

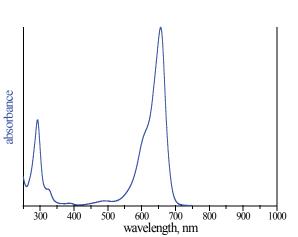


EDA-GTP-ATTO-MB2

2'/3'-O-(2-Aminoethyl-carbamoyl)-Guanosine-5'-triphosphate, labeled with ATTO-MB2, Triethylammonium salt

Cat. No.	Amount
NU-820-MB2	40 μl (1 mM)
	$\begin{array}{c} \begin{array}{c} & & \\ $

Structural formula of EDA-GTP-ATTO-MB2



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: $C_{32}H_{41}N_{10}O_{16}P_3S$ (free acid)

Molecular Weight: 946.71 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)

