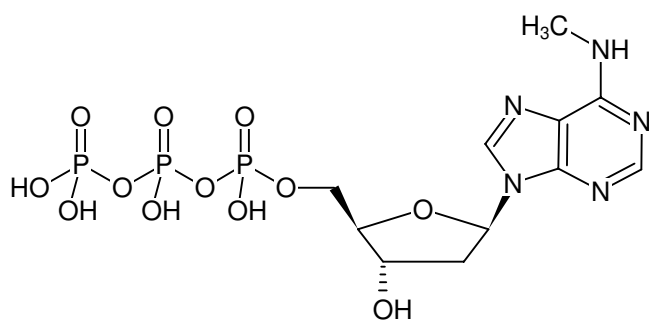


**N<sup>6</sup>-Methyl-dATP**

N<sup>6</sup>-Methyl-2'-deoxyadenosine-5'-triphosphate, Sodium salt  
N<sup>6</sup>-Methyl-2'-dATP

Cat. No.	Amount
NU-949S	10 µl (100 mM)
NU-949L	5 x 10 µl (100 mM)



Structural formula of N<sup>6</sup>-Methyl-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:** C<sub>11</sub>H<sub>18</sub>N<sub>5</sub>O<sub>12</sub>P<sub>3</sub> (free acid)

**Molecular Weight:** 505.21 g/mol (free acid)

**Exact Mass:** 505.02 g/mol (free acid)

**Purity:** ≥ 95 % (HPLC)

**Form:** solution in water

**Color:** colorless to slightly yellow

**Concentration:** 100 mM - 110 mM

**pH:** 7.5 ± 0.5

**Spectroscopic Properties:** λ<sub>max</sub> 269 nm, ε 15.4 L mmol<sup>-1</sup> cm<sup>-1</sup>  
(Tris-HCl pH 7.5)

**Selected References:**

Choi *et al.* (2016) The use of modified and non-natural nucleotides provide unique insights into pro-mutagenic replication catalyzed by polymerase eta. *Nucleic Acids Res.* **44** (3):1022.

Devadoss *et al.* (2007) Enhancing the 'A-rule' of translesion DNA synthesis: promutagenic DNA synthesis using modified nucleoside triphosphates. *Biochemistry* **46** (48):13752.

Ratel *et al.* (2006) N6-methyladenine: the other methylated base of DNA. *Bioessays*. **28** (3):309.

Kuwahara *et al.* (2003) Simultaneous incorporation of three different modified nucleotides during PCR. *Nucleic Acids Res. Suppl.* **3**:37.