

**5-AcOHg-dUTP**

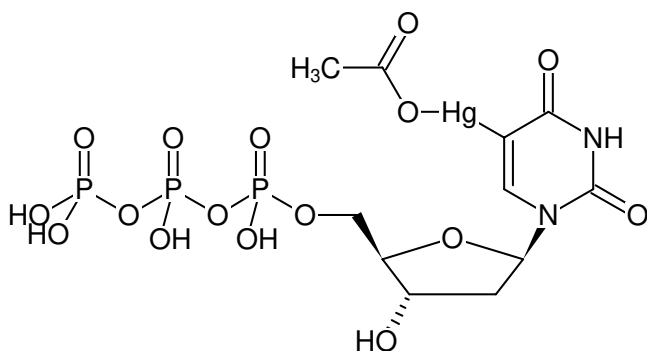
5-Mercurylacetate-2'-deoxyuridine-5'-triphosphate, Triethylammonium salt

Cat. No.	Amount
NU-910S	100 µl (10 mM)
NU-910L	5 x 100 µl (10 mM)

**Selected References:**

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

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-dUTP

**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:** C<sub>11</sub>H<sub>17</sub>HgN<sub>2</sub>O<sub>16</sub>P<sub>3</sub> (free acid)**Molecular Weight:** 726.77 g/mol (free acid)**Exact Mass:** 727.95 g/mol (free acid)**Purity:** ≥ 95 % (HPLC)**Form:** clear aqueous solution**Concentration:** 10 mM - 11 mM**pH:** 7.5 ±0.5**Spectroscopic Properties:** λ<sub>max</sub> 267 nm, ε 10.1 L mmol<sup>-1</sup> cm<sup>-1</sup> (Tris-HCl pH 7.5)