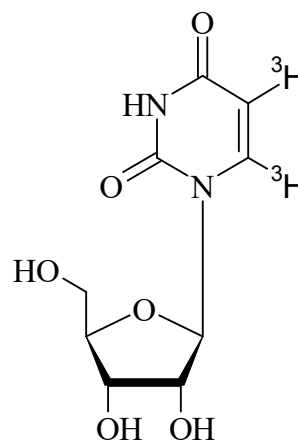


Uridine, [5,6- ^3H]-

Product Number: NET367

LOT SPECIFIC INFORMATION

Lot Number:	3386538
Specific Activity:	39.1 Ci/mmol
	1446.7 GBq/mmol
Production Date:	6 Mar 2025

M.W. 244
 $\text{C}_9\text{H}_{12}\text{N}_2\text{O}_6$ **PACKAGING:** 1.0 mCi/ml (37 MBq/ml) in aqueous solution, steri-packaged.**STABILITY AND STORAGE RECOMMENDATIONS:** When uridine, [5,6- ^3H]- is stored at 5°C in its original solvent and at its original concentration, the rate of decomposition is initially 4% per month from date of purification. Stability is nonlinear and not correlated to isotope half-life. Lot to lot variation may occur.It is recommended that aqueous solutions of uridine, [5,6- ^3H]- be used within one month of receipt. A 70% ethanol solution is recommended as a solvent if the compound is to be stored longer than one month, as the material has been found to be much more stable in this solvent. The rate of decomposition in 70% ethanol is initially 2% for 6 months.**SPECIFIC ACTIVITY RANGE:** >25 Ci/mmol (>925 GBq/mmol)**RADIOCHEMICAL PURITY:** This product was initially found to be greater than 97% when determined by the following method. The rate of decomposition can accelerate. It is advisable to check purity prior to use:

High pressure liquid chromatography on a C18-AQ column using the following mobile phases:

- 1% triethylamine acetate pH 4.
- water.

QUALITY CONTROL: The radiochemical purity of uridine, [5,6- ^3H]- is checked at appropriate intervals using the listed chromatography method.**SPECIAL INFORMATION:** To remove the 70% ethanol solution from stored material, an aliquot or the entire sample can be taken to dryness by directing a gentle stream of inert gas (nitrogen) over the surface of the solution or by rotary evaporation under vacuum. The temperature of the solution should not be allowed to exceed 20°C during the drying process, and the compound should not be permitted to remain in the solid state any longer than necessary.

HAZARD INFORMATION: WARNING: This product contains a chemical known to the state of California to cause cancer.

Limited Warranty: The information provided in this document is valid for the specified lot number and date of analysis. This information is for reference purposes only and does not constitute a warranty or guarantee of the product's suitability for any specific use. Revvity, Inc., its subsidiaries, and/or affiliates (collectively, "Revvity") do not assume any liability for any errors or damages arising from the use of this document or the product described herein. REVVITY EXPRESSLY DISCLAIMS ALL WARRANTIES, INCLUDING WARRANTIES OF MERCHANTABILITY OR FITNESS FOR A PARTICULAR PURPOSE, REGARDLESS OF WHETHER ORAL OR WRITTEN, EXPRESS OR IMPLIED, ALLEGEDLY ARISING FROM ANY USAGE OF ANY TRADE OR ANY COURSE OF DEALING, IN CONNECTION WITH THE USE OF INFORMATION CONTAINED HEREIN OR THE PRODUCT ITSELF.

NET367-R Rev-01

revvity