

# PRODUCT INFORMATION



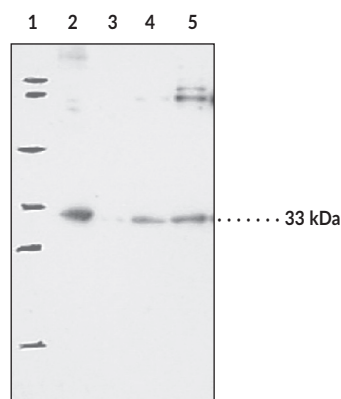
## Prostaglandin E Synthase-2 (microsomal) Polyclonal Antibody

Item No. 160145

### Overview and Properties

<b>Contents:</b>	This vial contains 500 µl of peptide affinity-purified IgG
<b>Synonyms:</b>	Membrane-Associated PGES-2, mPGES-2, mPGE Synthase-2, Prostaglandin H/E Isomerase-2
<b>Immunogen:</b>	Synthetic peptide from the internal region of human mPGES-2
<b>Species Reactivity:</b>	(+) Human, African green monkey, bovine, mouse, ovine, and rat; other species not tested
<b>Cross Reactivity:</b>	(+) mPGES-2; (-) mPGES-1 and cPGES
<b>Uniprot No.:</b>	Q9H7Z7
<b>Form:</b>	Liquid
<b>Storage:</b>	-20°C (as supplied)
<b>Stability:</b>	≥1 year
<b>Storage Buffer:</b>	TBS, pH 7.4, with 50% glycerol, 0.1% BSA, and 0.02% sodium azide
<b>Host:</b>	Rabbit
<b>Applications:</b>	Western blot (WB) and immunohistochemistry (IHC); the recommended starting dilution is 1:200 for WB and 1:500 for IHC. For IHC with formalin-fixed paraffin-embedded tissues the recommended starting dilution is 1:1,000. Other applications were not tested, therefore optimal working concentration/dilution should be determined empirically.

### Image



Lane 1: Protein MW standards  
Lane 2: Mouse heart homogenate (30 µg)  
Lane 3: Bovine aorta homogenate (30 µg)  
Lane 4: Ram seminal vesicle microsomes (30 µg)  
Lane 5: Cos-7 cell lysate (10 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, 04/12/2019

**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

---

Prostaglandin E synthase-2 (microsomal) (mPGES-2) is membrane-associated PGE<sub>2</sub>-producing enzyme.<sup>1,2</sup> Unlike mPGES-1, mPGES-2 does not require glutathione for catalytic activity and is activated in the presence of a broad range of reducing agents. Cayman Chemical's mPGES-2 Polyclonal Antibody can be used for Western blot and Immunohistochemistry (IHC) applications. The antibody recognizes mPGES-2 at 33 kDa from human, mouse, sheep, cow, and monkey samples.

## References

---

1. Watanabe, K., Kurihara, K., and Suzuki, T. Purification and characterization of membrane-bound prostaglandin E synthase from bovine heart. *Biochim. Biophys. Acta* **1439(3)**, 406-414 (1999).
2. Tanikawa, N., Ohmiya, Y., Ohkubo, H., *et al.* Identification and characterization of a novel type of membrane-associated prostaglandin E synthase. *Biochem. Biophys. Res. Commun.* **291(4)**, 884-889 (2002).

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