

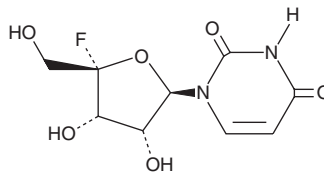
PRODUCT INFORMATION



EIDD-2749

Item No. 38249

CAS Registry No.: 1613589-24-4
Formal Name: 4'-C-fluoro-uridine
Synonyms: 4'-FIU, 4'-Fluorouridine
MF: C₉H₁₁FN₂O₆
FW: 262.2
Purity: ≥98%
UV/Vis.: λ_{max}: 261 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

EIDD-2749 is supplied as a solid. A stock solution may be made by dissolving the EIDD-2749 in the solvent of choice, which should be purged with an inert gas. EIDD-2749 is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of EIDD-2749 in these solvents is approximately 10 and 20 mg/ml, respectively.

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 EIDD-2749 can be prepared by directly dissolving the solid in aqueous buffers. The solubility of EIDD-2749 in PBS (pH 7.2) is approximately 10 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

EIDD-2749 is an antiviral ribonucleoside analog.¹ Incorporation of EIDD-2749 into RSV or severe acute respiratory syndrome coronavirus 2 (SARS-CoV-2) RNA inhibits viral RNA-dependent RNA polymerase (RdRp) and induces transcriptional stalling in cell-free assays. It reduces viral yield in HEp-2 cells infected with clinical isolates of respiratory syncytial virus (RSV; EC₅₀s = 0.61-1.2 μM). EIDD-2749 reduces apical virus shedding in RSV-infected isolated human airway epithelial (HAE) cells (EC₅₀ = 55 nM). *In vivo*, EIDD-2749 (5 mg/kg) reduces lung virus load in a mouse model of RSV infection. It also reduces nasal viral titers in a ferret model of SARS-CoV-2 infection when administered at doses of 15 and 50 mg/kg.

Reference

1. Sourimant, J., Lieber, C.M., Aggarwal, M., *et al.* 4'-Fluorouridine is an oral antiviral that blocks respiratory syncytial virus and SARS-CoV-2 replication. *Science* **375**(6577), 161-167 (2022).

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, 05/05/2023

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