

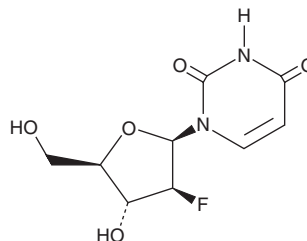
PRODUCT INFORMATION



2'-Fluoro-2'-deoxy-arabino-Uridine

Item No. 40303

CAS Registry No.: 69123-94-0
Formal Name: 1-(2-deoxy-2-fluoro-β-D-arabinofuranosyl)-2,4(1H,3H)-pyrimidinedione
Synonym: 2'-Arafluoro-2'-deoxyuridine
MF: C₉H₁₁FN₂O₅
FW: 246.2
Purity: ≥98%
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

2-Fluoro-2'-deoxy-arabino-uridine is supplied as a solid. A stock solution may be made by dissolving the fluoro-2'-deoxy-arabino-uridine in the solvent of choice, which should be purged with an inert gas. Fluoro-2'-deoxy-arabino-uridine is soluble in methanol.

Description

2'-Fluoro-2'-deoxy-arabino-uridine is an antiviral nucleoside and a derivative of the pyrimidine nucleoside uridine (Item No. 20300).¹ It inhibits the replication of herpes simplex virus 1 (HSV-1) in infected Vero cells when used at a concentration of 1 µg/ml and does not induce cytotoxicity in L5178Y or P815 cells (IC₅₀ = >10 µg/ml for both).

Reference

1. Watanabe, K.A., Reichman, U., Hirota, K., et al. Nucleosides. 110. Synthesis and antiherpes virus activity of some 2'-fluoro-2'-deoxyarabinofuranosylpyrimidine nucleosides. J. Med. Chem. 22(1), 21-24 (1979).

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, 03/19/2024

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