

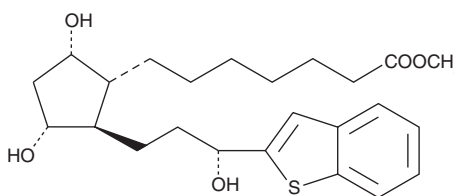
PRODUCT INFORMATION



AR-102 methyl ester

Item No. 9000700

CAS Registry No.: 1245699-42-6
Formal Name: (1R,2R,3R,5S)-2-[(3R)-3-benzo[b]thien-2-yl-3-hydroxypropyl]-3,5-dihydroxy-cyclopentaneheptanoic acid, methyl ester
Synonym: Dukeprost methyl ester
MF: C₂₄H₃₄O₅S
FW: 434.6
Purity: ≥98%
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥4 years



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

Laboratory Procedures

AR-102 methyl ester is supplied as a crystalline solid. A stock solution may be made by dissolving the AR-102 methyl ester in the solvent of choice, which should be purged with an inert gas. AR-102 methyl ester is soluble (≥10 mg/ml) in DMSO and sparingly soluble (1-10 mg/ml) in ethanol.

Description

AR-102 methyl ester is a derivative of AR-102 (Item No. 10008049), a prostaglandin F derivative.

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, 04/30/2025

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