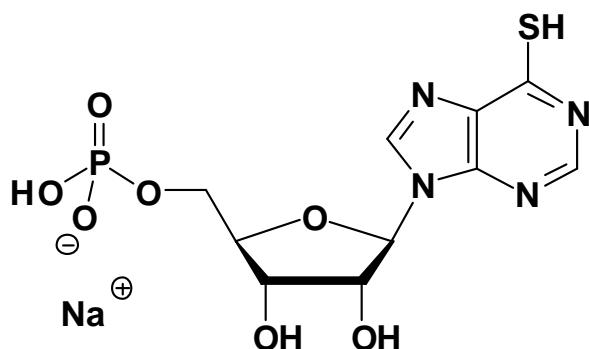


# 6-Mercaptopurine-riboside-5'-monophosphate

6-Mercaptopurine-riboside-5'-monophosphate, Sodium salt

Cat. No.	Amount
NU-1148S	100 Units
NU-1148L	500 Units



**Cat. No.:** NU-1148

**Molecular Formula:** C<sub>10</sub>H<sub>12</sub>N<sub>4</sub>O<sub>7</sub>PS (Anion)

**Molecular Weight:** 363.26 (Anion)

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

**Spectroscopic properties:** λ<sub>max</sub> 322 nm;  
ε 27300

### 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 μl of a 10 mM solution

### Selected References:

Wielinga *et al.* (2002) Thiopurine metabolism and identification of the thiopurine metabolites transported by MRP4 and MRP5 overexpressed in human embryonic kidney cells. *Mol Pharmacol.* **62(6)**:1321.

Szabados *et al.* (1994) 5-Aminoimidazole-4-carboxamide ribotide transformylase-IMP cyclohydrolase from human CCRFCM leukemia cells: purification, pH dependence, and inhibitors. *Biochemistry.* **33(47)**:14237.

Stet *et al.* (1993) A biochemical basis for synergism of 6-mercaptopurine and mycophenolic acid in Molt F4, a human malignant T-lymphoblastic cell line. *Biochim Biophys Acta.* **1180(3)**:277.

Drake *et al.* (1982) Metabolism and mechanisms of action of 9-(tetrahydro-2-furyl)-6-mercaptopurine in Chinese hamster ovary cells. *Chem Biol Interact.* **41(1)**:105.