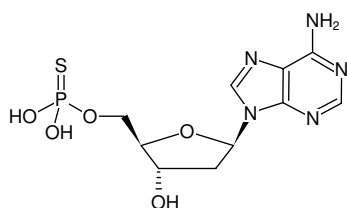


**dAMP α S**2'-Deoxyadenosine-5'-(α -thio)-monophosphate, Sodium salt

Cat. No.	Amount
NU-430-1	1 mg
NU-430-5	5 mg

Structural formula of dAMP α S**For research use only!****Shipping:** shipped on blue ice**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:** C₁₀H₁₄N₅O₅PS (free acid)**Molecular Weight:** 347.29 g/mol (free acid)**Exact Mass:** 347.05 g/mol (free acid)**Purity:** \geq 97 % (HPLC)**Form:** solid**Spectroscopic Properties:** λ_{\max} 259 nm, ϵ 15.4 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)