

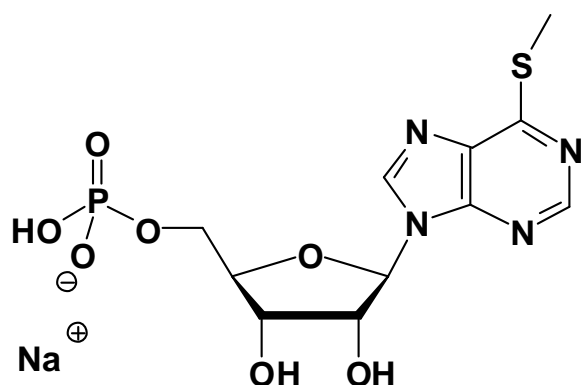
## 6-Methylthio-IMP

6-Methylthioinosine-5'-monophosphate, Sodium salt

Cat. No.	Amount
NU-1131S	150 Units
NU-1131L	750 Units

### 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.



**Cat. No.:** NU-1131

**Molecular Formula:** C<sub>11</sub>H<sub>14</sub>N<sub>4</sub>O<sub>7</sub>PS (Anion)

**Molecular Weight:** 377.28 (Anion)

**Purity:** > 95%, clear aqueous solution, pH 7.5

**Spectroscopic properties:**  $\lambda_{\max}$  224/292 nm;  
 $\epsilon$  11500/19000

### Storage conditions:

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

### For research use only!

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