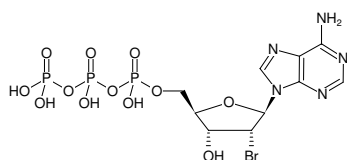


**2'-Bromo-dATP**

(2'Br-dATP)

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

Cat. No.	Amount
NU-102	150 µl (10 mM)



Structural formula of 2'-Bromo-dATP

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:** C₁₀H₁₅N₅O₁₂P₃Br (free acid)**Molecular Weight:** 570.08 g/mol (free acid)**Exact Mass:** 568.91 g/mol (free acid)**CAS#:** 107245-23-8**Purity:** ≥ 95 % (HPLC)**Form:** clear aqueous solution**Concentration:** 10 mM - 11 mM**pH:** 7.5 ±0.5**Spectroscopic Properties:** λ_{max} 259 nm, ε 15.1 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.0)**Selected References:**Gruen *et al.* (1999) 2'-Halo-ATP and -GTP analogues: Rational phasing tools for protein crystallography. *Protein Sci.* **8**:2524.Naber *et al.* (1995) A novel adenosine-triphosphate analog with a heavy-atom to target the nucleotide-binding site of proteins. *Protein Sci.* **4**:1824.