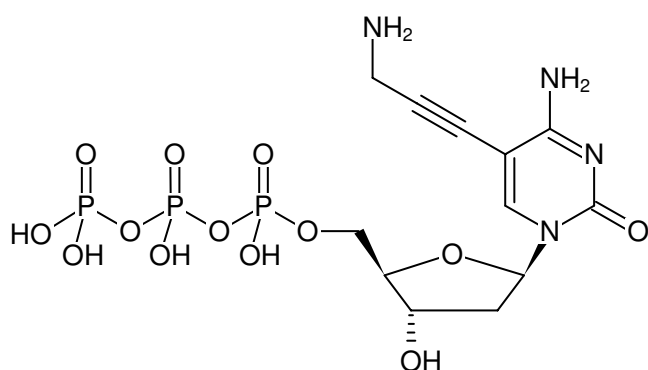




### 5-Propargylamino-dCTP - Solid

5-Propargylamino-2'-deoxycytidine-5'-triphosphate, Sodium salt

Cat. No.	Amount
NU-809-1	1 mg
NU-809-5	5 mg



Structural formula of 5-Propargylamino-dCTP - Solid

**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>12</sub>H<sub>19</sub>N<sub>4</sub>O<sub>13</sub>P<sub>3</sub> (free acid)

**Molecular Weight:** 520.22 g/mol (free acid)

**Exact Mass:** 520.02 g/mol (free acid)

**CAS#:** 115899-39-3

**Purity:** ≥ 93 % (HPLC)

**Form:** solid

**Spectroscopic Properties:** λ<sub>max</sub> 294 nm, ε 9.3 L mmol<sup>-1</sup> cm<sup>-1</sup> (Tris-HCl pH 7.5)