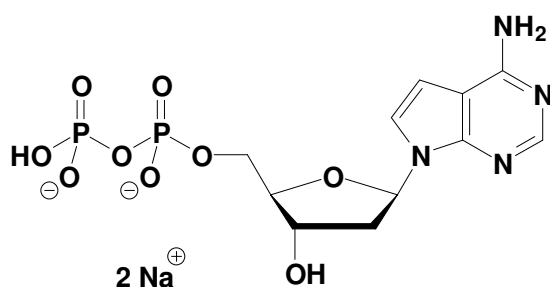


7-Deaza-dADP

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

Cat. No.	Amount
NU-1174S	100 Units
NU-1174L	500 Units



Cat. No.: NU-1174

Molecular Formula: C₁₁H₁₄N₄O₉P₂ (Anion)

Molecular Weight: 408.2 (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.