



5-AcOHg-dCTP

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

Cat. No.	Amount
NU-1149S	100 µl (10 mM)
NU-1149L	5 x 100 µl (10 mM)

Applications:

Selective isolation of Hg-DNA^[1]

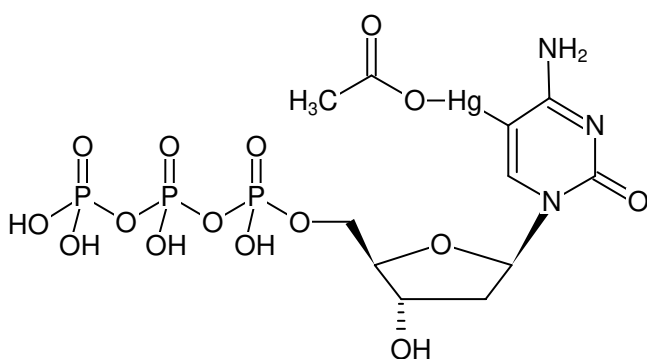
Substrate for polymerases^[2]

Crystallization with heavy atom^[2]

Selected References:

[1] Banfalvi *et al.* (1995) Effect of mercury substitution of DNA on its susceptibility to cleavage by restriction endonucleases. *DNA Cell Biol.* **14** (5):445.

[2] Dale *et al.* (1973) The Synthesis and Enzymatic Polymerization of Nucleotides Containing Mercury: Potential Tools for Nucleic Acid Sequencing and Structural Analysis. *Proc. Natl. Acad. Sci. USA* **70** (8):2238.



Structural formula of 5-AcOHg-dCTP

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: C₁₁H₁₈HgN₃O₁₅P₃ (free acid)

Molecular Weight: 725.78 g/mol (free acid)

Exact Mass: 726.97 g/mol (free acid)

Purity: ≥ 95 % (HPLC)

Form: solution in water

Color: colorless to slightly yellow

Concentration: 10 mM - 11 mM

pH: 7.5 ± 0.5

Spectroscopic Properties: λ_{max} 272 nm, ε 9.1 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)