



7-Propargylamino-7-deaza-dGTP-ATTO-594

7-Deaza-7-propargylamino-2'-deoxyguanosine-5'-triphosphate, labeled with ATTO 594, Triethylammonium salt

| Cat. No. | Amount |
|-------------|--------------|
| NU-1615-594 | 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: C₅₅H₆₅N₈O₂₂P₃S₂ (free acid)

Molecular Weight: 1347.20 g/mol (free acid)

Exact Mass: 1346.29 g/mol (free acid)

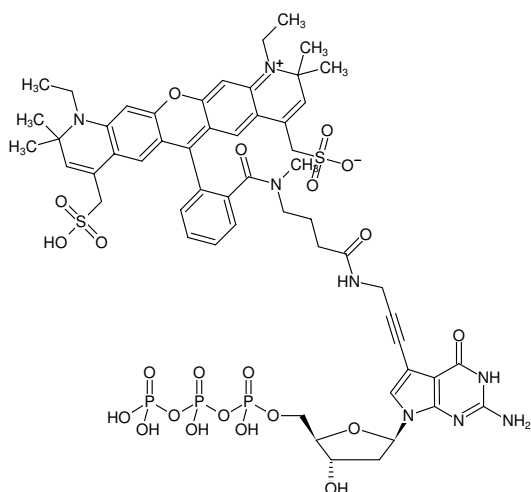
Purity: ≥ 95 % (HPLC)

Form: clear aqueous solution

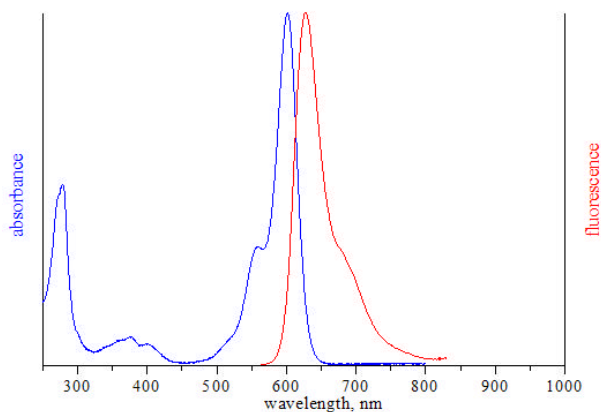
Concentration: 1.0 mM - 1.1 mM

pH: 7.5 ± 0.5

Spectroscopic Properties: λ_{exc} 603 nm, λ_{em} 626 nm, ε 120.0 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)



Structural formula of 7-Propargylamino-7-deaza-dGTP-ATTO-594



excitation and emission spectrum of ATTO 594