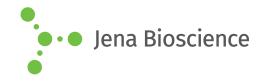
DATA SHEET



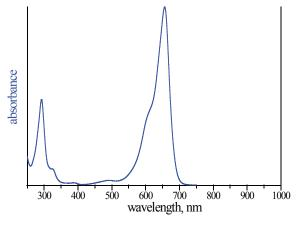


EDA-AppNHp (EDA-AMPPNP)-ATTO-MB2

2'/3'-O-(2-Aminoethyl-carbamoyl)-Adenosine-5'-[(β , γ)-imido] triphosphate, labeled with ATTO-MB2, Triethylammonium salt

Cat. No.	Amount
NU-810-MB2	40 μl (1 mM)

Structural formula of EDA-AppNHp (EDA-AMPPNP)-ATTO-MB2



absorption spectrum of ATTO MB2

For general laboratory use.

Shipping: shipped on dry ice

Storage Conditions: store at -20 °C

Shelf Life: 6 months after date of delivery

Molecular Formula: $C_{32}H_{42}N_{11}O_{14}P_3S$ (free acid)

Molecular Weight: 929.73 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)