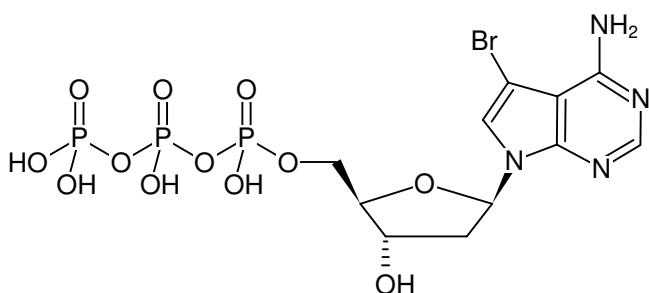


**7-Deaza-7-bromo-dATP**

(7-Deaza-7-Br-dATP)

7-Deaza-7-bromo-2'-deoxyadenosine-5'-triphosphate, Sodium salt

Cat. No.	Amount
NU-149S	50 µl (10 mM)
NU-149L	5 x 50 µl (10 mM)



Structural formula of 7-Deaza-7-bromo-dATP

**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:** C<sub>11</sub>H<sub>16</sub>N<sub>4</sub>O<sub>12</sub>P<sub>3</sub>Br (free acid)**Molecular Weight:** 569.09 g/mol (free acid)**Exact Mass:** 567.92 g/mol (free acid)**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<sub>max</sub> 277 nm, ε 9.5 L mmol<sup>-1</sup> cm<sup>-1</sup> (Tris-HCl pH 7.5)