

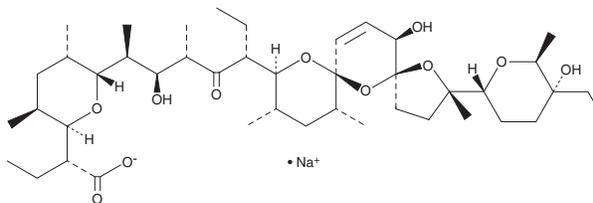
PRODUCT INFORMATION



Narasin (sodium salt)

Item No. 19447

CAS Registry No.: 58331-17-2
Formal Name: 4S-methyl-salinomycin, monosodium salt
Synonyms: 4-Methylsalinomycin, Monteban, Narasin A
MF: C₄₃H₇₁O₁₁ • Na
FW: 787.0
Purity: ≥95%
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

Narasin (sodium salt) is supplied as a solid. A stock solution may be made by dissolving the narasin (sodium salt) in the solvent of choice, which should be purged with an inert gas. Narasin (sodium salt) is soluble in organic solvents such as ethanol, methanol, DMSO, and dimethyl formamide.

Narasin (sodium salt) is sparingly soluble in aqueous solutions. To enhance aqueous solubility, dilute the organic solvent solution into aqueous buffers or isotonic saline. If performing biological experiments, ensure the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. We do not recommend storing the aqueous solution for more than one day.

Description

Narasin is an ionophore antibiotic isolated from certain *Streptomyces* sp. and used in veterinary practice as a coccidiostat for gastrointestinal parasites. It has also been shown to inhibit NF-κB signaling via inhibition of IκBα phosphorylation (IC₅₀ = 3.2 μM) and to stimulate tumor necrosis factor-related apoptosis-induced ligand (TRAIL)-mediated apoptosis in glioma cells via ER stress.^{1,2}

References

1. Miller, S.C., Huang, R., Sakamuru, S., *et al.* Identification of known drugs that act as inhibitors of NF-κB signaling and their mechanism of action. *Biochem. Pharmacol.* **79(9)**, 1272-1280 (2016).
2. Yoon, M.J. and Kang, Y.J. Monensin, a polyether ionophore antibiotic, overcomes TRAIL resistance in glioma cells via endoplasmic reticulum stress, DR5 upregulation and c-FLIP downregulation. *Carcinogenesis* **34(8)**, 1918-1928 (2013).

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, 11/17/2022

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