

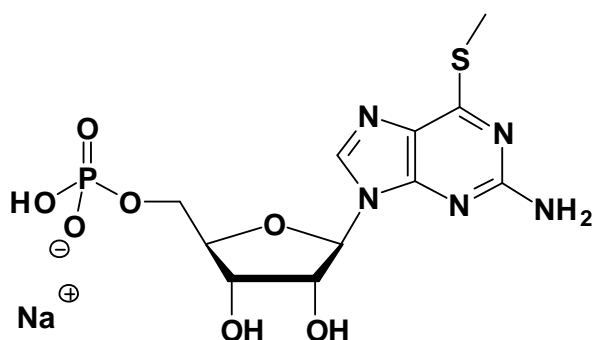
## 6-Methylthio-GMP

6-Methylthioguanosine-5'-monophosphate, Sodium salt

Cat. No.	Amount
NU-1128S	100 Units
NU-1128L	500 Units

### Selected References:

Kaminsky *et al.* (2007) The c-Myc Target Gene Rcl (C6orf108) Encodes a Novel Enzyme, Deoxynucleoside 5'-monophosphate N-Glycosidase. *J. Biol. Chem.* **282** (11): 8150.



**Cat. No.:** NU-1128

**Molecular Formula:** C<sub>11</sub>H<sub>15</sub>N<sub>5</sub>O<sub>7</sub>PS (Anion)

**Molecular Weight:** 392.30 (Anion)

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

**Spectroscopic properties:**  $\lambda_{\max}$  242 nm;  
 $\epsilon$  12800

### 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