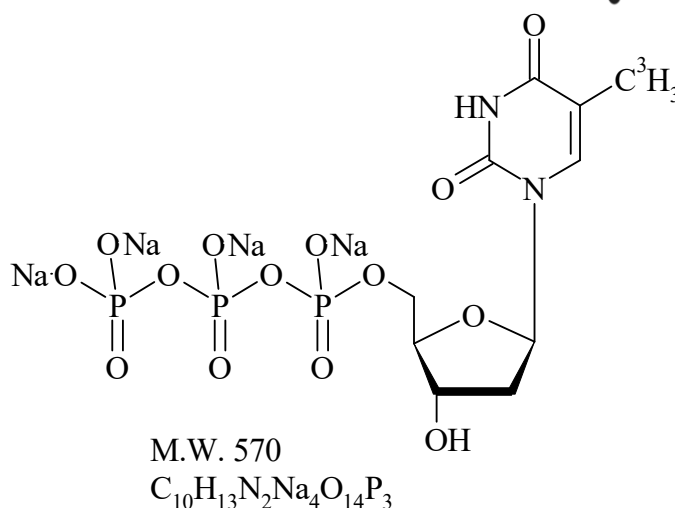


2'-Deoxythymidine-5'-Triphosphate, Tetrasodium Salt, [Methyl- ^3H]-

Product Number: NET221A

LOT SPECIFIC INFORMATION

Lot Number:	3393529
Specific Activity:	83 Ci/mmol
	3071 GBq/mmol
Production Date:	4-April-2025



PACKAGING: 2.5 mCi/ml (92.5 MBq/ml) in aqueous solution in 10mM tricine-NaOH buffer (1), pH 7.6. Shipped in dry ice.

STABILITY AND STORAGE RECOMMENDATIONS: When 2'-Deoxythymidine 5'-triphosphate, tetrasodium salt, [methyl- ^3H]- is stored at -80°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.

- At -80°C , the rate of decomposition gradually increases after 3 or 4 months.
- The compound should be stored at -80°C . Rapid freezing at dry ice temperature is recommended for thawed solutions. The purity should be checked before use. Lot to lot variation may occur.

SPECIFIC ACTIVITY RANGE: >70 Ci/mmol (>2590 GBq/mmol)

RADIOCHEMICAL PURITY: This product was initially found to be greater than 95% when determined by the following methods. The rate of decomposition can accelerate. It is advisable to check purity prior to use:

High pressure liquid chromatography on a SAX column using the following mobile phase:
0.6M ammonium phosphate, (pH 3.5).

SPECIAL INFORMATION:

1. Conversion to other salts:

The tetrasodium salt may be converted to that of another cation by passing it through a small column of exchange resin in the form of the desired cation. A 1-2 ml quantity of thoroughly washed resin is generally satisfactory. Necessary precautions should be exercised to minimize any breakdown from occurring during this procedure.

2. Removal of solvent:

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. 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. For additional information regarding stability and storage, see discussion above.

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.

NET221A-R Rev01

revvity