

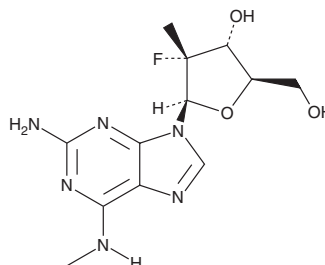
PRODUCT INFORMATION



HCV-IN-31

Item No. 36778

CAS Registry No.: 1998705-62-6
Formal Name: (2'R)-2-amino-2'-deoxy-2'-fluoro-N,2'-dimethyl-adenosine
Synonym: Hepatitis C Virus Inhibitor 31
MF: C₁₂H₁₇FN₆O₃
FW: 312.3
Purity: ≥98%
UV/Vis.: λ_{max}: 224, 282 nm
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

HCV-IN-31 is supplied as a solid. A stock solution may be made by dissolving the HCV-IN-31 in the solvent of choice, which should be purged with an inert gas. HCV-IN-31 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of HCV-IN-31 in DMSO and DMF is approximately 1 mg/ml. HCV-IN-31 is slightly soluble in ethanol.

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 HCV-IN-31 can be prepared by directly dissolving the solid in aqueous buffers. The solubility of HCV-IN-31 in PBS (pH 7.2) is approximately 0.2 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

HCV-IN-31 is a nucleoside with antiviral activity.^{1,2} It inhibits hepatitis C virus (HCV) replication in an HCV replicon assay (EC₅₀ = 15.7 μM).²

References

1. Good, S.S., Moussa, A., Zhou, X.-J., *et al.* Preclinical evaluation of AT-527, a novel guanosine nucleotide prodrug with potent, pan-genotypic activity against hepatitis C virus. *PLoS One* **15**(1), e0227104 (2020).
2. Sommadossi, J.-P. and Moussa, A. Beta-D-2'-deoxy-2'-alpha-fluoro-2'-beta-C-substituted-2'-modified-N6-substituted purine nucleotides for HCV treatment. *Atea Pharmaceuticals, Inc.* **US20160257706A1** (2016).

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, 12/21/2022

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