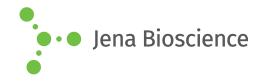
# **DATA SHEET**





## ■ 3'-O-Azidomethyl-dGTP

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

Cat. No.	Amount
NU-938-1	1 mg

# HO OH OH OH OH NH2

Structural formula of 3'-O-Azidomethyl-dGTP

## For general laboratory use.

**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>17</sub>N<sub>8</sub>O<sub>13</sub>P<sub>3</sub> (free acid)

Molecular Weight: 562.22 g/mol (free acid)

Exact Mass: 562.01 g/mol (free acid)

**Purity:** ≥ 95 % (HPLC)

Form: solid

Color: white to off-white

Spectroscopic Properties: λ<sub>max</sub> 252 nm, ε 13.7 L mmol<sup>-1</sup> cm<sup>-1</sup> (Tris-HCl

pH 7.5)

### 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.