

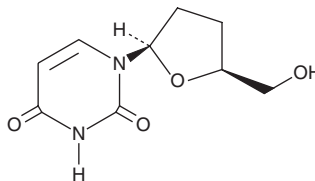
PRODUCT INFORMATION



2',3'-Dideoxyuridine

Item No. 36494

CAS Registry No.: 5983-09-5
Formal Name: 2',3'-dideoxy-uridine
Synonym: ddU
MF: $C_9H_{12}N_2O_4$
FW: 212.2
Purity: $\geq 98\%$
UV/Vis.: λ_{max} : 262 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

2',3'-Dideoxyuridine is supplied as a solid. A stock solution may be made by dissolving the 2',3'-dideoxyuridine in the solvent of choice, which should be purged with an inert gas. 2',3'-Dideoxyuridine is soluble in organic solvents such as ethanol, DMSO and dimethyl formamide (DMF). The solubility of 2',3'-dideoxyuridine in DMSO and DMF is approximately 10 mg/ml. 2',3'-Dideoxyuridine is slightly soluble in ethanol.

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 2',3'-dideoxyuridine can be prepared by directly dissolving the solid in aqueous buffers. The solubility of 2',3'-dideoxyuridine in PBS (pH 7.2) is approximately 5 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

2',3'-Dideoxyuridine is an inactive nucleoside analog.^{1,2}

References

1. Mehellou, Y., Balzarini, J., and McGuigan, C. An investigation into the anti-HIV activity of 2',3'-didehydro-2',3'-dideoxyuridine (d4U) and 2',3'-dideoxyuridine (ddU) phosphoramidate 'ProTide' derivatives. *Org. Biomol. Chem.* **7**(12), 2548-2553 (2009).
2. Pertenbreiter, F., Balzarini, J., and Meier, C. Nucleoside mono- and diphosphate prodrugs of 2',3'-dideoxyuridine and 2',3'-dideoxy-2',3'-didehydrouridine. *ChemMedChem* **10**(1), 94-106 (2015).

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, 12/19/2022

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