

**3'-O-Azidomethyl-dATP**

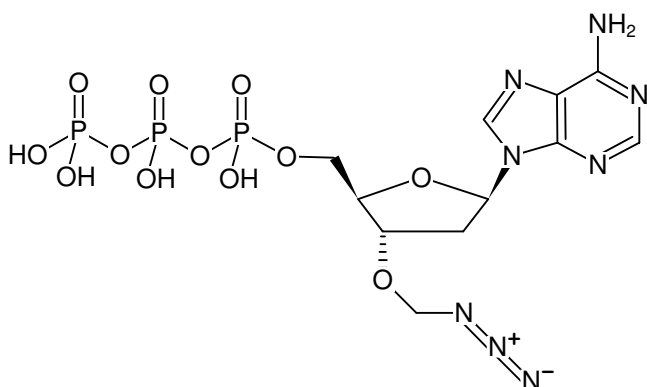
3'-O-Azidomethyl-2'-deoxyadenosine-5'-triphosphate, Sodium salt

Cat. No.	Amount
NU-937-1	1 mg

**Selected References:**

Palla *et al.* (2014) DNA sequencing by synthesis using 3'-O-azidomethyl nucleotide reversible terminators and surface-enhanced Raman spectroscopic detection. *RSC Adv.* **4** (90):49342.

Zavgorodny *et al.* (1991) Alkylthioalkylation of Nucleoside Hydroxyl Functions and Its Synthetic Applications: A New Versatile Method in Nucleoside Chemistry. *Tetrahedron Letters* **32** (51):7593.



Structural formula of 3'-O-Azidomethyl-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:** C<sub>11</sub>H<sub>17</sub>N<sub>8</sub>O<sub>12</sub>P<sub>3</sub> (free acid)**Molecular Weight:** 546.22 g/mol (free acid)**Exact Mass:** 546.02 g/mol (free acid)**Purity:** ≥ 95 % (HPLC)**Form:** solid**Color:** white to off-white**Spectroscopic Properties:** λ<sub>max</sub> 259 nm, ε 15.4 L mmol<sup>-1</sup> cm<sup>-1</sup> (Tris-HCl pH 7.5)