

PRODUCT INFORMATION



AR-102-d₄
Item No. 10009682

Formal Name: 3-hydroxy-2,2-bis(hydroxymethyl)propyl
7-((1R,2R,3R,5S)-2-((R)-3-(benzo[b]
thiophen-2-yl)-3-hydroxypropyl)-3,5-
dihydroxycyclopentyl)heptanoate-3,3,4,4-d₄

Synonym: Dukeprost-d₄ pentaerythritol ester

MF: C₂₈H₃₈D₄O₈S

FW: 542.7

Chemical Purity: ≥95% (AR-102)

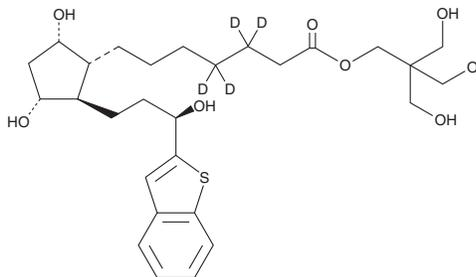
Deuterium

Incorporation: ≥99% deuterated forms (d₁-d₄); ≤1% d₀

Supplied as: A solution in methanol

Storage: -20°C

Stability: ≥2 years



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

Laboratory Procedures

AR-102-d₄ is intended for use as an internal standard for the quantification of AR-102 (Item No. 9000362) 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).

Description

AR-102 is an FP receptor agonist, a derivative of prostaglandin F (PGF), and a prodrug form of AR-102 (free acid) (Item Nos. 10008049 | 43439).^{1,2}

References

1. Fogagnolo, P. and Rossetti, L. Medical treatment of glaucoma: Present and future. *Expert Opin. Investig. Drugs* **20(7)**, 947-959 (2011).
2. Wierzbowska, J., Robaszkiewicz, J., Figurska, M., *et al.* Future possibilities in glaucoma therapy. *Med. Sci. Monit.* **16(11)**, RA252-RA259 (2010).

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, 05/08/2025

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