

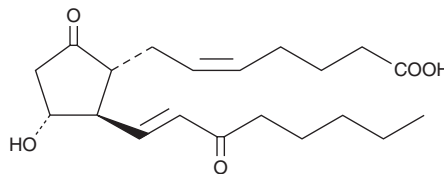
PRODUCT INFORMATION



15-keto Prostaglandin E₂

Item No. 14720

CAS Registry No.: 26441-05-4
Formal Name: 9,15-dioxo-11 α -hydroxy-prosta-5Z,13E-dien-1-oic acid
Synonyms: 15-keto PGE₂, 15-oxo PGE₂
MF: C₂₀H₃₀O₅
FW: 350.5
Purity: \geq 98%
UV/Vis.: λ_{max} : 229 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: \geq 4 years



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

Laboratory Procedures

15-keto Prostaglandin E₂ (15-keto PGE₂) is supplied as a crystalline solid. 15-keto PGE₂ is sparingly soluble in water but freely soluble in organic solvents such as ethanol, DMSO, or dimethyl formamide. The solubility of 15-keto PGE₂ in these solvents is at least 100 mg/ml.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of 15-keto PGE₂ can be prepared by directly dissolving the crystalline compound in aqueous buffers. The solubility of 15-keto PGE₂ in PBS, pH 7.2, is approximately 5 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

15-keto PGE₂ is a metabolite of PGE₂ (Item No. 14010) formed by 15-hydroxy prostaglandin dehydrogenase (15-PGDH).¹ Unlike PGE₂, 15-keto PGE₂ does not bind effectively to the PGE₂ receptors EP₂ and EP₄ expressed in CHO cells (K_is = 2.6 and 15 μ M, respectively) or induce adenylate cyclase activity in the same cells (EC₅₀s = 1.8 and >33 μ M, respectively). However, it does bind to EP₂ and EP₄ in HEK cells expressing these receptors (IC₅₀s = 0.117 and 2.82 μ M, respectively), as well as induces cAMP formation (EC₅₀s = 0.137 and 0.426 μ M, respectively) and the transcriptional activity of β -catenin/TCF in the same cells.² 15-keto PGE₂ inhibits CD3-CD28-MHC-I-induced proliferation of isolated human CD4⁺ T cells in a concentration-dependent manner.³ It also reduces mortality in a mouse model of LPS-induced sepsis when administered at a dose of 15 mg/kg.⁴

References

1. Nishigaki, N., Negishi, M., and Ichikawa, A. Two G_s-coupled prostaglandin E receptor subtypes, EP2 and EP4, differ in desensitization and sensitivity to the metabolic inactivation of the agonist. *Mol. Pharmacol.* **50**(4), 1031-1037 (1996).
2. Endo, S., Suganami, A., Fukushima, K., *et al.* 15-Keto-PGE₂ acts as a biased/partial agonist to terminate PGE₂-evoked signaling. *J. Biol. Chem.* **295**(38), 13338-13352 (2020).
3. Schmidleithner, L., Thabet, Y., Schönfeld, E., *et al.* Enzymatic activity of HPGD in Treg cells suppresses Tconv cells to maintain adipose tissue homeostasis and prevent metabolic dysfunction. *Immunity* **50**(5), 1232-1248 (2019).
4. Chen, I.-J., Hee, S.-W., Liao, C.-H., *et al.* Targeting the 15-keto-PGE₂-PTGR2 axis modulates systemic inflammation and survival in experimental sepsis. *Free Radic. Biol. Med.* **115**, 113-126 (2018).

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, 01/16/2024

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