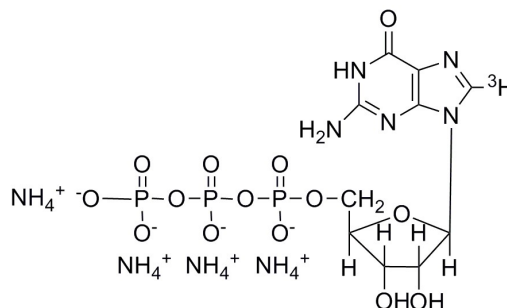


Guanosine 5'-Triphosphate, Ammonium Salt, [8-³H(N)]- (GTP, Ammonium Salt, [8-³H(N)]-)

Product Number: NET1201

LOT SPECIFIC INFORMATION

Lot Number:	3294056
Specific Activity:	3.7 Ci/mmol
	136.9 GBq/mmol
Production Date:	12-Jun-2024



M.W. 593.31
C₁₀H₁₂N₅O₁₄P₃ (4 NH₄)

PACKAGING: 1.0 mCi/ml (37 MBq/ml) in ethanol : water (1:1). Shipped in dry ice.

STABILITY AND STORAGE RECOMMENDATIONS: When guanosine 5'-triphosphate, ammonium salt, [8-³H(N)]- is stored at -20°C in its original solvent and at its original concentration, the rate of decomposition is initially 1-2% per month from date of purification. The rate of decomposition can accelerate. Stability is nonlinear and not correlated to isotope half-life. Lot to lot variation may occur.

- Ethanol is employed to reduce the rate of decomposition. If it is necessary to store aqueous solutions of this compound, the solution should be rapidly frozen by rotating the vessel in a dry-ice acetone bath.

SPECIFIC ACTIVITY RANGE: >3 Ci/mmol (>111 GBq/mmol)

RADIOCHEMICAL PURITY: This product was initially found to be greater than 97% when determined by the following methods. It is advisable to check purity prior to use:

High pressure liquid chromatography on a SAX column using the following mobile phase:
0.3M ammonium phosphate, pH 3.5.

QUALITY CONTROL: The radiochemical purity of guanosine 5'-triphosphate, ammonium salt, [8-³H(N)]- is checked at appropriate intervals using the listed chromatography method.

SPECIAL INFORMATION:

1. Conversion to other salts:

The ammonium 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 1-2ml 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: