

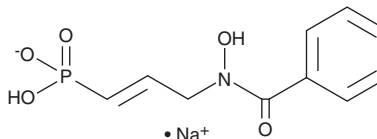
PRODUCT INFORMATION



DXR Inhibitor 11a (sodium salt)

Item No. 39305

Formal Name: (E)-hydrogen (3-(N-hydroxybenzamido)prop-1-en-1-yl)phosphonate, monosodium salt
Synonym: 1-Deoxy-D-Xylulose-5-Phosphate Reductoisomerase Inhibitor 11a
MF: C₁₀H₁₁NO₅P • Na
FW: 279.2
Purity: ≥98%
Supplied as: A solution in methanol
Storage: -20°C
Stability: ≥1 year



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

Description

DXR inhibitor 11a is an inhibitor of *P. falciparum* 1-deoxy-D-xylulose-5-phosphate reductoisomerase (DXR; IC₅₀ = 0.11 μM).¹ It inhibits the growth of *P. falciparum* in isolated human erythrocytes (IC₅₀ = 0.37 μM).

Reference

1. Wang, X., Edwards, R.L., Ball, H.S., et al. MEPicides: α,β-Unsaturated fosmidomycin N-acyl analogs as efficient inhibitors of *Plasmodium falciparum* 1-deoxy-D-xylulose-5-phosphate reductoisomerase. *ACS Infect. Dis.* **9**(7), 1387-1395 (2023).

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, 10/11/2023

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