

PRODUCT INFORMATION



AR-102 free acid (sodium salt)

Item No. 43439

Formal Name: (1R,2R,3R,5S)-2-[(3R)-3-benzo[b]thien-2-yl-3-hydroxypropyl]-3,5-dihydroxy-cyclopentaneheptanoic acid, monosodium salt

Synonym: Dukeprost

MF: $C_{23}H_{31}O_5S \cdot Na$

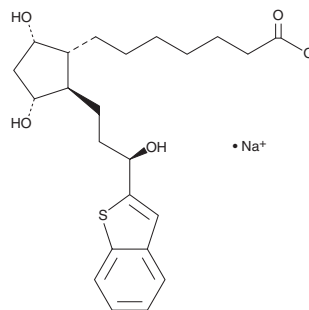
FW: 442.5

Purity: $\geq 98\%$

Supplied as: A solid

Storage: $-20^\circ C$

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

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of AR-102 free acid (sodium salt) can be prepared by directly dissolving the solid in aqueous buffers. AR-102 free acid (sodium salt) is soluble (≥ 10 mg/ml) in PBS (pH 7.2). We do not recommend storing the aqueous solution for more than one day.

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