

# PRODUCT INFORMATION



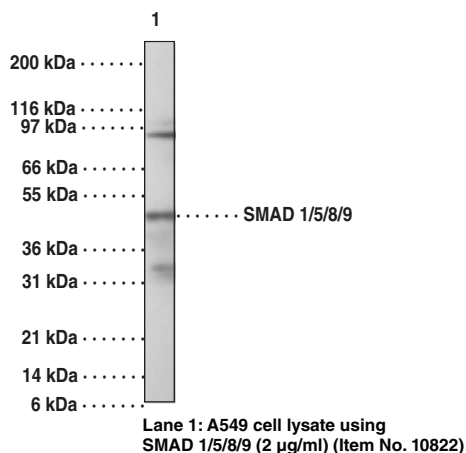
## SMAD1/5/8/9 Polyclonal Antibody

Item No. 10822

### Overview and Properties

<b>Contents:</b>	This vial contains 100 µg of protein G-purified polyclonal antibody.
<b>Synonym:</b>	Mothers Against Decapentaplegic Homolog 1
<b>Immunogen:</b>	Two synthetic peptides corresponding to amino acids 18-33 and amino acids 315-330 of human SMAD1
<b>Cross Reactivity:</b>	(+) Human, mouse, and rat SMAD1/5/8/9; predicted to react with with chicken, chimpanzee, invertebrates, mammals, porcine, <i>Xenopus</i> , and zebrafish
<b>Form:</b>	Liquid
<b>Storage:</b>	-20°C (as supplied); avoid freeze/thaw cycles by aliquoting protein
<b>Stability:</b>	≥9 months
<b>Storage Buffer:</b>	200 µl PBS, with 0.05% BSA and 0.05% sodium azide
<b>Concentration:</b>	0.5 mg/ml
<b>Host:</b>	Rabbit
<b>Isotype:</b>	IgG
<b>Applications:</b>	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



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, 06/22/2021

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 beta superfamily.<sup>1</sup> 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. This antibody is expected to detect SMAD 1, 5, 8, and 9.

## Reference

---

1. Topper, J.N., Cai, J., Qui, Y., *et al.* 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