

# 6-Methylthio-Guanosine Kit

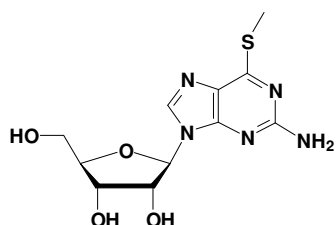
Cat.No.: NK-104

## Kit Contents

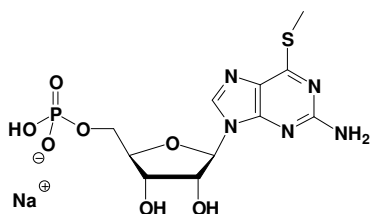
6-Methylthio-Guanosine Analogs	Cat. No.	Amount
6-Methylthio-G	N-1042	3 mg
6-Methylthio-GMP	NU-1128	100 Units
6-Methylthio-GDP	NU-1129	100 Units
6-Methylthio-GTP	NU-1130	100 Units

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

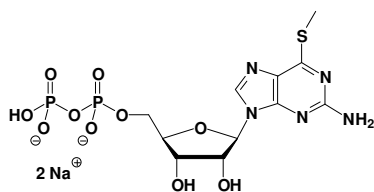
## Structures



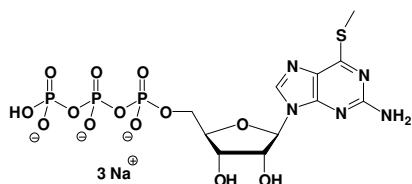
6-Methylthio-G



6-Methylthio-GMP



6-Methylthio-GDP



6-Methylthio-GTP

## Introduction

6-Methylthio-Guanosines are structurally related to 6-Thio-GTP. 6-Thioguanosines are metabolites of Azathioprine, a drug used in the therapy of chronic inflammatory and auto-immune diseases. The immunosuppressive properties of Azathioprine are very likely caused by binding of 6-Thio-GTP to the GTPase Rac.

## Kit Description

The 6-Methylthio-Guanosine Kit contains a set of 4 typical 6-Methylthio-Guanosine analogs (Nucleoside, 5'-Mono, Di and Triphosphate).

## Selected References:

Nelson *et al.* (1977) Oxypurine and 6-thiopurine nucleoside triphosphate formation in human erythrocytes. *Adv. Exp. Med. Biol.* **76A**:121.

## Storage and Stability

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