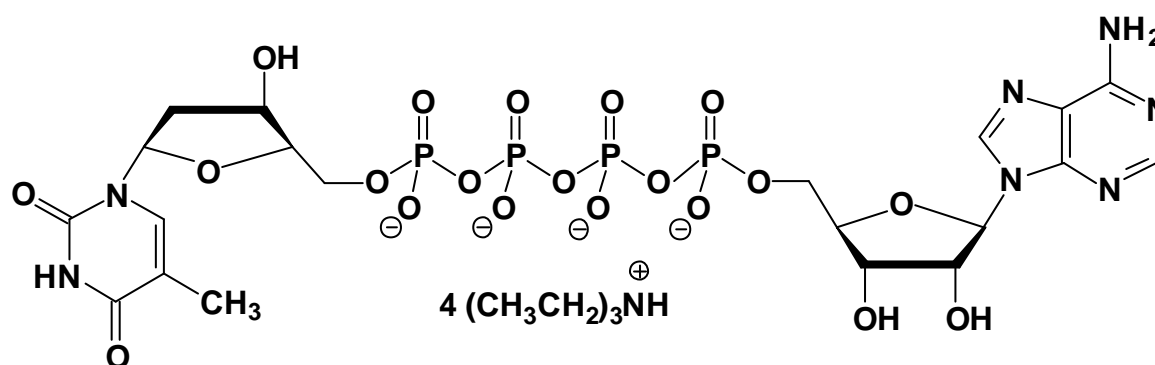


AP₄dT

P¹-(5'-adenosyl) P⁴-(5'-(2'-deoxy-thymidyl)) tetraphosphate, Triethylammonium salt

Cat. No.	Amount
NU-501S	50 Units
NU-501L	250 Units



Cat. No.: NU-501

Molecular Formula: C₂₀H₂₅N₇O₂OP₄ (Anion)

Molecular Weight: 807.34 (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