRODUCT INFORMATION



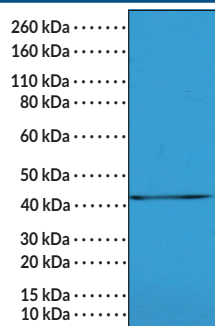
α -SMA (N-Term) Rabbit Monoclonal Antibody (Clone RM253)

Item No. 32204

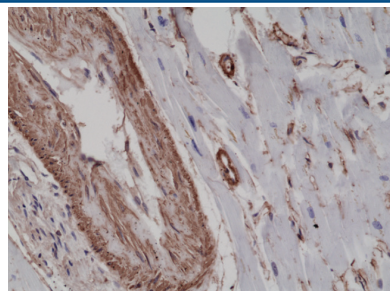
Overview and Properties

Contents:	This vial contains 100 μ l of protein A-affinity purified monoclonal antibody.
Synonyms:	α -Actin 2, Smooth Muscle, ACTA2, α -Smooth Muscle Actin
Immunogen:	Synthetic peptide corresponding to the N-terminus of human α -SMA
Species Reactivity:	(+) Human, mouse
Cross Reactivity:	(+) α -SMA
Form:	Liquid
Storage:	-20°C (as supplied)
Stability:	\geq 1 year
Storage Buffer:	PBS with 50% glycerol, 1% BSA, and 0.09% sodium azide
Concentration:	50 μ g/ml
Clone:	RM253
Host:	Rabbit
Isotype:	IgG
Applications:	Immunohistochemistry (IHC) and Western blot (WB); the recommended starting dilution for IHC is 1:1,100-1:2,500 and 1:1,100-1:2,000 for WB. Other applications were not tested, therefore optimal working concentration/dilution should be determined empirically.

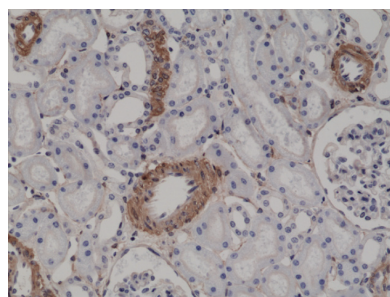
Images



WB of mouse heart tissue lysate using α -SMA (N-term) Rabbit Monoclonal Antibody (Clone RM253) at a dilution of 1:1,000.



Immunohistochemical staining of formalin-fixed and paraffin-embedded human heart tissue using α -SMA (N-term) Rabbit Monoclonal Antibody (Clone RM253) at a dilution of 1:2,500.



Immunohistochemical staining of formalin-fixed and paraffin-embedded human kidney tissue using α -SMA (N-term) Rabbit Monoclonal Antibody (Clone RM253) at a dilution of 1:2,500.

WARNING
THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA
This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the [complete](#) Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY
Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 01/25/2024

CAYMAN CHEMICAL
1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA
PHONE: [800] 364-9897
[734] 971-3335
FAX: [734] 971-3640
CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM

PRODUCT INFORMATION



Description

α -Smooth muscle actin (α -SMA) is a cytoskeletal protein encoded by *ACTA2* in humans and a component of the smooth muscle cell (SMC) contractile unit.¹ It is expressed in the cytosol of SMCs throughout the vascular system where it polymerizes to form the thin filaments of the SMC contractile unit. α -SMA is also expressed in pericytes and myofibroblasts.² *Acta2* expression is upregulated in, and positively correlates with enhanced migration and contraction of, hepatic stellate cells, which are precursors to myofibroblasts, in a rat model of carbon tetrachloride-induced liver injury. Missense mutations in *ACTA2* cause approximately 14% of inherited ascending thoracic aortic aneurysms and dissections (TAAD) via induction of defective SMC contraction, aortic medial degeneration, and SMC hyperplasia or disarray.³ Mutations in *ACTA2* are also positively correlated with a predisposition for various additional vascular diseases including premature coronary artery disease, stroke, and Moyamoya disease.⁴ Cayman's α -SMA (N-Term) Rabbit Monoclonal Antibody (Clone RM253) can be used for immunohistochemistry (IHC) and Western blot (WB) applications.

References

1. Karimi, A. and Milewicz, D.M. Structure of the elastin-contractile units in the thoracic aorta and how genes that cause thoracic aortic aneurysms and dissections disrupt this structure. *Can. J. Cardiol.* **32(1)**, 26-34 (2016).
2. Rockey, D.C., Weymouth, N., and Shi, Z. Smooth muscle α actin (*Acta2*) and myofibroblast function during hepatic wound healing. *PLoS One* **8(10)**, e77166 (2013).
3. Cho, D.-C., Pannu, H., Tran-Fadulu, V., *et al.* Mutations in smooth muscle alpha-actin (*ACTA2*) lead to thoracic aortic aneurysms and dissections. *Nat. Genet.* **39(12)**, 1448-1493 (2007).
4. Guo, D.-C., Papke, C.L., Tran-Fadulu, V., *et al.* Mutations in smooth muscle alpha-actin (*ACTA2*) cause coronary artery disease, stroke, and Moyamoya disease, along with thoracic aortic disease. *Am. J. Hum. Genet.* **84(5)**, 617-627 (2009).

CAYMAN CHEMICAL
1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA
PHONE: [800] 364-9897
[734] 971-3335
FAX: [734] 971-3640
CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM