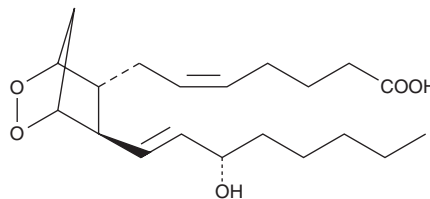


PRODUCT INFORMATION



Prostaglandin H₂ Item No. 17020

CAS Registry No.: 42935-17-1
Formal Name: 9 α ,11 α -epidioxy-15S-hydroxy-prosta-5Z,13E-dien-1-oic acid
Synonym: PGH₂
MF: C₂₀H₃₂O₅
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 H₂ (PGH₂) is a very unstable compound and therefore, stock solutions should be kept on ice, at all times, while performing experiments in the lab. PGH₂ is supplied as a solution in acetone. To change the solvent, first place the vial of PGH₂ on ice. Next, evaporate the acetone under a gentle stream of nitrogen and immediately add the solvent of choice. PGH₂ is soluble in the organic solvent ethanol at a concentration of approximately 100 mg/ml.

PGH₂ is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, the acetic solution of PGH₂ should be diluted with the aqueous buffer of choice. The solubility of PGH₂ in PBS (pH 7.2) is approximately 2 mg/ml. Store aqueous solutions of PGH₂ on ice and use immediately as the half-life of PGH₂ in aqueous solutions is approximately 10 minutes.

Description

PGH₂ was first isolated from incubations of arachidonic acid with ovine seminal vesicle microsomes, and was described as a potent vasoconstrictor.¹ PGH₂ is the precursor for all 2-series PGs and thromboxanes (TXs), and is a TP receptor agonist which irreversibly aggregates human platelets at 50-100 ng/ml.²

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, 10/24/2023

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