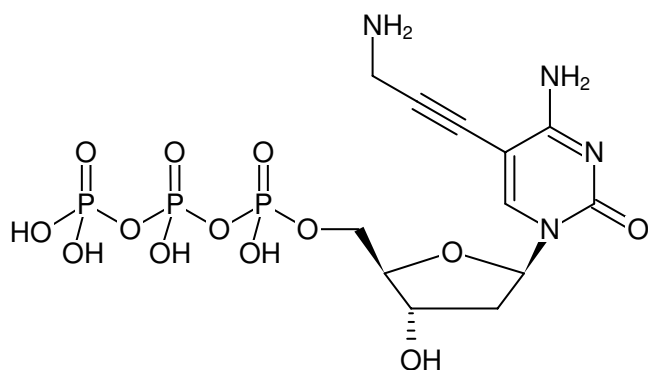




5-Propargylamino-dCTP - Solid

5-Propargylamino-2'-deoxycytidine-5'-triphosphate, Sodium salt

Cat. No.	Amount
NU-809-1	1 mg
NU-809-5	5 mg



Structural formula of 5-Propargylamino-dCTP - Solid

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: C₁₂H₁₉N₄O₁₃P₃ (free acid)

Molecular Weight: 520.22 g/mol (free acid)

Exact Mass: 520.02 g/mol (free acid)

CAS#: 115899-39-3

Purity: ≥ 93 % (HPLC)

Form: solid

Color: white to off-white

Spectroscopic Properties: λ_{max} 294 nm, ε 9.3 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)