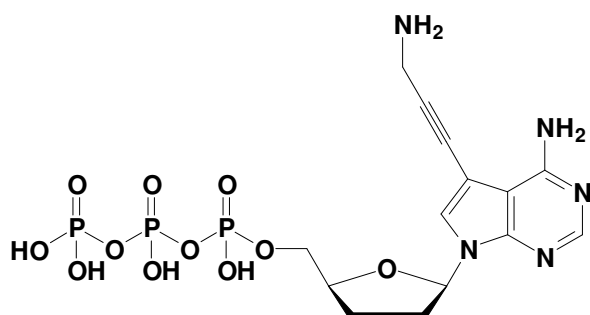


## 7-Deaza-7-propargylamino-ddATP

7-Deaza-7-propargylamino-2'/3'-dideoxy-adenosine-5'-triphosphate, sodium salt

Cat. No.	Amount
NU-1612-1	1 mg



**Cat. No.:** NU-1612

**Molecular Formula:** C<sub>14</sub>H<sub>20</sub>N<sub>5</sub>O<sub>11</sub>P<sub>3</sub> (free acid)

**Molecular Weight:** 527.26 (free acid)

**Purity:** > 95%

### 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!**