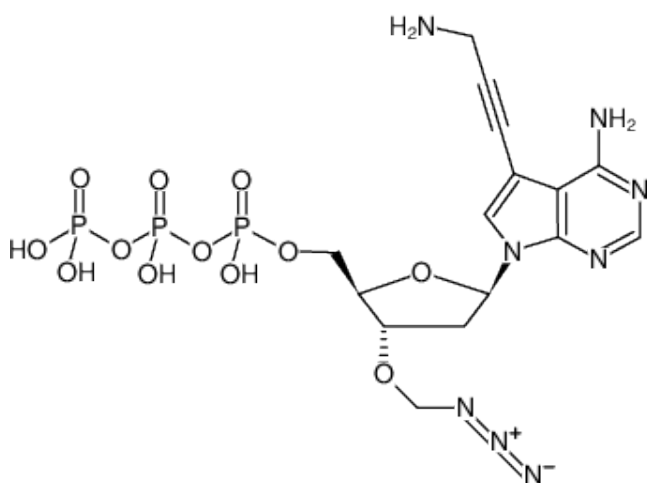




### 3'-O-Azidomethyl-7-Deaza-7-Propargylamino-dATP

3'-O-Azidomethyl-7-Deaza-7-Propargylamino-2'-deoxyadenosine-5'-triphosphate, Sodium salt  
7-Deaza-7-Propargylamino-3'-Azidomethyl-dATP

Cat. No.	Amount
NU-287-1	1 mg



Structural formula of 3'-O-Azidomethyl-7-Deaza-7-Propargylamino-dATP

#### 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>15</sub>H<sub>21</sub>N<sub>8</sub>O<sub>12</sub>P<sub>3</sub> (free acid)

**Molecular Weight:** 598.30 g/mol (free acid)

**Exact Mass:** 598.05 g/mol (free acid)

**Purity:** ≥ 95 % (HPLC)

**Form:** solid

**Color:** white to off-white

**Spectroscopic Properties:** λ<sub>max</sub> 280 nm, ε 12.7 L mmol<sup>-1</sup> cm<sup>-1</sup> (Tris-HCl pH 7.5)