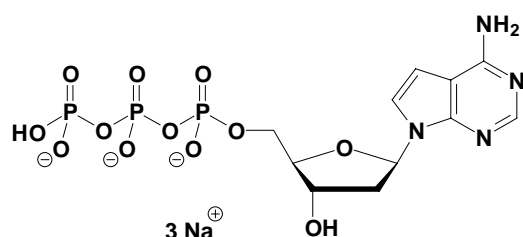


7-Deaza-dATP

7-Deaza-2'-Deoxy-Adenosine-5'-triphosphate, Sodium salt

Cat. No.	Amount
NU-1175S	100 Units
NU-1175L	500 Units



Cat. No.: NU-1175

Molecular Formula: C₁₁H₁₄N₄O₁₂P₃ (Anion)

Molecular Weight: 487.17 (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.

* 1 unit = 1 µl of a 10 mM solution

For research use only!

Selected References:

Fletcher *et al.* (1996) Human Telomerase Inhibition by 7'-Deaza-2'-deoxypurine Nucleoside Triphosphates. *Biochemistry* **35** (49):15611.

Yamakawa *et al.* (1997) A DNA cycle sequencing reaction that minimizes compression on automated fluorescent sequencers. *Nucleic Acids Research* **25** (6):1311.