

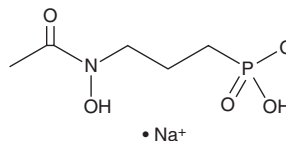
PRODUCT INFORMATION



FR900098 (sodium salt)

Item No. 10012681

CAS Registry No.: 73226-73-0
Formal Name: P-[3-(acetylhydroxyamino)propyl]-phosphonic acid, monosodium salt
MF: C₅H₁₁NO₅P • Na
FW: 219.1
Purity: ≥95%
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥4 years
Item Origin: Synthetic



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

Laboratory Procedures

FR900098 (sodium salt) is supplied as a crystalline solid. Aqueous solutions of FR900098 (sodium salt) can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of FR900098 (sodium salt) in PBS, pH 7.2, is approximately 5 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

FR900098 is a derivative of fosmidomycin (Item No. 16000) that has antimalarial activity (IC₅₀s = 170, 170, and 90 nM for HB3, A2, and Dd2 *P. falciparum* strains, respectively).¹ FR900098 inhibits DOXP reductoisomerase, a key enzyme in the non-mevalonate pathway of isoprenoid biosynthesis, *in vitro* in a dose-dependent manner.¹ FR900098 (>10 mg/kg, i.p.) eradicated *P. vinckei* parasites within four days in mice.

Reference

1. Jomaa, H., Wiener, J., Sanderbrand, S., *et al.* Inhibitors of the nonmevalonate pathway of isoprenoid biosynthesis as antimalarial drugs. *Science* **285(5433)**, 1573-1576 (1999).

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, 12/12/2022

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