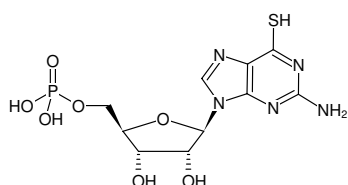


**6-Thio-GMP**

6-Thio-guanosine-5'-monophosphate, Sodium salt

Cat. No.	Amount
NU-1121S	150 µl (10 mM)
NU-1121L	5 x 150 µl (10 mM)



Structural formula of 6-Thio-GMP

**For general laboratory use.****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>14</sub>N<sub>5</sub>O<sub>7</sub>PS (free acid)**Molecular Weight:** 379.28 g/mol (free acid)**Exact Mass:** 379.04 g/mol (free acid)**CAS#:** 15867-02-4**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> 257 nm, ε 8.8 L mmol<sup>-1</sup> cm<sup>-1</sup> (Tris-HCl pH 7.5), λ<sub>max</sub> 342 nm, ε 24.8 L mmol<sup>-1</sup> cm<sup>-1</sup> (Tris-HCl pH 7.5)**Selected References:**

Karner *et al.* (2010) Determination of 6-thioguanosine diphosphate and triphosphate and nucleoside diphosphate kinase activity in erythrocytes: novel targets for thiopurine therapy? *Ther. Drug Monit.* **32** (2):119.

Courtois *et al.* (1969) Interaction entre pigments et acides nucleiques Etude sur l'interaction de l'Actinomycine D avec les nucléosides par Dichroïsme Circulaire. *FEBS Lett.* **3**:330.