

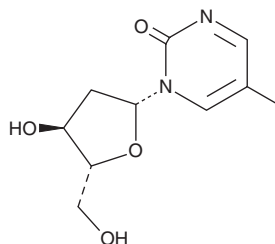
PRODUCT INFORMATION



Ropidoxuridine

Item No. 41220

CAS Registry No.: 93265-81-7
Formal Name: 1-(2-deoxy-β-D-erythro-pentofuranosyl)-5-iodo-2(1H)-pyrimidinone
Synonyms: IPdR, 5-Iodo-2-pyrimidinone-2'-deoxyribose
MF: C₉H₁₁IN₂O₄
FW: 338.1
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

Ropidoxuridine is supplied as a solid. A stock solution may be made by dissolving the ropidoxuridine in the solvent of choice, which should be purged with an inert gas. Ropidoxuridine is sparingly soluble (1-10 mg/ml) in DMSO.

Description

Ropidoxuridine is a radiosensitizer and a prodrug form of the antiviral compound idoxuridine (Item No. 20222).¹ It also inhibits viral plaque formation in HeLa cells infected by herpes simplex virus 1 (HSV-1) or HSV-2 when used at a concentration of 100 μM.² Ropidoxuridine (1,000 mg/kg per day for six days) enhances radiation-induced decreases in tumor growth without reducing body weight in an HT-29 colon cancer mouse xenograft model.¹ It increases animal survival in a mouse model of lethal HSV-2 infection when administered at a dose of 100 mg/kg every five days.²

References

1. Kinsella, T.J., Kunugi, K.A., Vielhuber, K.A., *et al.* Preclinical evaluation of 5-iodo-2-pyrimidinone-2'-deoxyribose as a prodrug for 5-iodo-2'-deoxyuridine-mediated radiosensitization in mouse and human tissues. *Clin. Cancer Res.* **4**(1), 99-109 (1998).
2. Lewandowski, G.A., Grill, S.P., Fisher, M.H., *et al.* Anti-herpes simplex virus activity of 5-substituted 2-pyrimidinone nucleosides. *Antimicrob. Agents Chemother.* **33**(3), 340-344 (1989).

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, 08/09/2024

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