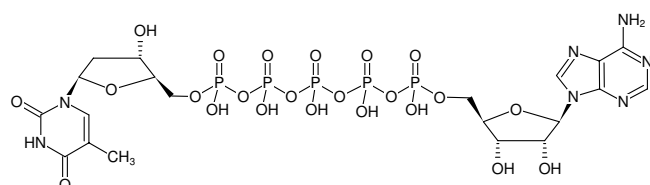


**AP₅dT**

P¹-(5'-Adenosyl) P⁵-[5'-(2'-deoxy-thymidyl)] pentaphosphate, Triethylammonium salt

| Cat. No. | Amount |
|----------|-------------------|
| NU-502S | 50 µl (10 mM) |
| NU-502L | 5 x 50 µl (10 mM) |



Structural formula of AP₅dT

For research use only!

Shipping: shipped on gel packs

Storage Conditions: store at -20 °C

Short term exposure (up to 1 week cumulative) to ambient temperature possible.

Shelf Life: 12 months after date of delivery

Molecular Formula: C₂₀H₃₀N₇O₂₃P₅ (free acid)

Molecular Weight: 891.36 g/mol (free acid)

Exact Mass: 891.01 g/mol (free acid)

CAS#: 103137-88-8

Purity: ≥ 95 % (HPLC)

Form: solution in water

Color: colorless to slightly yellow

Concentration: 10 mM - 11 mM

pH: 7.5 ± 0.5

Spectroscopic Properties: λ_{max} 262 nm, ε 22.3 L mmol⁻¹ cm⁻¹
(Tris-HCl pH 7.5)

Selected References:

Haouz *et al.* (2003) Enzymatic and Structural Analysis of Inhibitors Designed against *Mycobacterium tuberculosis* Thymidylate Kinase. *J. Biol. Chem.* **278** (7):4963.

Lavie *et al.* (1998) 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.