

# PRODUCT INFORMATION



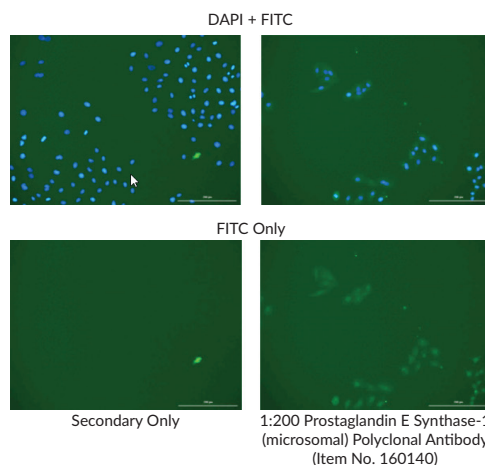
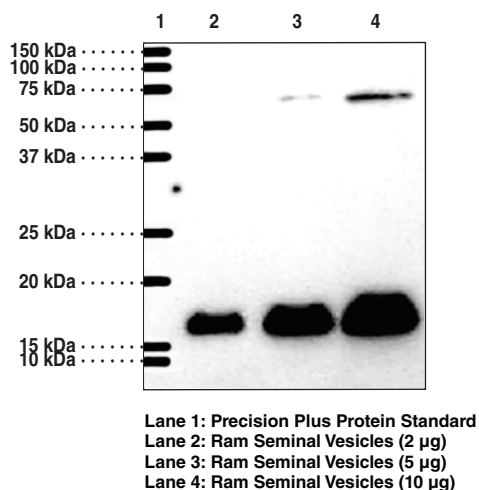
## Prostaglandin E Synthase-1 (microsomal) Polyclonal Antibody

Item No. 160140

### Overview and Properties

<b>Contents:</b>	This vial contains 500 µl of peptide-affinity purified polyclonal antibody.
<b>Synonyms:</b>	Membrane-Associated PGES-1, MGST1-L1, mPGES-1, mPGE Synthase-1, PIG12, Prostaglandin H/E Isomerase
<b>Immunogen:</b>	Peptide from an internal region of human mPGES-1
<b>Species Reactivity:</b>	(+) Human, mouse, ovine, and rat; other species not tested
<b>Uniprot No.:</b>	O14684
<b>Form:</b>	Liquid
<b>Storage:</b>	-20°C (as supplied)
<b>Stability:</b>	≥3 years
<b>Host:</b>	Rabbit
<b>Storage Buffer:</b>	PBS, pH 7.2, with 50% glycerol and 0.02% sodium azide
<b>Application:</b>	Immunofluorescence (IF), Western blot (WB); the recommended starting dilution for is 1:200. Other applications were not tested, therefore optimal working concentration/dilution should be determined empirically.

### Images



Immunofluorescence analysis of paraformaldehyde-fixed, A549 cells. After incubation with Prostaglandin E Synthase-1 (microsomal) Polyclonal Antibody (Item No. 160140), at a 1:200 dilution (or negative control) cells were incubated with FITC labeled anti-rabbit IgG (Item No. 10006588), followed by DAPI nuclear stain. Images show FITC alone or both fluorescence channels to highlight nuclear staining (where applicable).

**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, 11/16/2023

**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-1 (PGES-1) catalyzes the isomerization of  $\text{PGH}_2$  (Item No. 17020) to  $\text{PGE}_2$  (Item No. 14010).  $\text{PGE}_2$  is one of the primary COX products of arachidonic acid (item No. 90010) and one of the most widely investigated PGs. Its activity influences inflammation, fertility and parturition, gastric mucosal integrity, and immune modulation.<sup>1-4</sup> PGES-1 is a 16 kDa membrane-associated protein with glutathione-dependent activity.<sup>5</sup> The mRNA for PGES-1 is expressed in a variety of tissues including prostate, testes, and small intestine, as well as in A549 and HeLa cells.<sup>5</sup> PGES-1 protein expression is increased in A549 cells following treatment with IL-1 $\beta$ .<sup>5</sup>

## References

1. Willis, A.L. and Cornelsen, M. Repeated injection of prostaglandin  $\text{E}_2$  in rat paws induces chronic swelling and a marked decrease in pain threshold. *Prostaglandins* **3**(3), 353-357 (1973).
2. Jackson, G.M., Sharp, H.T., and Varner, M.W. Cervical ripening before induction of labor: A randomized trial of prostaglandin  $\text{E}_2$  gel versus low-dose oxytocin. *Am. J. Obstet. Gynecol.* **171**(4), 1092-1096 (1994).
3. Robert, A., Schultz, J.R., Nezamis, J.E., et al. Gastric antisecretory and antiulcer properties of  $\text{PGE}_2$ , 15-methyl  $\text{PGE}_2$ , and 16,16-dimethyl  $\text{PGE}_2$ . Intravenous, oral and intrajejunal administration. *Gastroenterology* **70**(3), 359-370 (1976).
4. Arvind, P., Papavassiliou, E.D., Tsioulas, G.J., et al. Prostaglandin  $\text{E}_2$  down-regulates the expression of HLA-DR antigen in human colon adenocarcinoma cell lines. *Biochemistry* **34**(16), 5604-5609 (1995).
5. Jakobsson, P.-J., Thorén, S., Morgenstern, R., et al. Identification of human prostaglandin E synthase: A microsomal, glutathione-dependent, inducible enzyme, constituting a potential novel drug target. *Proc. Natl. Acad. Sci. USA* **96**(13), 7220-7225 (1999).

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