

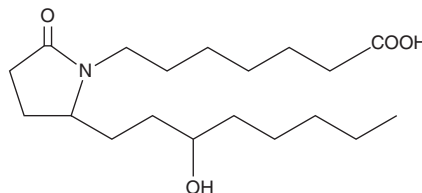
PRODUCT INFORMATION



CAY10580

Item No. 16835

CAS Registry No.: 64054-40-6
Formal Name: 2-(3-hydroxyoctyl)-5-oxo-1-pyrrolidineheptanoic acid
MF: C₁₉H₃₅NO₄
FW: 341.5
Purity: ≥96%
Supplied as: A solution in ethanol
Storage: -20°C
Stability: ≥2 years



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

Laboratory Procedures

CAY10580 is supplied as a solution in ethanol. To change the solvent, simply evaporate the ethanol under a gentle stream of nitrogen and immediately add the solvent of choice. Solvents such as DMSO and dimethyl formamide purged with an inert gas can be used. The solubility of CAY10580 in these solvents is approximately 10 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. If an organic solvent-free solution of CAY10580 is needed, it can be prepared by evaporating the ethanol and directly dissolving the neat oil in aqueous buffers. The solubility of CAY10580 in PBS (pH 7.2) is approximately 2.5 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Prostaglandin E₂ (PGE₂) activates four E prostanoid (EP) receptors, EP₁₋₄. EP₄ is a G_s protein-coupled receptor that, by elevating the second messenger cAMP, plays important roles in bone formation and resorption, cancer, and atherosclerosis.¹⁻³ CAY10580 is an 8-aza-9-oxo-15-hydroxy saturated analog of PGE₂. It selectively binds the EP₄ receptor (K_i = 35 nM) relative to the EP₁, EP₂, and EP₃ receptors (K_i = 3,000, 2,000, and >3,000 nM, respectively).⁴ CAY10580 stimulates cAMP formation in excised mouse ovaries.⁵

References

1. Li, M., Thompson, D.D., and Paralkar, V.M. Prostaglandin E₂ receptors in bone formation. *Int. Orthop.* **31(6)**, 767-772 (2007).
2. Hawcroft, G., Ko, C.W.S., and Hull, M.A. Prostaglandin E₂-EP₄ receptor signalling promotes tumorigenic behaviour of HT-29 human colorectal cancer cells. *Oncogene* **26(21)**, 3006-3019 (2007).
3. Babaev, V.R., Chew, J.D., Ding, L., et al. Macrophage EP₄ deficiency increases apoptosis and suppresses early atherosclerosis. *Cell Metab.* **8(6)**, 492-501 (2008).
4. Billot, X., Chateauneuf, A., Chauret, N., et al. Discovery of a potent and selective agonist of the prostaglandin EP₄ receptor. *Bioorg. Med. Chem. Lett.* **13(6)**, 1129-1132 (2003).
5. Smith, R.L., Lee, T., Gould, N.P., et al. Prostaglandin isosteres. 1. (8-Aza-, 8,10-Diaza-, and 8-Aza-11-thia)-9-oxoprostanoic acids and their derivatives. *J. Med. Chem.* **20(10)**, 1292-1299 (1977).

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, 02/02/2024

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