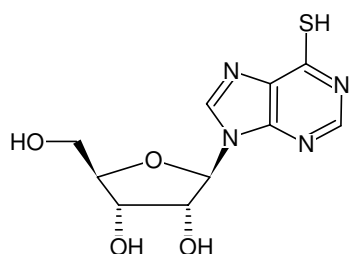


**6-Mercaptopurine-riboside**

6-Thio-Inosine

Cat. No.	Amount
N-1063-100	100 mg



Structural formula of 6-Mercaptopurine-riboside

For general laboratory use.**Shipping:** shipped at ambient temperature**Storage Conditions:** store at 4 °C

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

Shelf Life: 24 months after date of delivery**Molecular Formula:** C₁₀H₁₂N₄O₄S**Molecular Weight:** 284.29 g/mol**Exact Mass:** 284.06 g/mol**CAS#:** 4988-64-1, 574-25-4**EC number:** 209-371-5**Purity:** ≥ 95 % (HPLC)**Form:** solid**Color:** white to off-white