

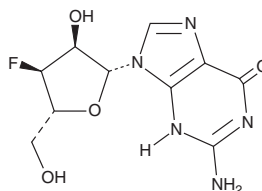
PRODUCT INFORMATION



3'-Deoxy-3'-fluoroguanosine

Item No. 35699

CAS Registry No.: 123402-21-1
MF: $C_{10}H_{12}FN_5O_4$
FW: 285.2
Purity: $\geq 98\%$
Supplied as: A solid
Storage: $-20^{\circ}C$
Stability: ≥ 4 years



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

Laboratory Procedures

3'-Deoxy-3'-fluoroguanosine is supplied as a solid. A stock solution may be made by dissolving the 3'-deoxy-3'-fluoroguanosine in the solvent of choice, which should be purged with an inert gas. 3'-Deoxy-3'-fluoroguanosine is soluble in DMSO and slightly soluble in acetonitrile and methanol.

3'-Deoxy-3'-fluoroguanosine is slightly 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

3'-Deoxy-3'-fluoroguanosine is a nucleoside analog and an inhibitor of hepatitis C virus (HCV) RNA polymerase non-structural protein 5B (NS5B; $IC_{50} = 1.8 \mu M$).¹

Reference

1. Eldrup, A.B., Allerson, C.R., Bennett, C.F., *et al.* Structure-activity relationship of purine ribonucleosides for inhibition of hepatitis C virus RNA-dependent RNA polymerase. *J. Med. Chem.* **47**(9), 2283-2295 (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, 02/28/2024

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