



3'-O-Azidomethyl-dTTP

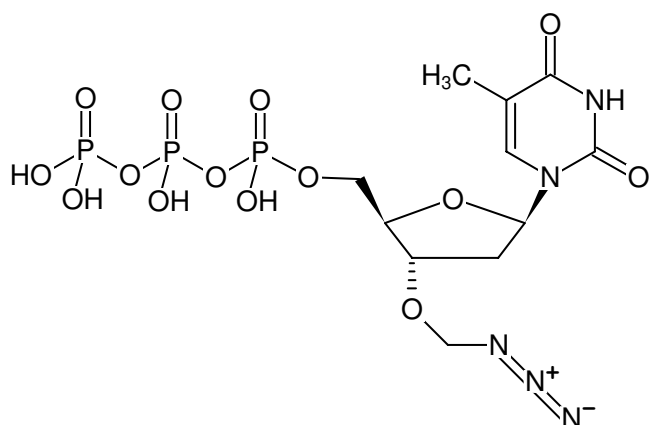
3'-O-Azidomethyl-2'-deoxythymidine-5'-triphosphate, Sodium salt

Cat. No.	Amount
NU-939-1	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-dTTP

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: C₁₁H₁₈N₅O₁₄P₃ (free acid)

Molecular Weight: 537.21 g/mol (free acid)

Exact Mass: 537.01 g/mol (free acid)

Purity: ≥ 95 % (HPLC)

Form: solid

Color: white to off-white

Spectroscopic Properties: λ_{max} 267 nm, ε 9.6 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)