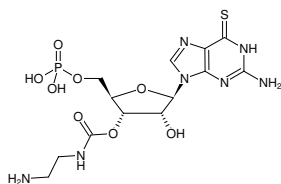
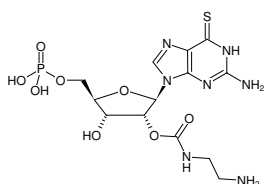


**EDA-6-Thio-GMP**

2'/3'-O-(2-Aminoethyl-carbamoyl)-6-Thio-guanosine-5'-monophosphate, Sodium salt

Cat. No.	Amount
NU-1614S	50 µl (10 mM)
NU-1614L	5 x 50 µl (10 mM)



Structural formula of EDA-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:** C₁₃H₂₀N₇O₈PS (free acid)**Molecular Weight:** 465.38 g/mol (free acid)**Exact Mass:** 465.08 g/mol (free acid)**Purity:** ≥ 95 % (HPLC)**Form:** solution in water**Concentration:** 10 mM - 11 mM**pH:** 7.5 ±0.5**Spectroscopic Properties:** λ_{max} 257/342 nm, ε 8.8/24.8 L mmol⁻¹ cm⁻¹
(Tris-HCl pH 7.5)