

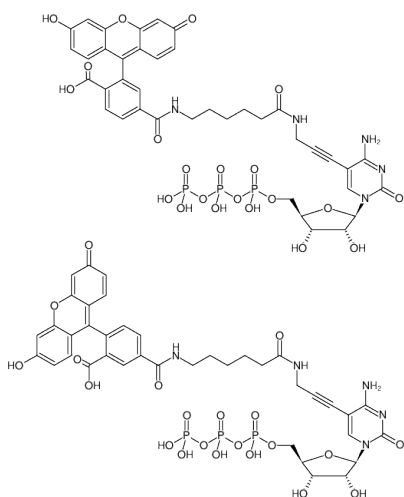


## Fluorescein-12-CTP

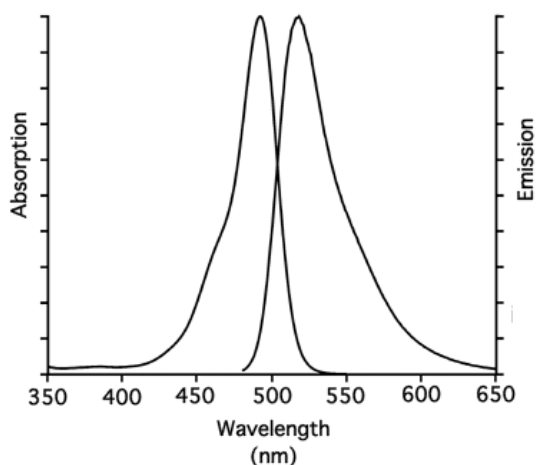
Fluorescein-X-5-Propargylamino-CTP

5/6-Fluorescein-X-5-Propargylamino-cytidine-5'-triphosphate, Triethylammonium salt

Cat. No.	Amount
NU-831-FAMX-S	10 µl (5 mM)
NU-831-FAMX-L	5 x 10 µl (5 mM)



Structural formula of Fluorescein-12-CTP



excitation and emission spectrum of 5/6-FAM

**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>39</sub>H<sub>40</sub>N<sub>5</sub>O<sub>21</sub>P<sub>3</sub> (free acid)

**Molecular Weight:** 1007.68 g/mol (free acid)

**Exact Mass:** 1007.14 g/mol (free acid)

**Purity:** ≥ 95 % (HPLC)

**Form:** solution in 10 mM Tris-HCl

**Color:** yellow-orange

**Concentration:** 5.0 mM - 5.5 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)