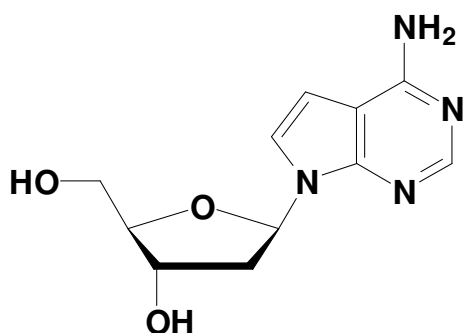


7-Deaza-2'-deoxyadenosine

Cat. No.	Amount
N-1046	100 mg



Cat. No.: N-1046

Molecular Formula: C₁₁H₁₄N₄O₃

Molecular Weight: 250.25

Purity: > 98%

Storage conditions:

Short term exposure (up to 1 week cumulative) to ambient temperature possible. Long term storage at +4°C. If stored as recommended, Jena Bioscience guarantees optimal performance of this product for 12 months after date of delivery.

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.