

7-Deaza-2'-Deoxy-Adenosine Kit

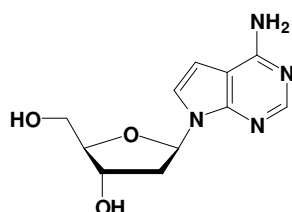
Cat.No.: NK-111

Kit Contents

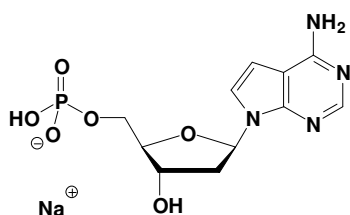
7-Deaza-2'-Deoxy-Adenosine Analogs	Cat. No.	Amount
7-Deaza-dA	N-1046	3 mg
7-Deaza-dAMP	NU-1176	100 Units
7-Deaza-dADP	NU-1174	100 Units
7-Deaza-dATP	NU-1175	100 Units

1 Unit = 1 μ l of a 10 mM solution

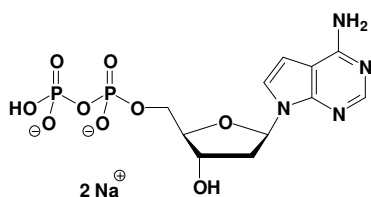
Structures



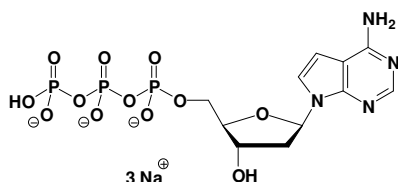
7-Deaza-dA



7-Deaza-dAMP



7-Deaza-dADP



7-Deaza-dATP

Introduction

7-Deaza-2'-Deoxy-Adenosines are able to replace their natural counterparts in PCR reactions, which leads to modified DNA with improved properties, such as lower melting temperature of the double strand. They are therefore used in DNA sequencing reactions, for PCR of GC-rich templates and to improve the DNA bands in polyacrylamide gel electrophoresis. Furthermore, they lead to increased stability of the helical DNA structure.

Kit Description

The 7-Deaza-2'-Deoxy-Adenosine Kit contains a set of 4 typical 7-Deaza-2'-Deoxy-Adenosine analogs (Nucleoside, 5'-Mono, Di and Triphosphate).

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 compressions on automated fluorescent sequencers. *Nucleic Acids Research* **25** (6):1311.

Storage and Stability

Short term exposure (up to 1 week cumulative) to ambient temperature is possible. Long term storage is recommended at < -20°C. If properly stored, Jena Bioscience guarantees optimal performance of the compounds for 12 months after date of delivery.