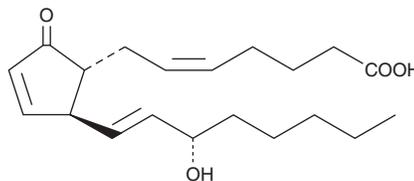


PRODUCT INFORMATION



Prostaglandin A₂ Item No. 10210

CAS Registry No.: 13345-50-1
Formal Name: 15S-hydroxy-9-oxo-prosta-5Z,10,13E-trien-1-oic acid
Synonyms: Medullin, PGA₂
MF: C₂₀H₃₀O₄
FW: 334.5
Purity: ≥98%
UV/Vis.: λ_{max}: 216 nm
Supplied as: A solution in methyl acetate
Storage: -20°C
Stability: ≥2 years



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

Laboratory Procedures

Prostaglandin A₂ (PGA₂) is supplied as a solution in methyl acetate. To change the solvent, simply evaporate the methyl acetate under a gentle stream of nitrogen and immediately add the solvent of choice. Solvents such as ethanol, DMSO, and dimethyl formamide (DMF) purged with an inert gas can be used. The solubility of PGA₂ in these solvents is approximately 100, 50, and 75 mg/ml, respectively.

PGA₂ is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, the methyl acetate solution of PGA₂ should be diluted with the aqueous buffer of choice. The solubility of PGA₂ in PBS (pH 7.2) is approximately 2.4 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

PGA₂ is a naturally occurring prostaglandin in gorgonian corals where it may function in self defense. It is generally not present in mammals. PGA₂ has low biological potency in most bioassays, but it does show some anti-viral/anti-tumor activity.¹ At a 25 μM concentration, PGA₂ blocks the cell cycle progression of NIH 3T3 cells at the G₁ and G₂/M phase.² It has also been shown to act as a vasodilator with natriuretic properties.³

References

1. Fukushima, M., Kato, T., Narumiya, S., *et al.* Prostaglandin A and J: Antitumor and antiviral prostaglandins. *Adv. Prostaglandin Thromboxane Leukotriene Res.* **19**, 415-418 (1989).
2. Hitomi, M., Shu, J., Strom, D., *et al.* Prostaglandin A₂ blocks the activation of G₁ phase cyclin-dependant kinase without altering mitogen-activated protein kinase stimulation. *J. Biol. Chem.* **271(16)**, 9376-9383 (1996).
3. Frolich, J.C., Sweetman, B.J., Carr, K., *et al.* Assessment of the levels of PGA₂ in human plasma by gas chromatography-mass spectrometry. *Prostaglandins* **10(1)**, 185-195 (1975).

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, 02/29/2024

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