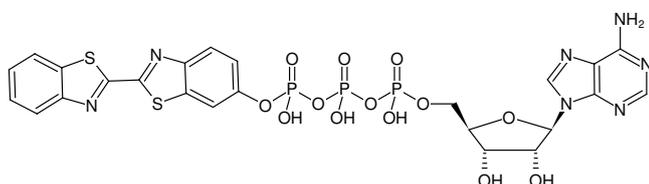


**γ-(BBT)-ATP**

ATTO-ATP

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

Cat. No.	Amount
NU-1700	1 mg



Structural formula of γ-(BBT)-ATP

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:** C₂₄H₂₂N₇O₁₃P₃S₂ (free acid)**Molecular Weight:** 773.51 g/mol (free acid)**Exact Mass:** 772.99 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:**Niyomrattanakit *et al.* (2015) Stabilization of dengue virus polymerase in de novo initiation assay provides advantages for compound screening. *Antiviral Res.* **119**:36.Takahashi *et al.* (2012) Establishment of a robust dengue virus NS3-NS5 binding assay for identification of protein-protein interaction inhibitors. *Antiviral Res.* **96 (3)**:305.Niyomrattanakit *et al.* (2011) A Fluorescence-Based Alkaline Phosphatase-Coupled Polymerase Assay for Identification of Inhibitors of Dengue Virus RNA-Dependent RNA Polymerase. *J. Biomol. Screen.* **16 (2)**:201.