

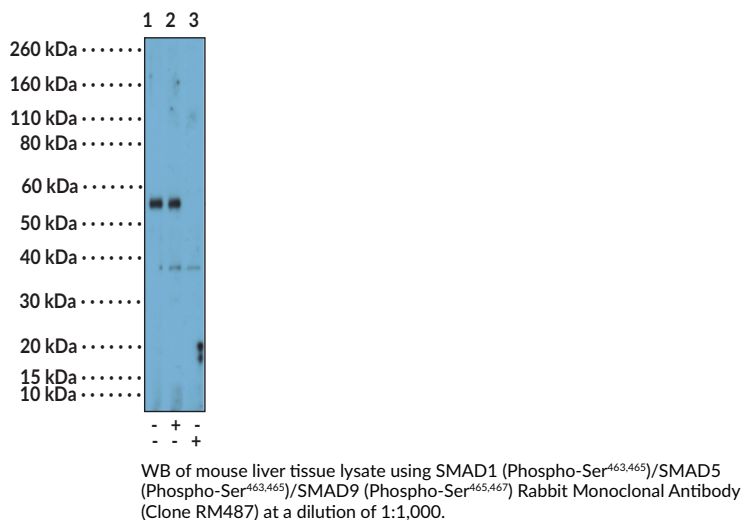
# PRODUCT INFORMATION

**SMAD1 (Phospho-Ser<sup>463,465</sup>)/SMAD5 (Phospho-Ser<sup>463,465</sup>)/SMAD9 (Phospho-Ser<sup>465,467</sup>) Rabbit Monoclonal Antibody (Clone RM487)**  
Item No. 42817

## Overview and Properties

|                            |                                                                                                                                                                                                   |
|----------------------------|---------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------|
| <b>Contents:</b>           | This vial contains 100 µl of protein A-affinity purified monoclonal antibody.                                                                                                                     |
| <b>Immunogen:</b>          | Phosphopeptide corresponding to of SMAD1 (phospho-Ser <sup>463,465</sup> )/SMAD5 (phospho-Ser <sup>463,465</sup> )/SMAD9 (phospho-Ser <sup>465,467</sup> )                                        |
| <b>Cross Reactivity:</b>   | (+) SMAD1 (phospho-Ser <sup>463,465</sup> )/SMAD5 (phospho-Ser <sup>463,465</sup> )/SMAD9 (phospho-Ser <sup>465,467</sup> ); (-) SMAD1/SMAD5/SMAD9 without phosphorylation                        |
| <b>Species Reactivity:</b> | (+) Human, mouse, rat                                                                                                                                                                             |
| <b>Form:</b>               | Liquid                                                                                                                                                                                            |
| <b>Storage:</b>            | -20°C (as supplied)                                                                                                                                                                               |
| <b>Stability:</b>          | ≥1 year                                                                                                                                                                                           |
| <b>Storage Buffer:</b>     | PBS with 50% glycerol, 1% BSA, and 0.09% sodium azide                                                                                                                                             |
| <b>Concentration:</b>      | 1.0 mg/ml                                                                                                                                                                                         |
| <b>Clone:</b>              | RM487                                                                                                                                                                                             |
| <b>Host:</b>               | Rabbit                                                                                                                                                                                            |
| <b>Isotype:</b>            | IgG                                                                                                                                                                                               |
| <b>Application:</b>        | Western blot (WB); the recommended starting dilution is 1:1,000 - 1:2,000. 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, 03/03/2025

**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

SMAD proteins are intracellular transcription factors and mediators of TGF- $\beta$  signaling.<sup>1,3</sup> They are composed of an N-terminal MAD homology 1 (MH1) domain containing nuclear localization and DNA-binding sequences and a C-terminal MH2 domain that is essential for SMAD-receptor interactions, as well as SMAD-SMAD interactions. Binding of TGF- $\beta$  or bone morphogenetic protein (BMP) to their respective cell surface receptors initiates recruitment and activation of SMAD proteins. TGF- $\beta$  induces SMAD2 and SMAD3 activation, while BMP binding induces activation of SMAD1, SMAD5, and SMAD9 via phosphorylation of serine 463 and 465 (Ser<sup>463,465</sup>) in SMAD1 and SMAD5 and Ser<sup>465,467</sup> in SMAD9.<sup>1-3</sup> SMAD1, -5, and -9 activation is associated with hypertrophy in patients with osteoarthritis, but associated with inhibition of fibrotic gene expression in various fibrotic diseases, including chronic kidney disease (CKD).<sup>2,3</sup> Cayman's SMAD1 (Phospho-Ser<sup>463,465</sup>)/SMAD5 (Phospho-Ser<sup>463,465</sup>)/SMAD9 (Phospho-Ser<sup>465,467</sup>) Rabbit Monoclonal Antibody (Clone RM487) can be used for Western blot (WB).

## References

1. Lai, L.Y.S., Gracie, N.P., Gowripalan, A., *et al.* SMAD proteins: Mediators of diverse outcomes during infection. *Eur. J. Cell Biol.* **101**(2), 151204 (2022).
2. Thielen, N.G.M., van Caam, A.P.M., v. Beuningen, H.M., *et al.* Separating friend from foe: Inhibition of TGF- $\beta$ -induced detrimental SMAD1/5/9 phosphorylation while maintaining protective SMAD2/3 signaling in OA chondrocytes. *Osteoarthritis Cartilage* **31**(11), 1481-1490 (2023).
3. Hu, H.H., Chen, D.Q., Wang, Y.N., *et al.* New insights into TGF- $\beta$ /Smad signaling in tissue fibrosis. *Chem. Biol. Interact.* **292**, 76-83 (2018).

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