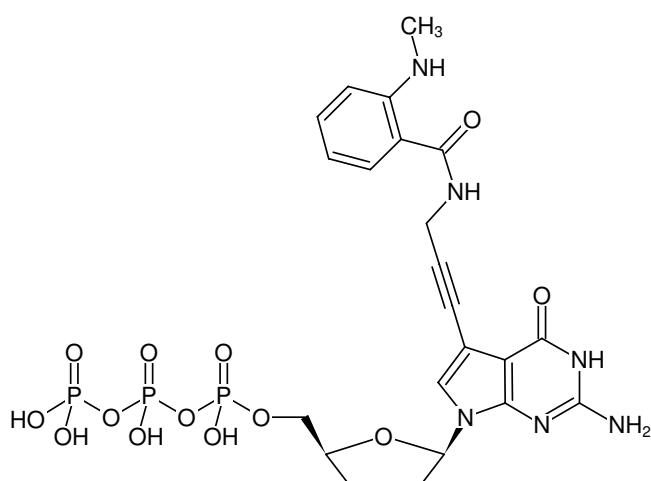




## 7-Propargylamino-7-deaza-ddGTP-MANT

7-Deaza-7-propargylamino-2',3'-dideoxyguanosine-5'-triphosphate, labeled with MANT, Triethylammonium salt

Cat. No.	Amount
NU-1618-MNT	20 $\mu$ l (5 mM)



Structural formula of 7-Propargylamino-7-deaza-ddGTP-MANT

### 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>22</sub>H<sub>27</sub>N<sub>6</sub>O<sub>13</sub>P<sub>3</sub> (free acid)

**Molecular Weight:** 676.41 g/mol (free acid)

**Exact Mass:** 676.08 g/mol (free acid)

**Purity:**  $\geq$  95 % (HPLC)

**Form:** solution in water

**Color:** colorless to slightly yellow

**Concentration:** 5.0 mM - 5.5 mM

**pH:** 7.5  $\pm$  0.5

**Spectroscopic Properties:**  $\lambda_{exc}$  335 nm,  $\lambda_{em}$  440 nm,  $\epsilon$  2.7 L mmol<sup>-1</sup> cm<sup>-1</sup> (Tris-HCl pH 7.5)