

PRODUCT INFORMATION



Prostaglandin D₂ Dopamine-d₄ Item No. 19104

Formal Name: 9 α ,15S-dihydroxy-11-oxo-prosta-5Z,13E-dien-1-oic N-[2-(3,4-dihydroxyphenyl)ethyl-d₄] amide

Synonyms: PGD₂DA-d₄, PGD₂ Dopamine-d₄

MF: C₂₈H₃₇D₄NO₆

FW: 491.7

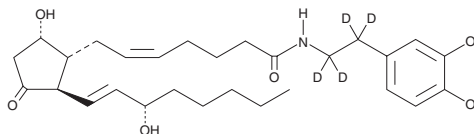
Chemical Purity: \geq 98% Prostaglandin D₂ Dopamine Deuterium

Incorporation: \geq 99% deuterated forms (d₁-d₄); \leq 1% d₀

Stability: \geq 1 year at -20°C

Supplied as: A solution in ethanol

UV/Vis.: λ_{max} : 283 nm



Laboratory Procedures

Prostaglandin D₂ dopamine-d₄ (PGD₂ dopamine-d₄) contains four deuterium atoms. It is intended for use as an internal standard for the quantification of PGD₂ dopamine (Item No. 9002504) by GC- or LC-mass spectrometry (MS). For long term storage, we suggest that PGD₂ dopamine-d₄ be stored as supplied at -20°C. It should be stable for at least one year.

PGD₂ dopamine-d₄ is supplied as a solution in ethanol. To change the solvent, simply evaporate the PGD₂ dopamine-d₄ under a gentle stream of nitrogen and immediately add the solvent of choice. Solvents such as ethanol, DMSO, and dimethyl formamide (DMF) purged with an inert gas can be used. The solubility of PGD₂ dopamine-d₄ in these solvents is approximately 50, 20, and 30 mg/ml respectively.

PGD₂ dopamine-d₄ is used as an internal standard for the quantification of PGD₂ dopamine by stable isotope dilution MS. The accuracy of the sample weight in this vial is between 5% over and 2% under the amount shown on the vial. If better precision is required, the deuterated standard should be quantitated against a more precisely weighed unlabeled standard by constructing a standard curve of peak intensity ratios (deuterated *versus* unlabeled).

Description

PGD₂ dopamine is a conjugate of the neurotransmitter dopamine and PGD₂ that can be used to study the biological function of PGD₂ in the brain and periphery.

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, 02/26/2016

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