

# PRODUCT INFORMATION



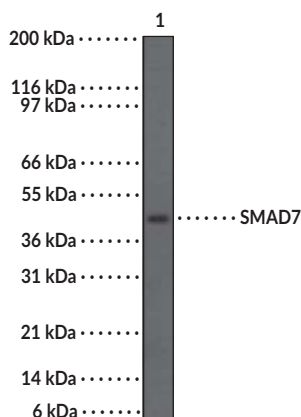
## SMAD7 Polyclonal Antibody

Item No. 10845

### Overview and Properties

<b>Contents:</b>	This vial contains 100 µg of protein G-purified IgG in 200 µl PBS, with 0.05% sodium azide.
<b>Immunogen:</b>	Synthetic peptides corresponding to amino acids 12-29 and 36-50 of human SMAD7
<b>Species Reactivity:</b>	(+) Human, mouse, ovine, and rat SMAD7
<b>Storage:</b>	-20°C (as supplied)
<b>Stability:</b>	≥6 months
<b>Concentration:</b>	1.0 mg/ml
<b>Host:</b>	Rabbit
<b>Applications:</b>	Western blot (WB); the recommended starting concentration is 2-5 µg/ml. Other applications were not tested, therefore optimal working concentration/dilution should be determined empirically.

### Image



Lane 1: HepG2 cell lysate

**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  $\beta$  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.

## 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