

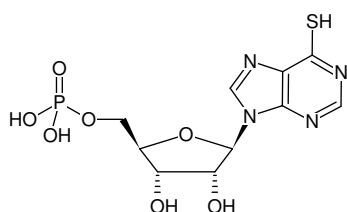


## 6-Mercaptopurine-riboside-5'-monophosphate

6-Thio-Inosine-5'-monophosphate

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

Cat. No.	Amount
NU-1148S	100 µl (10 mM)
NU-1148L	5 x 100 µl (10 mM)



Structural formula of 6-Mercaptopurine-riboside-5'-monophosphate

**For research use only!**

**Shipping:** shipped on gel packs

**Storage Conditions:** store at -20 °C

Short term exposure (up to 1 week cumulative) to ambient temperature possible.

**Shelf Life:** 12 months after date of delivery

**Molecular Formula:** C<sub>10</sub>H<sub>13</sub>N<sub>4</sub>O<sub>7</sub>PS (free acid)

**Molecular Weight:** 364.26 g/mol (free acid)

**Exact Mass:** 364.02 g/mol (free acid)

**CAS#:** 53-83-8

**Purity:** ≥ 95 % (HPLC)

**Form:** solution in water

**Color:** colorless to slightly yellow

**Concentration:** 10 mM - 11 mM

**pH:** 7.5 ± 0.5

**Spectroscopic Properties:** λ<sub>max</sub> 322 nm, ε 27.3 L mmol<sup>-1</sup> cm<sup>-1</sup> (Tris-HCl pH 7.5)

### Selected References:

Neurath *et al.* (2005) 6-Thioguanosine Diphosphate and Triphosphate Levels in Red Blood Cells and Response to Azathioprine Therapy in Crohn's Disease. *Clinical Gastroenterology and Hepatology* **3** (10): 1007.

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.