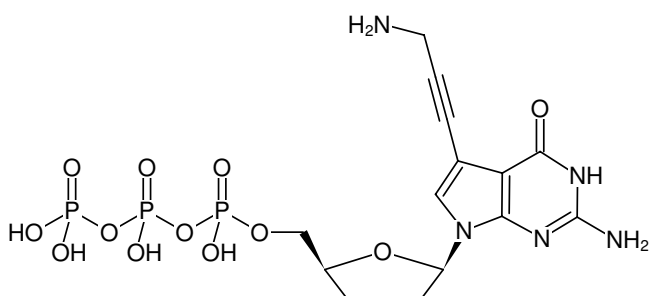




## 7-Deaza-7-propargylamino-ddGTP

7-Deaza-7-propargylamino-2',3'-dideoxyguanosine-5'-triphosphate, Lithium salt

Cat. No.	Amount
NU-994	1 mg



Structural formula of 7-Deaza-7-propargylamino-ddGTP

**For research use only!**

**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>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:** ≥ 93 % (HPLC)

**Form:** solid

**Color:** white to off-white

**Spectroscopic Properties:** λ<sub>max</sub> 272 nm, ε 11.9 L mmol<sup>-1</sup> cm<sup>-1</sup> (Tris-HCl pH 7.5)