



### 3'-O-Azidomethyl-dTTP

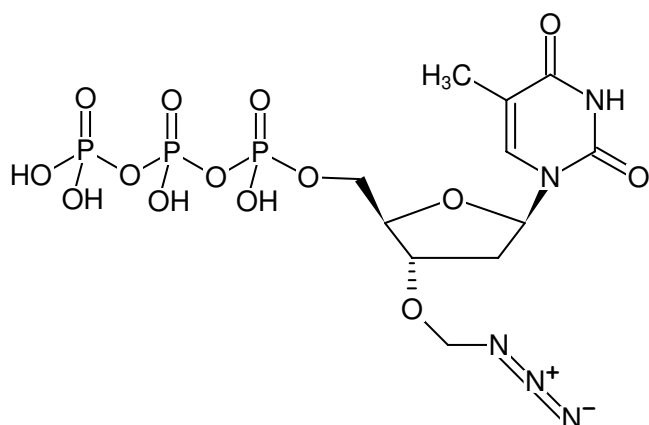
3'-O-Azidomethyl-2'-deoxythymidine-5'-triphosphate, Sodium salt

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:** C<sub>11</sub>H<sub>18</sub>N<sub>5</sub>O<sub>14</sub>P<sub>3</sub> (free acid)

**Molecular Weight:** 537.21 g/mol (free acid)

**Exact Mass:** 537.01 g/mol (free acid)

**Purity:** ≥ 95 % (HPLC)

**Form:** solid

**Spectroscopic Properties:** λ<sub>max</sub> 267 nm, ε 9.6 L mmol<sup>-1</sup> cm<sup>-1</sup> (Tris-HCl pH 7.5)