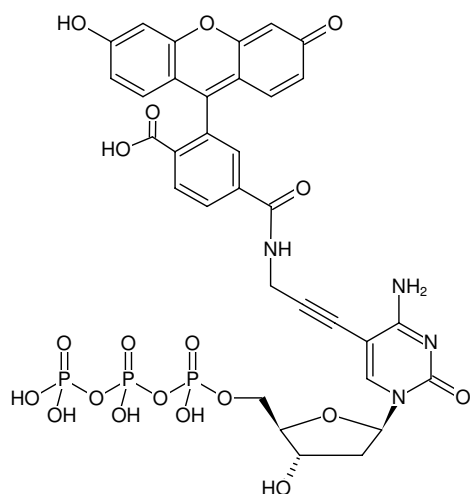




### 5-Propargylamino-dCTP-6-FAM

5-Propargylamino-2'-deoxycytidine-5'-triphosphate, labeled with 6 FAM, Triethylammonium salt

Cat. No.	Amount
NU-809-6FM	30 µl (1 mM)



Structural formula of 5-Propargylamino-dCTP-6-FAM

**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>33</sub>H<sub>29</sub>N<sub>4</sub>O<sub>19</sub>P<sub>3</sub> (free acid)

**Molecular Weight:** 878.53 g/mol (free acid)

**Exact Mass:** 878.06 g/mol (free acid)

**Purity:** ≥ 95 % (HPLC)

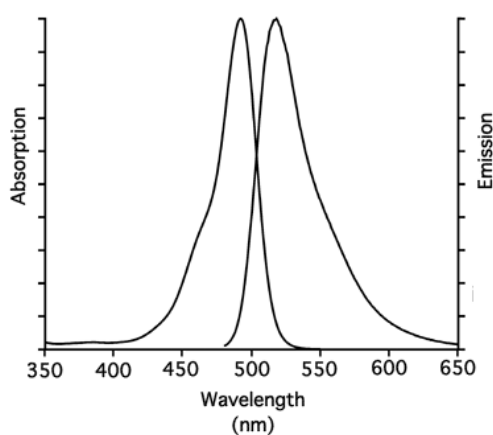
**Form:** solution in water

**Color:** yellow

**Concentration:** 1.0 mM - 1.1 mM

**pH:** 7.5 ± 0.5

**Spectroscopic Properties:** λ<sub>exc</sub> 492 nm, λ<sub>em</sub> 517 nm, ε 83.0 L mmol<sup>-1</sup> cm<sup>-1</sup> (Tris-HCl pH 7.5)



excitation and emission spectrum of 6-FAM