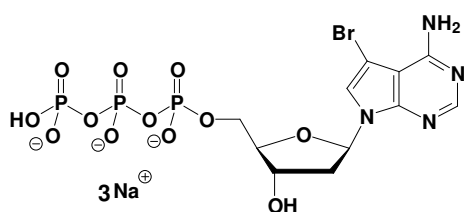


7-Deaza-7-Br-dATP

7-Deaza-7-Bromo-2'-Deoxy-Adenosine-5'-triphosphate, Sodium salt

Cat. No.	Amount
NU-149S	50 Units
NU-149L	250 Units



Cat. No.: NU-149

Molecular Formula: $C_{11}H_{13}N_4O_{12}P_3Br$ (Anion)

Molecular Weight: 566.06 (Anion)

Purity: > 95%, clear aqueous solution, pH 7.5

Storage conditions:

Short term exposure (up to 1 week cumulative) to ambient temperature possible. Long term storage at < -20°C. If stored as recommended, Jena Bioscience guarantees optimal performance of this product for 12 months after date of delivery.

* 1 unit = 1 μ l of a 10 mM solution

For research use only!