

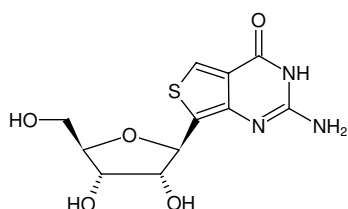


Thienoguanosine

Thieno-G, thG

2-Aminothienu[3,4-d]pyrimidine-ribonucleoside

Cat. No.	Amount
NU-317	10 μ l (100 mM)



Structural formula of Thienoguanosine

For general laboratory use.

Please centrifuge briefly before opening (volume \leq 2 ml).

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: C₁₁H₁₃N₃O₅S (free acid)

Molecular Weight: 299.30 g/mol (free acid)

Exact Mass: 299.06 g/mol (free acid)

Purity: \geq 95 % (HPLC)

Form: solution in DMSO

Color: colorless

Concentration: 100 mM - 110 mM

Spectroscopic Properties: λ_{exc} 321 nm, λ_{em} 453 nm, ϵ 4.15 L mmol⁻¹ cm⁻¹ (water)