

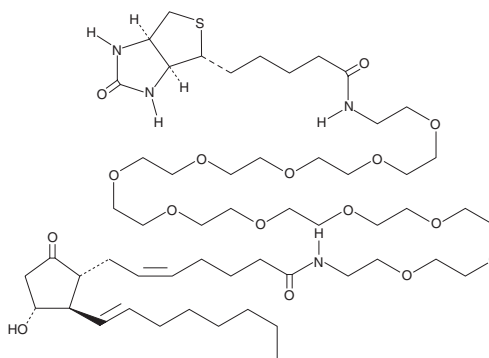
PRODUCT INFORMATION



Prostaglandin E₂-PEG₁₁-biotin

Item No. 9000376

Formal Name: 11 α ,15S-dihydroxy-9-oxo-prosta-52,13E-dien-1-oyl-polyethylene glycol 11-N'-biotinoyl-1,5-diaminopentane
Synonym: PGE₂-PEG₁₁-biotin
MF: C₅₄H₉₆N₄O₁₇S
FW: 1,105.4
Purity: \geq 98%
Supplied as: A solution in ethanol
Storage: -20°C
Stability: \geq 2 years



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

Laboratory Procedures

Prostaglandin E₂-PEG₁₁-biotin (PGE₂-PEG₁₁-biotin) 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 ethanol, DMSO, and dimethyl formamide purged with an inert gas can be used. The solubility of PGE₂-PEG₁₁-biotin in these solvents is approximately 50 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 PGE₂-PEG₁₁-biotin is needed, it can be prepared by evaporating the ethanol and directly dissolving the neat oil in aqueous buffers. The solubility of PGE₂-PEG₁₁-biotin in PBS (pH 7.2) is approximately 1 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

PGE₂ is the major eicosanoid product of cyclooxygenase-1 and 2 in epithelial cells, leukocytes, reproductive tissues, and many others. PGE₂ contracts smooth muscle, protects the gastric mucosa, modulates immune function, and regulates fertility and implantation. PGE₂ acts through at least four G-protein coupled receptors, EP₁-EP₄. PGE₂-PEG₁₁-biotin is an affinity probe which allows PGE₂ to be detected or immobilized using the biotin ligand. The PEG₁₁ moiety serves to separate the biotin linker from PGE₂ with a hydrophilic spacer. It is thus a tool to be used in the elucidation of PGE₂ actions or interactions in aqueous solutions.

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, 02/07/2024

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