

PRODUCT INFORMATION



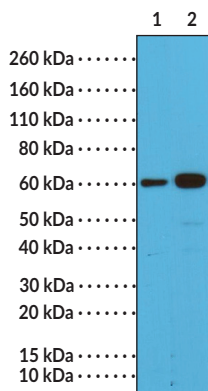
SMAD4 (C-Term) Rabbit Monoclonal Antibody (Clone RM277)

Item No. 32227

Overview and Properties

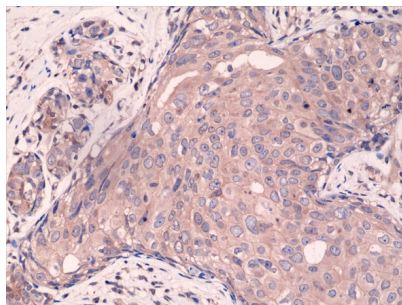
Contents:	This vial contains 100 µl of protein A-affinity purified monoclonal antibody.
Synonyms:	Deleted in Pancreatic Carcinoma Locus 4, DPC4, MADH4, Mothers Against Decapentalplegic Homolog 4
Immunogen:	Peptide from the C-terminal region of human SMAD4
Cross Reactivity:	(+) SMAD4
Species Reactivity:	(+) Human
Form:	Liquid
Storage:	-20°C (as supplied)
Stability:	≥1 year
Storage Buffer:	PBS with 50% glycerol, 1% BSA, and 0.09% sodium azide
Clone:	RM277
Host:	Rabbit
Isotype:	IgG
Applications:	Immunohistochemistry (IHC) and Western blot (WB); the recommended starting dilution is 1:1,000-1:2,000 and 1:500-1:1,000, respectively. Other applications were not tested, therefore optimal working concentration/dilution should be determined empirically.

Images

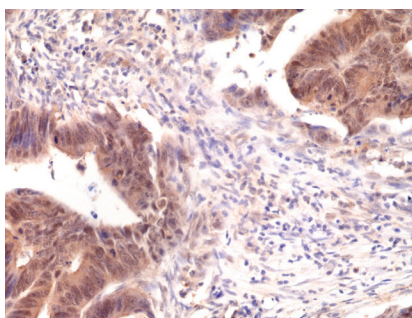


Lane 1: Jurkat cell lysates
Lane 2: IMR32 cell lysates

WB of Jurkat and IMR32 cell lysates using SMAD4 (C-Term) Rabbit Monoclonal Antibody (Clone RM277) at a dilution of 1:500.



Immunohistochemical staining of formalin-fixed and paraffin-embedded human breast cancer tissue using SMAD4 (C-Term) Rabbit Monoclonal Antibody (Clone RM277) at a 1:2,000 dilution.



Immunohistochemical staining of formalin-fixed and paraffin-embedded human colon cancer tissue using SMAD4 (C-Term) Rabbit Monoclonal Antibody (Clone RM277) at a 1:5,000 dilution.

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, 01/29/2024

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

SMAD4, also known as deleted in pancreatic carcinoma locus 4 (DPC4), is a tumor suppressor and a signal transduction protein that functions as a central mediator of the TGF- β and bone morphogenic protein (BMP) signaling pathways.^{1,2} It is comprised of an N-terminal MH1 domain that binds DNA, a linker region containing a C-terminal SMAD activation domain essential to SMAD4 transcriptional activity, and a C-terminal MH2 domain that interacts with the MH1 domain of other SMAD proteins to facilitate homo- and heterodimerization.² SMAD4 is ubiquitously expressed and localized to both the nucleus and cytosol.¹ Upon TGF- β or BMP receptor activation, SMAD4 forms complexes with phosphorylated SMAD2 and SMAD3 or SMAD1, SMAD5, and SMAD8, respectively, which are translocated to the nucleus to induce cell cycle arrest and apoptosis.² Homozygous deletion of, or intragenic inactivation mutations in, SMAD4 have been found in pancreatic and colorectal cancers. Heterozygous mutations in SMAD4 have been found in approximately 20% of patients with juvenile polyposis syndrome and are positively correlated with manifestation of hereditary hemorrhagic telangiectasia.³ Cayman's SMAD4 (C-Term) Rabbit Monoclonal Antibody (Clone RM277) can be used for immunohistochemistry (IHC) and Western blot (WB) applications.

References

1. McCarthy, A.J. and Chetty, R. Smad4/DPC4. *J. Clin. Pathol.* **71(8)**, 661-664 (2018).
2. Zhao, M., Mishra, L., and Deng, C.-X. The role of TGF- β /SMAD4 signaling in cancer. *Int. J. Biol. Sci.* **14(2)**, 111-123 (2018).
3. Andrabi, S., Bekheirnia, M.R., Robbins-Furman, P., et al. SMAD4 mutation segregating in a family with juvenile polyposis, aortopathy, and mitral valve dysfunction. *Am. J. Med. Genet. A.* **155A(5)**, 1165-1169 (2011).

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