

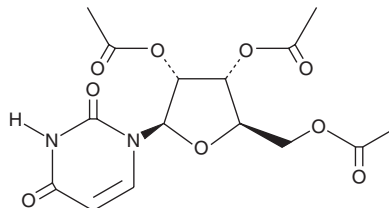
PRODUCT INFORMATION



2',3',5'-Triacetyluridine

Item No. 27445

CAS Registry No.: 4105-38-8
Formal Name: 2',3',5'-triacetate uridine
Synonyms: PN401, RG2133,
2',3',5'-Tri-O-acetyluridine,
Uridine Triacetate
MF: $C_{15}H_{18}N_2O_9$
FW: 370.3
Purity: $\geq 98\%$
UV/Vis.: λ_{max} : 257 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',5'-Triacetyluridine is supplied as a solid. A stock solution may be made by dissolving the 2',3',5'-triacetyluridine in the solvent of choice, which should be purged with an inert gas. 2',3',5'-Triacetyluridine is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of 2',3',5'-triacetyluridine in these solvents is approximately 30 mg/ml.

2',3',5'-Triacetyluridine is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, 2',3',5'-triacetyluridine should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. 2',3',5'-Triacetyluridine has a solubility of approximately 0.25 mg/ml in a 1:3 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

2',3',5'-Triacetyluridine is a prodrug of uridine (Item No. 20300).¹ It is more lipid soluble than uridine and resistant to degradation by uridine phosphorylase. It is cleaved by plasma esterases *in vivo* to release uridine. 2',3',5'-Triacetyluridine (6% in the diet) decreases neurodegeneration in the piriform cortex and striatum, as well as reduces the amount of huntingtin-positive aggregates and increases BDNF protein levels in the piriform cortex in a transgenic mouse model of Huntington's disease.² It also improves rotarod performance and increases survival in transgenic mouse models of Huntington's disease. 2',3',5'-Triacetyluridine reverses toxicity and increases survival in a mouse model of dihydropyrimidine dehydrogenase (DPD) deficiency-induced 5-fluorouracil (5-FU; Item No. 14416) overdose when used at a concentration of 2,000 mg/kg three times per day beginning within 24 hours of 5-FU administration.³ Formulations containing 2',3',5'-triacetyluridine have been used in the treatment of hereditary orotic aciduria and of overdose or life-threatening toxicity due to fluorouracil or capecitabine.

References

1. Ashour, O.M., Naguib, F.N.M., and el Kouni, M.H. 5-(*m*-Benzyloxybenzyl)barbituric acid acyclonucleoside, a uridine phosphorylase inhibitor, and 2',3',5'-tri-O-acetyluridine, a prodrug of uridine, as modulators of plasma uridine concentration: Implications for chemotherapy. *Biochem. Pharmacol.* **51(12)**, 1601-1611 (1996).
2. Saydoff, J.A., Garcia, R.A.G., Browne, S.E., *et al.* Oral uridine pro-drug PN401 is neuroprotective in the R6/2 and N171-82Q mouse models of Huntington's disease. *Neurobiol. Dis.* **24(3)**, 455-465 (2006).
3. von Borstel, R.W., O'Neil, J.D., Saydoff, J.A., *et al.* Uridine triacetate for lethal 5-FU toxicity due to dihydropyrimidine dehydrogenase (DPD) deficiency. *J. Clin. Oncol.* **28(15_suppl)**, e13505-e13505 (2010).

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, 11/04/2022

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