

PRODUCT INFORMATION



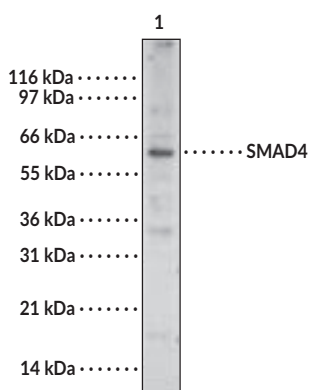
SMAD4 Polyclonal Antibody

Item No. 10838

Overview and Properties

- Contents:** This vial contains 100 µg of protein G-purified IgG in 200 µl PBS, with 0.2% gelatin and 0.05% sodium azide.
- Immunogen:** Synthetic peptides corresponding to amino acids 186-199 and 509-523 of human SMAD4
- Species Reactivity:** (+) Human, mouse, new world monkey, and rat SMAD4
- Storage:** -20°C (as supplied)
- Concentration:** 1.0 mg/ml
- Stability:** ≥6 months
- Host:** Rabbit
- Applications:** Western blot (WB); the recommended starting concentration is 1-3 µg/ml. Other applications were not tested, therefore optimal working concentration/dilution should be determined empirically.

Image



Lane 1: A549 cell lysate (15 µg)

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, 11/18/2024

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

SMADs are a family of intracellular proteins that are essential components in the signaling pathways of the serine/threonine kinase receptors of the transforming growth factor β superfamily.¹ SMADs can be divided into receptor-regulated SMADs (R-SMADs: SMAD1, SMAD2, SMAD3, SMAD5, SMAD8, and SMAD9), common-mediator SMAD (co-SMAD: SMAD4), and inhibitory SMADs (I-SMADs: SMAD6 and SMAD7). SMAD1, SMAD5, SMAD8, and SMAD9 have high degrees of homology and antibodies are available that recognize sequences common to all of them. SMAD8 and SMAD9 are typically used as alternate names for one another in the literature.

Reference

1. Topper, J.N., Cai, J., and Qui, Y. Vascular MADs: Two novel MAD-related genes selectively inducible by flow in human vascular endothelium. *Proc. Natl. Acad. Sci. USA* **94**(17), 9314-9319 (1997).

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