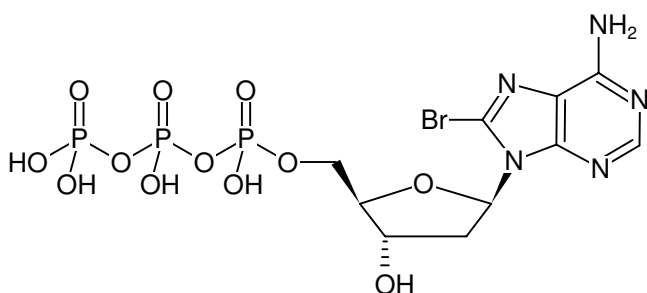


**8-Bromo-dATP**

(8Br-dATP)

8-Bromo-2'-deoxyadenosine-5'-triphosphate, Sodium salt

Cat. No.	Amount
NU-117S	50 µl (10 mM)
NU-117L	5 x 50 µl (10 mM)



Structural formula of 8-Bromo-dATP

For research use only!**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:** C₁₀H₁₅N₅O₁₂P₃Br (free acid)**Molecular Weight:** 570.08 g/mol (free acid)**Exact Mass:** 568.91 g/mol (free acid)**CAS#:** 92410-99-6**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} 264 nm, ε 17.0 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)