

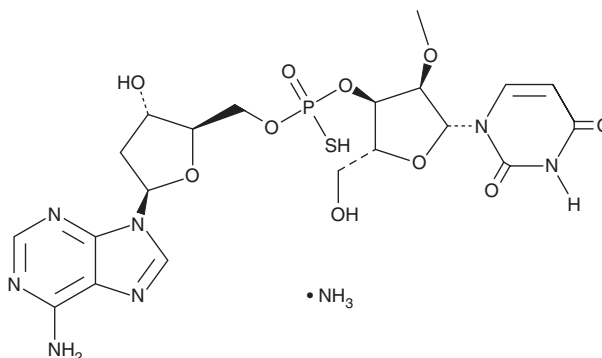
PRODUCT INFORMATION



Inarigivir (ammonium salt)

Item No. 42216

CAS Registry No.: 2172788-92-8
Formal Name: 2'-O-methyl-P-thiouridylyl-(3'→5')-2'-deoxy-adenosine, monoammonium salt
Synonym: ORI-9020
MF: C₂₀H₂₆N₇O₁₀PS • NH₃
FW: 604.5
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

Inarigivir (ammonium salt) is supplied as a solid. A stock solution may be made by dissolving the inarigivir (ammonium salt) in the solvent of choice, which should be purged with an inert gas. Inarigivir (ammonium salt) is sparingly soluble (1-10 mg/ml) in 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 inarigivir (ammonium salt) can be prepared by directly dissolving the solid in aqueous buffers. Inarigivir (ammonium 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

Inarigivir is a dinucleotide antiviral agent.¹ It decreases extracellular hepatitis B virus (HBV) DNA and intracellular HBV RNA levels in HepG2 2.2.15 cells (EC₅₀s = 0.258 and 1.4 μM, respectively) without inducing cytotoxicity with a 50% cytotoxic concentration (CC₅₀) value of >1,000 μM. Inarigivir (100 mg/kg per day) reduces hepatic HBV DNA levels in HBV-expressing transgenic mice.²

References

1. Iyer, R.P., Jin, Y., Roland, A., *et al.* Phosphorothioate di- and trinucleotides as a novel class of anti-hepatitis B virus agents. *Antimicrob. Agents Chemother.* **48(6)**, 2199-2205 (2004).
2. Iyer, R.P., Roland, A., Jin, Y., *et al.* Anti-hepatitis B virus activity of ORI-9020, a novel phosphorothioate dinucleotide, in a transgenic mouse model. *Antimicrob. Agents Chemother.* **48(6)**, 2318-2320 (2004).

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, 01/28/2025

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