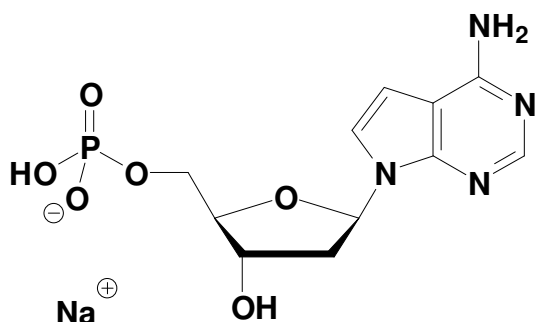


## 7-Deaza-dAMP

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

Cat. No.	Amount
NU-1176S	100 Units
NU-1176L	500 Units



### 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.

**Cat. No.:** NU-1176

**Molecular Formula:** C<sub>11</sub>H<sub>14</sub>N<sub>4</sub>O<sub>6</sub>P (Anion)

**Molecular Weight:** 329.2 (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!**