

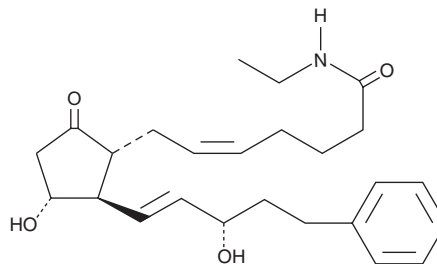
PRODUCT INFORMATION



17-phenyl trinor Prostaglandin E₂ ethyl amide

Item No. 13532

CAS Registry No.: 1219032-20-8
Formal Name: N-ethyl-9-oxo-11 α ,15S-dihydroxy-17-phenyl-18,19,20-trinor-prosta-5Z,13E-dien-1-amide
Synonym: 17-phenyl trinor PGE₂ ethyl amide
MF: C₂₅H₃₅NO₄
FW: 413.6
Purity: ≥98%
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

17-phenyl trinor Prostaglandin E₂ ethyl amide (17-phenyl trinor PGE₂ ethyl amide) 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 17-phenyl trinor PGE₂ ethyl amide in these solvents is approximately 100 mg/ml.

17-phenyl trinor PGE₂ ethyl amide is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, the ethanolic solution of 17-phenyl trinor PGE₂ ethyl amide should be diluted with the aqueous buffer of choice. The solubility of 17-phenyl trinor PGE₂ ethyl amide in PBS (pH 7.2) is approximately 0.5 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

17-phenyl trinor PGE₂ ethyl amide is derived from 17-phenyl trinor PGE₂, a synthetic analog of PGE₂ that acts as an agonist of EP₁ and EP₃ receptors in mice (K_i = 14 and 3.7 nM, respectively) and EP₁, EP₃, and EP₄ in rats (K_i = 25, 4.3, and 54 nM, respectively).^{1,2} 17-phenyl trinor PGE₂ causes contraction of guinea pig ileum at a concentration of 11 μ M and is 4.4 times more potent than PGE₂ as an antifertility agent in hamsters.^{1,2} Modification of the C-1 carboxyl group to an ethyl amide serves to increase lipid solubility, thereby improving uptake into tissues and further lowering the effective concentration. Ethyl amide groups are then removed by amidases, regenerating the active free acid.

References

1. Kiriya, M., Ushikubi, F., Kobayashi, T., *et al.* Ligand binding specificities of the eight types and subtypes of the mouse prostanoid receptors expressed in Chinese hamster ovary cells. *Br. J. Pharmacol.* **122(2)**, 217-224 (1997).
2. Boie, Y., Stocco, R., Sawyer, N., *et al.* Molecular cloning and characterization of the four rat prostaglandin E₂ prostanoid receptor subtypes. *Eur. J. Pharmacol.* **340(2-3)**, 227-241 (1997).
3. Lawrence, R.A., Jones, R.L., and Wilson, N.H. Characterization of receptors involved in the direct and indirect actions of prostaglandins E and I on the guinea-pig ileum. *Br. J. Pharmacol.* **105(2)**, 271-278 (1992).
4. Miller, W.L., Weeks, J.R., Lauderdale, J.W., *et al.* Biological activities of 17-phenyl-18,19,20-trinorprostaglandins. *Prostaglandins* **9(1)**, 9-18 (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, 02/23/2024

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