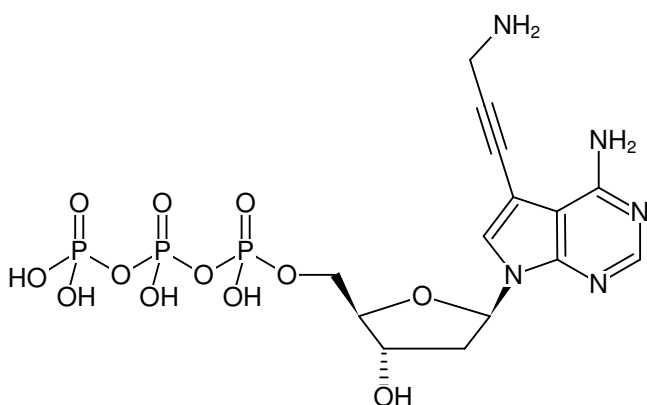


**7-Deaza-7-propargylamino-dATP**

7-Deaza-7-propargylamino-2'-deoxyadenosine-5'-triphosphate, Triethylammonium salt

Cat. No.	Amount
NU-1611-1	1 mg (1,84 µmol)



Structural formula of 7-Deaza-7-propargylamino-dATP

**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>14</sub>H<sub>20</sub>N<sub>5</sub>O<sub>12</sub>P<sub>3</sub> (free acid)**Molecular Weight:** 543.26 g/mol (free acid)**Exact Mass:** 543.03 g/mol (free acid)**Purity:** ≥ 95 % (HPLC)**Form:** solid**Spectroscopic Properties:** λ<sub>abs</sub> 280 nm, ε 12.7 L mmol<sup>-1</sup> cm<sup>-1</sup> (Tris-HCl pH 7.5)