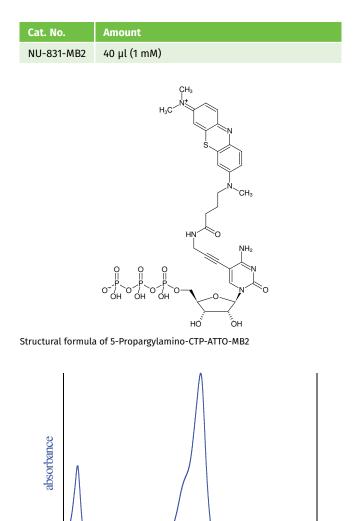




5-Propargylamino-CTP-ATTO-MB2

5-Propargylamino-cytidine-5'-triphosphate, labeled with ATTO-MB2, Triethylammonium salt



300 500 600 700 900 400 800 1000 wavelength, nm

absorption spectrum of ATTO MB2

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: C₃₁H₃₈N₇O₁₅P₃S (free acid)

Molecular Weight: 873.66 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_{abs} 668 nm, ϵ 100.0 L mmol $^{-1}$ cm $^{-1}$ (Tris-HCl pH 7.5)

