

# 6-Methylthio-Inosine Kit

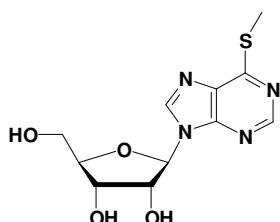
Cat.No.: NK-105

## Kit Contents

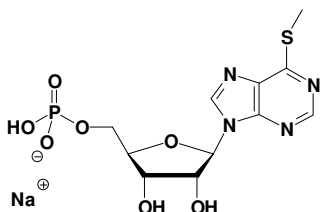
6-Methylthio-Inosine Analogs	Cat. No.	Amount
6-Methylthio-I	N-1053	3 mg
6-Methylthio-IMP	NU-1131	100 Units
6-Methylthio-IDP	NU-1132	100 Units
6-Methylthio-ITP	NU-1133	100 Units

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

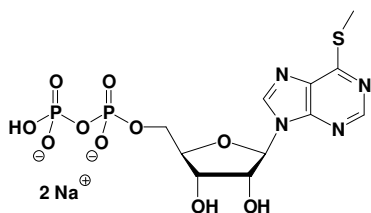
## Structures



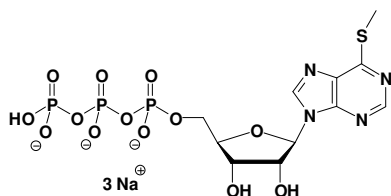
6-Methylthio-I



6-Methylthio-IMP



6-Methylthio-IDP



6-Methylthio-ITP

## Introduction

6-Methylthio-Inosines are structurally related to 6-Thio-GTP. 6-Thio-Guanosines 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-Inosine Kit contains a set of 4 typical 6-Methylthio-Inosine analogs. (Nucleoside, 5'-Mono, Di and Triphosphate).

## Selected References:

Lavi *et al.* (1985) A rapid and sensitive high-performance liquid chromatographic assay for 6-mercaptopurine metabolites in red blood cells. *Anal. Biochem.* **144** (2):514.

## 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.