

PRODUCT INFORMATION

SMAD2 Polyclonal Antibody

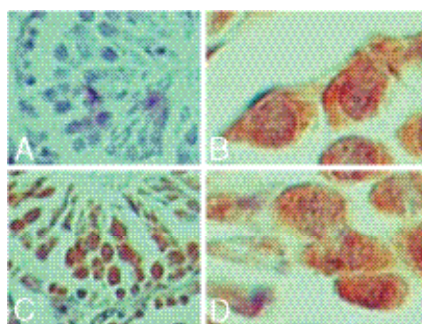
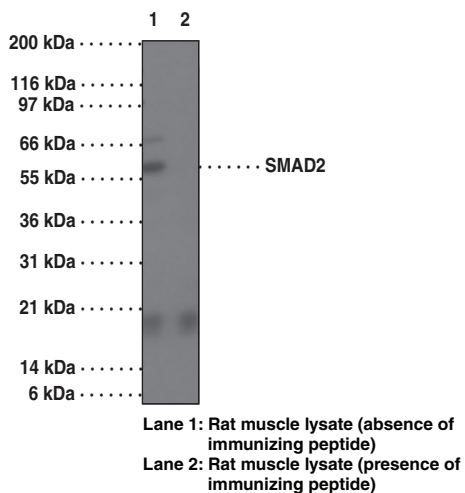
Item No. 10823



Overview and Properties

Contents:	This vial contains 100 µg of protein G-purified IgG
Immunogen:	Human SMAD2
Cross Reactivity:	(+) Human, chicken, mouse, and rat SMAD2
Form:	Liquid
Storage:	4°C (as supplied); Do not freeze
Stability:	≥6 months
Storage Buffer:	200 µl PBS with 0.05% BSA and 0.05% sodium azide.
Concentration:	1.0 mg/ml
Host:	Rabbit
Isotype:	IgG
Applications:	Immunohistochemistry (IHC) and Western blot (WB); the recommended starting concentration for IHC is 5 µg/ml and 2-4 µg/ml for WB.

Images



Immunohistochemical analysis of SMAD2 in formalin-fixed, paraffin-embedded mouse testis tissue using an isotype control (Panel A) and SMAD2 (Panel C, Panel D) at 5 µg/ml.

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, 03/06/2020

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

SMAD2 is an intracellular protein belonging to the Dwarf (DWA/B)/SMAD family. SMAD2 associates with SMAD4 for translocation to the nucleus. It acts as an intracellular mediator of the TGF β family of cytokines and activin type 1 receptor (ACVR1C). Hence it regulates multiple cellular processes like cell growth, proliferation, differentiation, and apoptosis and also cooperates with transcription factors to regulate expression of defined genes. Apart from playing an important role in TGF β 1-induced CTGF expression, SMADs also act as markers of epithelial-mesenchymal transition. SMAD2 also acts as an integrator of multiple signals in the regulation of NOS3 expression. It is ubiquitously expressed in most of tissues and may contribute to colorectal, hepatocellular, lung, and breast carcinomas as well as cervical cancer.

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