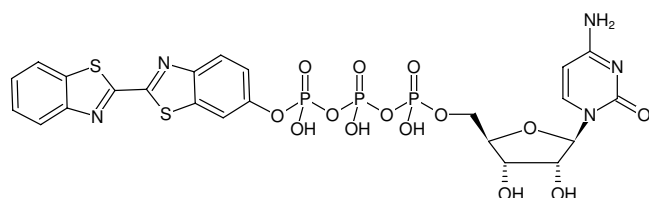


**γ-(BBT)-CTP**

ATTO-CTP

γ-[2'(2-Benzothiazoyl)-6'-hydroxybenzothiazole]-cytidine-5'-triphosphate, Sodium salt

Cat. No.	Amount
NU-933	1 mg



Structural formula of γ-(BBT)-CTP

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:** C₂₃H₂₂N₅O₁₄P₃S₂ (free acid)**Molecular Weight:** 749.50 g/mol (free acid)**Exact Mass:** 748.89 g/mol (free acid)**Purity:** ≥ 95 % (HPLC)**Form:** solid**Color:** white to off-white**Spectroscopic Properties:** λ_{max} 355 nm, ε 26.5 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)**Selected References:**

Tay *et al.* (2016) The C-terminal 18 Amino Acid Region of Dengue Virus NS5 Regulates its Subcellular Localization and Contains a Conserved Arginine Residue Essential for Infectious Virus Production. *PLoS Pathog.* **12** (9):e1005886.

Niyomrattanakit *et al.* (2015) Stabilization of dengue virus polymerase in de novo initiation assay provides advantages for compound screening. *Antiviral Res.* **119**:36.