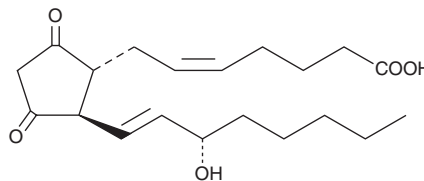


PRODUCT INFORMATION



Prostaglandin K₂ Item No. 18900

CAS Registry No.: 275816-51-8
Formal Name: 9,11-dioxo-15S-hydroxy-prosta-5Z,13E-dien-1-oic acid
Synonym: PGK₂
MF: C₂₀H₃₀O₅
FW: 350.5
Purity: ≥98%
UV/Vis.: λ_{max}: 245 nm
Supplied as: A solution in methyl acetate
Storage: -20°C
Stability: ≥1 year



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

Laboratory Procedures

Prostaglandin K₂ (PGK₂) 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 purged with an inert gas can be used. The solubility of PGK₂ in these solvents is approximately 75, 50, and 100 mg/ml, respectively.

PGK₂ is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, the methyl acetate solution of PGK₂ should be diluted with the aqueous buffer of choice. The solubility of PGK₂ in PBS (pH 7.2) is approximately 2.7 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

PGK₂ is the 9,11-diketone formed by the oxidation of PGE₂ or PGD₂. Whether this compound exists biologically is uncertain; it is known to be resistant to metabolism by 15-hydroxy PGDH *in vitro*.¹

Reference

1. Ohno, H., Morikawa, Y., and Hirata, F. Studies on 15-hydroxyprostaglandin dehydrogenase with various prostaglandin analogues. *J. Biochem.* **84**(6), 1485-1494 (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, 05/15/2023

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