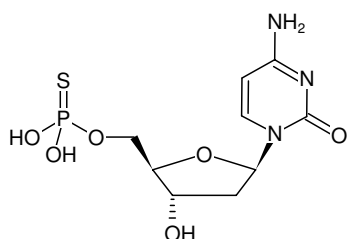


**dCMPαS**

2'-Deoxycytidine-5'-(α-thio)-monophosphate, Sodium salt
2'-Deoxycytidine-5'-O-monophosphorothioate

Cat. No.	Amount
NU-447-1	1 mg
NU-447-5	5 mg



Structural formula of dCMPαS

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: C₉H₁₄N₃O₆PS (free acid)

Molecular Weight: 323.26 g/mol (free acid)

Exact Mass: 323.03 g/mol (free acid)

CAS#: 63225-09-2 (component: 64145-27-3)

Purity: ≥ 97 % (HPLC)

Form: white powder

Spectroscopic Properties: λ_{max} 271 nm, ε 8.9 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)