

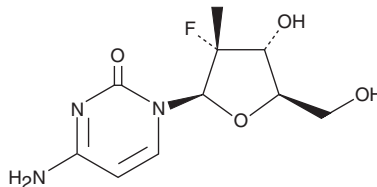
PRODUCT INFORMATION



PSI-6130

Item No. 13575

CAS Registry No.: 817204-33-4
Formal Name: (2'R)-2'-deoxy-2'-fluoro-2'-methyl-cytidine
MF: C₁₀H₁₄FN₃O₄
FW: 259.2
Purity: ≥98%
UV/Vis.: λ_{max}: 273 nm
Supplied as: A crystalline 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

PSI-6130 is supplied as a crystalline solid. A stock solution may be made by dissolving the PSI-6130 in the solvent of choice, which should be purged with an inert gas. PSI-6130 is soluble in the organic solvent DMSO at a concentration of approximately 50 mg/ml.

Description

PSI-6130 is a nucleoside inhibitor of the NS5B RNA polymerase of hepatitis C virus (HCV; $K_i = 4.3 \mu\text{M}$).¹ It inhibits HCV genotype 1b (GT-1b) Con1 and GT-1a H77 viral replication in Huh7 replicon cells using a luciferase-based assay ($\text{EC}_{50\text{s}} = 0.51$ and $0.30 \mu\text{M}$, respectively). PSI-6130 also inhibits replication of HCV GT-1b and GT-1a replicons from clinical isolates, with EC_{50} values ranging from 0.60 to $1.41 \mu\text{M}$, and 0.20 to $0.43 \mu\text{M}$, respectively, in the same assay. It has little or no activity against West Nile Virus (WNV), Dengue type 2 virus (DV), human immunodeficiency virus (HIV), or hepatitis B virus (HBV; $\text{EC}_{90\text{s}} = 46.3, >100, >100, \text{ and } >10 \mu\text{M}$, respectively).²

References

1. Murakami, E., Bao, H., Ramesh, M., *et al.* Mechanism of activation of β -D-2'-deoxy-2'-fluoro-2'-C-methylcytidine and inhibition of hepatitis C virus NS5B RNA polymerase. *Antimicrob. Agents Chemother.* **51**(2), 503-509 (2007).
2. Stuyver, L.J., McBrayer, T.R., Tharnish, P.M., *et al.* Inhibition of hepatitis C replicon RNA synthesis by β -D-2'-deoxy-2'-fluoro-2'-C-methylcytidine: A specific inhibitor of hepatitis C virus replication. *Antivir. Chem. Chemother.* **17**(2), 79-87 (2006).

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, 12/13/2022

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