

# 3'-Deoxy-Guanosine Kit

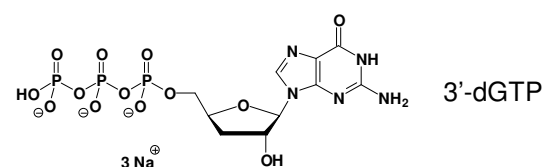
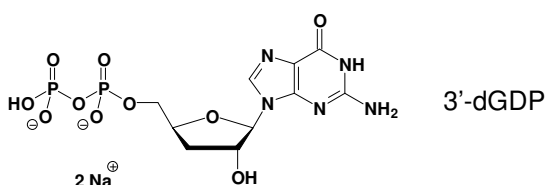
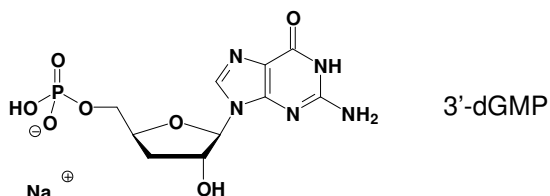
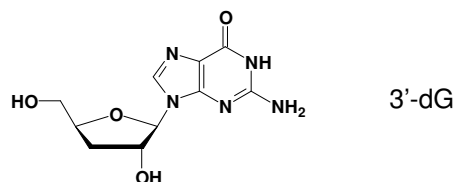
Cat.No.: NK-114

## Kit Contents

3'-Deoxy-Guanosine Analogs	Cat-No.	Amount
3'-dG	N-1048	1 mg
3'-dGMP	NU-1143	50 Units
3'-dGDP	NU-1144	50 Units
3'-dGTP	NU-1145	50 Units

1 Unit = 1  $\mu$ l of a 10 mM solution

## Structures



## Introduction

3'-Deoxy Guanosines are used in structure/activity investigations and as chain terminators in PCR reactions.

## Kit Description

The 3'-Deoxy-Guanosine Kit contains a set of 4 typical 3'-Deoxy-Guanosine analogs (Nucleoside, 5'-Mono, Di and Triphosphate).

## Selected References:

Shunji Izuta *et al.* (1996) 3'-Deoxyribonucleotides Inhibit Eukaryotic DNA Primase1. *J. Biochem*, **119** (6):1038.

Rasmussen *et al.* (1993) In vivo transcriptional pausing and cap formation on three *Drosophila* heat shock genes. *Proc. Natl. Acad. Sci. USA* **90**:7923.

Hamel *et al.* (1976) Interactions of Guanosine Triphosphate Analogues with Elongation Factor G of *Escherichia coli*. *Eur. J. Biochem.* **63**:431.

## 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.