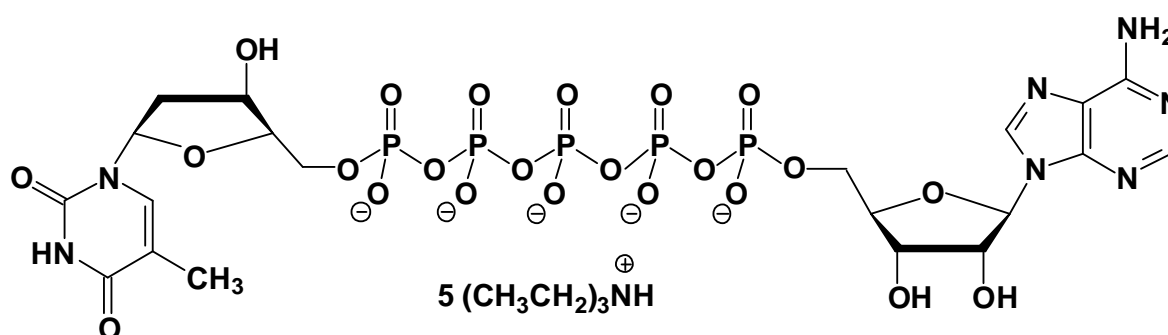


AP₅dT

P¹-(5'-adenosyl) P⁵-(5'-(2'-deoxy-thymidyl)) pentaphosphate, Triethylammonium salt

Cat. No.	Amount
NU-502S	50 Units
NU-502L	250 Units



Cat. No.: NU-502

Molecular Formula: C₂₀H₂₅N₇O₂₃P₅ (Anion)

Molecular Weight: 886.32 (Anion)

Purity: > 95%, clear aqueous solution, pH 7.5

Storage conditions:

Short term exposure (up to 1 week cumulative) to ambient temperature possible. Long term storage at < -20°C. If stored as recommended, Jena Bioscience guarantees optimal performance of this product for 12 months after date of delivery.

For research use only!

1 unit = 1 μl of a 10 mM solution

Selected References:

Lavie *et al.* Structural basis for efficient phosphorylation of 3'-azidothymidine monophosphate by Escherichia coli thymidylate kinase. *Proc. Natl. Acad. Sci. USA* **95** (24):14045.

Lavie *et al.* (1998) Crystal structure of yeast thymidylate kinase complexed with the bisubstrate inhibitor P-1-(5'-adenosyl) P-5-(5'-thymidyl) pentaphosphate (TP(5)A) at 2.0 angstrom resolution: Implications for catalysis and AZT activation. *Biochemistry* **37**:3677.