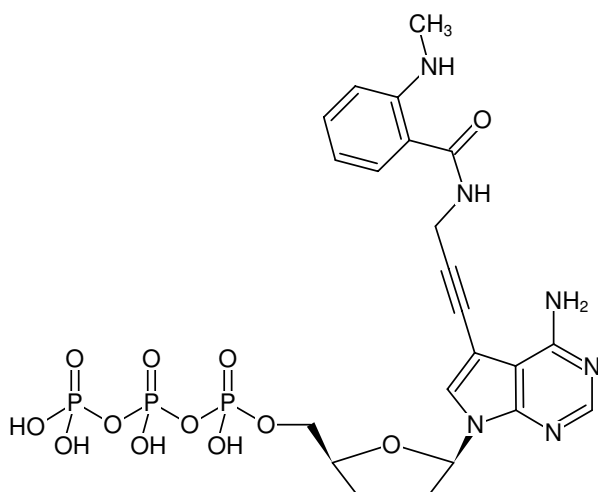




### 7-Propargylamino-7-deaza-ddATP-MANT

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

Cat. No.	Amount
NU-1612-MNT	25 µl (5 mM)



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

**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>22</sub>H<sub>27</sub>N<sub>6</sub>O<sub>12</sub>P<sub>3</sub> (free acid)

**Molecular Weight:** 660.41 g/mol (free acid)

**Purity:** ≥ 95 % (HPLC)

**Form:** solution in water

**Color:** colorless to slightly yellow

**Concentration:** 5.0 mM - 5.5 mM

**pH:** 7.5 ± 0.5

**Spectroscopic Properties:** λ<sub>exc</sub> 335 nm, λ<sub>em</sub> 440 nm, ε 2.7 L mmol<sup>-1</sup> cm<sup>-1</sup> (Tris-HCl pH 7.5)