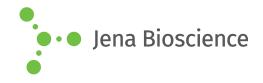
## **DATA SHEET**



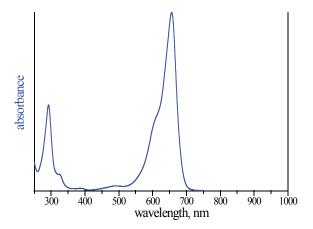


## γ-(6-Aminohexyl)-GTP-ATTO-MB2

γ-(6-Aminohexyl)-guanosine-5'-triphosphate, labeled with ATTO-MB2, Triethylammonium salt

Cat. No.	Amount
NU-834-MB2	40 μl (1 mM)

Structural formula of y-(6-Aminohexyl)-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<sub>35</sub>H<sub>48</sub>N<sub>9</sub>O<sub>15</sub>P<sub>3</sub>S (free acid) **Molecular Weight:** 959.79 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)