## **DATA SHEET**

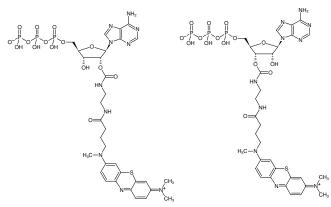




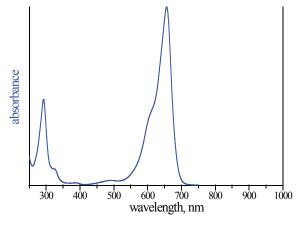
## **EDA-ATP-ATTO-MB2**

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

Cat. No.	Amount
NU-808-MB2	40 μl (1 mM)



Structural formula of EDA-ATP-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:**  $C_{32}H_{41}N_{10}O_{15}P_3S$  (free acid) **Molecular Weight:** 930.71 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:  $\lambda_{abs}$  668 nm,  $\epsilon$  100.0 L mmol  $^{\text{-}1}$  cm  $^{\text{-}1}$ 

(Tris-HCl pH 7.5)