

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)

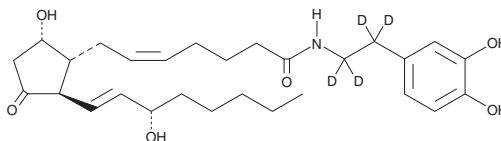
Deuterium Incorporation: $\geq 99\%$ deuterated forms (d₁-d₄); $\leq 1\%$ d₀

UV/Vis.: λ_{max} : 283 nm

Supplied as: A 100 $\mu\text{g/ml}$ solution in ethanol

Storage: -20°C

Stability: ≥ 1 year



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

Laboratory Procedures

Prostaglandin D₂ dopamine-d₄ (PGD₂ dopamine-d₄) is intended for use as an internal standard for the quantification of PGD₂ dopamine (Item No. 9002504) by GC- or LC-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).

PGD₂ dopamine-d₄ 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 PGD₂ dopamine-d₄ in these solvents is approximately 20 and 30 mg/ml respectively.

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, 01/05/2026

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