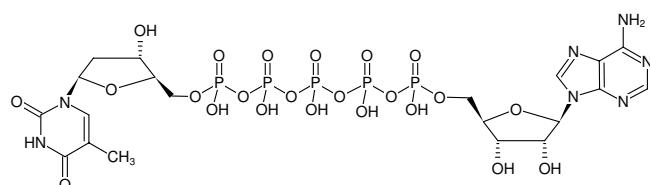


**AP<sub>5</sub>dT**

P<sup>1</sup>-(5'-Adenosyl) P<sup>5</sup>-[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 AP<sub>5</sub>dT

**For research use only!**

**Shipping:** shipped on blue ice

**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:** C<sub>20</sub>H<sub>30</sub>N<sub>7</sub>O<sub>23</sub>P<sub>5</sub> (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<sub>max</sub> 262 nm, ε 22.3 L mmol<sup>-1</sup> cm<sup>-1</sup> (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.