



### 7-Propargylamino-7-deaza-ddGTP-ATTO-647N

7-Deaza-7-propargylamino-2',3'-dideoxyguanosine-5'-triphosphate, labeled with ATTO 647N, Triethylammonium salt

Cat. No.	Amount
NU-1618-647N	80 µl (0,5 mM)

**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>56</sub>H<sub>69</sub>N<sub>8</sub>O<sub>14</sub>P<sub>3</sub> (free acid)

**Molecular Weight:** 1171.11 g/mol (free acid)

**Exact Mass:** 1170.41 g/mol (free acid)

**Purity:** ≥ 95 % (HPLC)

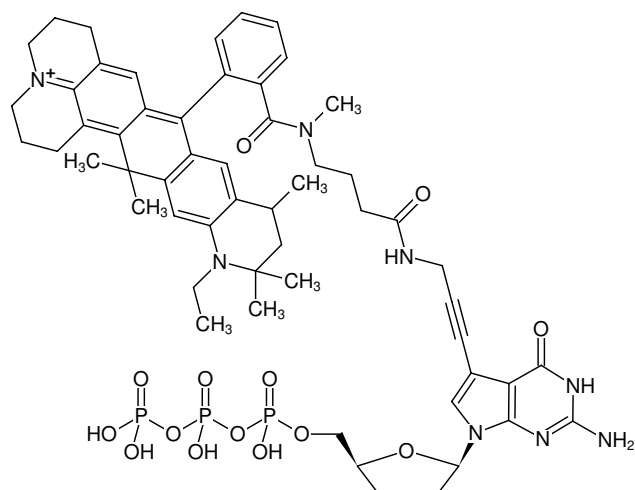
**Form:** solution in water

**Color:** blue

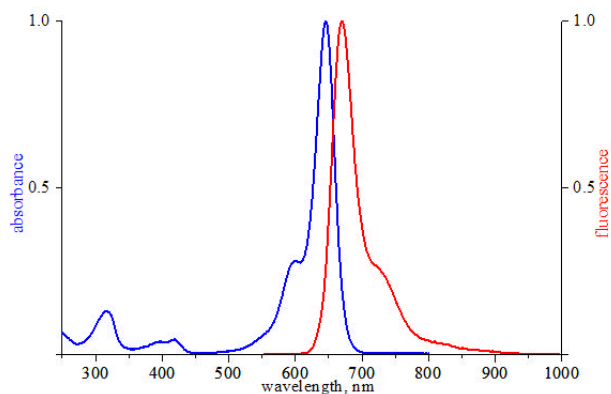
**Concentration:** 0.50 mM - 0.55 mM

**pH:** 7.5 ± 0.5

**Spectroscopic Properties:** λ<sub>exc</sub> 646 nm, λ<sub>em</sub> 664 nm, ε 150.0 L mmol<sup>-1</sup> cm<sup>-1</sup> (Tris-HCl pH 7.5)



Structural formula of 7-Propargylamino-7-deaza-ddGTP-ATTO-647N



excitation and emission spectrum of ATTO 647N