



### 7-Propargylamino-7-deaza-dGTP - ATTO-680

7-Deaza-7-propargylamino-2'-deoxyguanosine-5'-triphosphate, labeled with ATTO 680, Triethylammonium salt

Cat. No.	Amount
NU-1615-680	40 µl (1 mM)

**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>14</sub>H<sub>20</sub>N<sub>5</sub>O<sub>13</sub>P<sub>3</sub> - ATTO 680 (free acid)

**Molecular Weight:** 1066.25 g/mol (free acid)

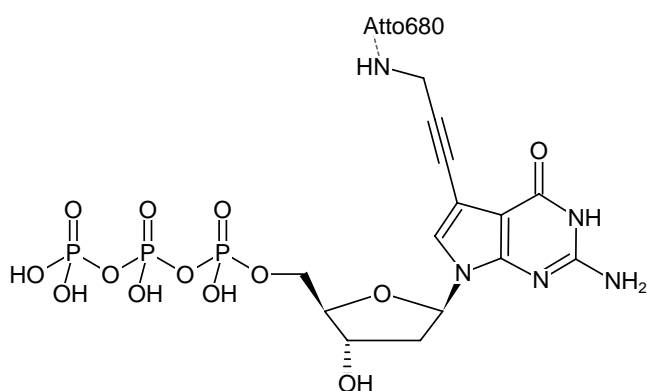
**Purity:** ≥ 95 % (HPLC)

**Form:** clear aqueous solution

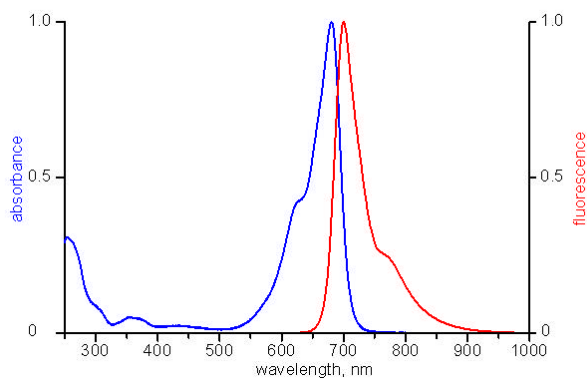
**Concentration:** 1.0 mM - 1.1 mM

**pH:** 7.5 ± 0.5

**Spectroscopic Properties:** λ<sub>exc</sub> 681 nm, λ<sub>em</sub> 698 nm, ε 125.0 L mmol<sup>-1</sup> cm<sup>-1</sup> (Tris-HCl pH 7.5)



Structural formula of 7-Propargylamino-7-deaza-dGTP - ATTO-680



excitation and emission spectrum of ATTO 680