



## 5-Propargylamino-ddCTP-ATTO-MB2

5-Propargylamino-2',3'-dideoxycytidine-5'-triphosphate, labeled with ATTO-MB2, Triethylammonium salt

Cat. No.	Amount
NU-850-MB2	40 µl (1 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>31</sub>H<sub>38</sub>N<sub>7</sub>O<sub>13</sub>P<sub>3</sub>S<sub>1</sub> (free acid)

**Molecular Weight:** 841.66 g/mol (free acid)

**Exact Mass:** 841.15 g/mol (free acid)

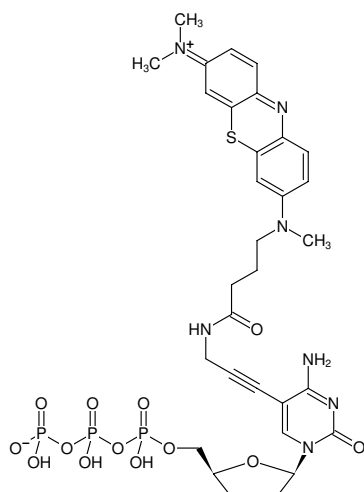
**Purity:** ≥ 95 % (HPLC)

**Form:** solution in water

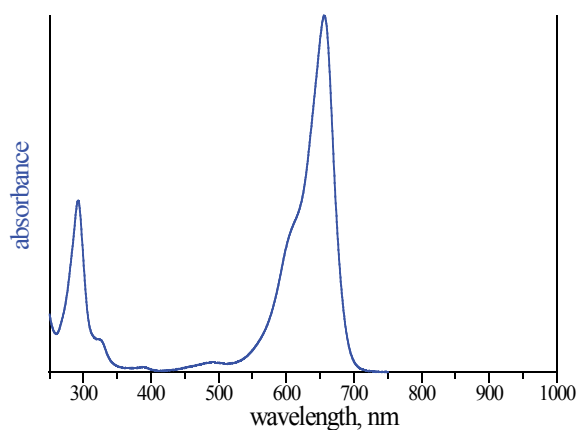
**Concentration:** 1.0 mM - 1.1 mM

**pH:** 7.5 ± 0.5

**Spectroscopic Properties:** λ<sub>abs</sub> 668 nm, ε 100.0 L mmol<sup>-1</sup> cm<sup>-1</sup> (Tris-HCl pH 7.5)



Structural formula of 5-Propargylamino-ddCTP-ATTO-MB2



absorption spectrum of ATTO MB2