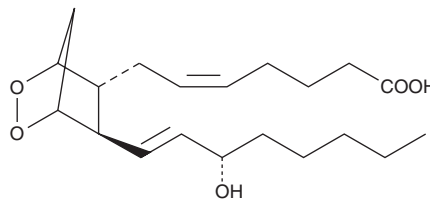


# PRODUCT INFORMATION



## Prostaglandin H<sub>2</sub> Item No. 17020

**CAS Registry No.:** 42935-17-1  
**Formal Name:** 9 $\alpha$ ,11 $\alpha$ -epidioxy-15S-hydroxy-prosta-5Z,13E-dien-1-oic acid  
**Synonyms:** PGH<sub>2</sub>  
**MF:** C<sub>20</sub>H<sub>32</sub>O<sub>5</sub>  
**FW:** 352.5  
**Purity:**  $\geq$ 95%  
**Supplied as:** A solution in acetone  
**Storage:** -80°C  
**Stability:**  $\geq$ 6 months



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

### Laboratory Procedures

Prostaglandin H<sub>2</sub> (PGH<sub>2</sub>) is supplied as a solution in acetone. To change the solvent, simply evaporate the acetone under a gentle stream of nitrogen and immediately add the solvent of choice. PGH<sub>2</sub> is soluble in the organic solvent ethanol at a concentration of approximately 100 mg/ml.

PGH<sub>2</sub> is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, the acetonic solution of PGH<sub>2</sub> should be diluted with the aqueous buffer of choice. The solubility of PGH<sub>2</sub> in PBS (pH 7.2) is approximately 2 mg/ml. We do not recommend storing the aqueous solution for more than one day.

### Description

PGH<sub>2</sub> was first isolated from incubations of arachidonic acid with ovine seminal vesicle microsomes, and was described as a potent vasoconstrictor.<sup>1</sup> PGH<sub>2</sub> is the precursor for all 2-series PGs and thromboxanes (TXs),<sup>2</sup> and is a TP receptor agonist which irreversibly aggregates human platelets at 50-100 ng/ml.

### References

1. Hamberg, M., Svensson, J., Wakabayashi, T., *et al.* Isolation and structure of two prostaglandin endoperoxides that cause platelet aggregation. *Proc. Natl. Acad. Sci. USA* **71**(2), 345-349 (1974).
2. Samuelsson, B., Goldyne, M., Granström, E., *et al.* Prostaglandins and thromboxanes. *Annu. Rev. Biochem.* **47**, 997-1029 (1978).

#### 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, 05/24/2023

#### 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