

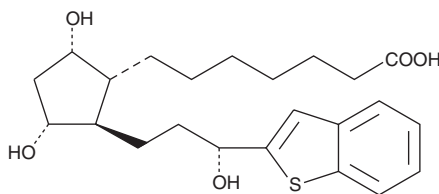
PRODUCT INFORMATION



AR-102 (free acid)

Item No. 10008049

CAS Registry No.: 955007-30-4
Formal Name: 7-((1R,2R,3R,5S)-2-((R)-3-(benzo[b]thiophen-2-yl)-3-hydroxypropyl)-3,5-dihydroxycyclopentyl)heptanoic acid
Synonym: Dukeprost
MF: C₂₃H₃₂O₅S
FW: 420.6
Purity: ≥98%
Supplied as: A 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 (free acid) is supplied as a crystalline solid. A stock solution may be made by dissolving the AR-102 (free acid) in the solvent of choice, which should be purged with an inert gas. AR-102 (free acid) is soluble (≥10 mg/ml) in ethanol and DMSO.

Description

AR-102 (free acid) is a derivative of prostaglandin F (PGF) and an active metabolite of AR-102 (Item No. 9000362).

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, 05/07/2025

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