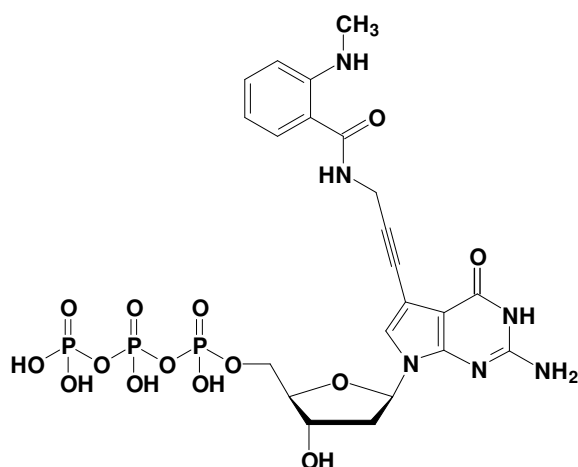


## 7-Propargylamino-7-deaza-dGTP - MANT

7-Deaza-7-propargylamino-2'-deoxy-guanosine-5'-triphosphate, labeled with MANT, Triethylammonium salt

Cat. No.	Amount
NU-1615-MNT	100 $\mu$ l (1 mM)



structural formula of 7-Propargylamino-7-deaza-dGTP - MANT

**Cat. No.:** NU-1615-MNT

**Molecular Formula:**  $C_{22}H_{27}N_6O_{14}P_3$  (free acid)

**Molecular Weight:** 692.40 (free acid)

**Purity:** > 95%, clear aqueous solution, pH 7.5

**Spectroscopic properties:**

$\lambda_{exc}$  335 nm;  $\lambda_{em}$  440 nm;  $\epsilon$  2,700  $cm^{-1} M^{-1}$

**Storage conditions:**

Short term exposure (up to 1 week cumulative) to ambient temperature possible. Long term storage at < -20°C. If stored as recommended, Jena Bioscience guarantees optimal performance of this product for 12 months after date of delivery.

**For research use only!**