



### 8-(6-Aminoethyl)-amino-GTP-ATTO-MB2

8-(6-Aminoethyl)-amino-guanosine-5'-triphosphate, labeled with ATTO-MB2, Triethylammonium salt

Cat. No.	Amount
NU-830-MB2	40 µl (1 mM)

For general laboratory use.

Please centrifuge briefly before opening (volume ≤2 ml).

**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>49</sub>N<sub>10</sub>O<sub>15</sub>P<sub>3</sub>S (free acid)

**Molecular Weight:** 974.81 g/mol (free acid)

**Purity:** ≥ 95 % (HPLC)

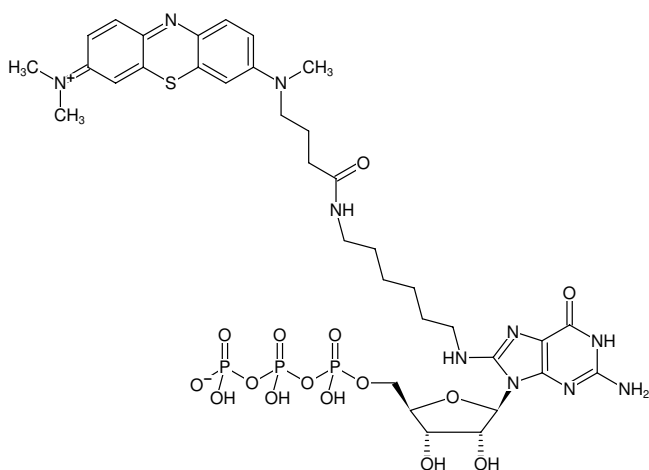
**Form:** solution in water

**Color:** blue

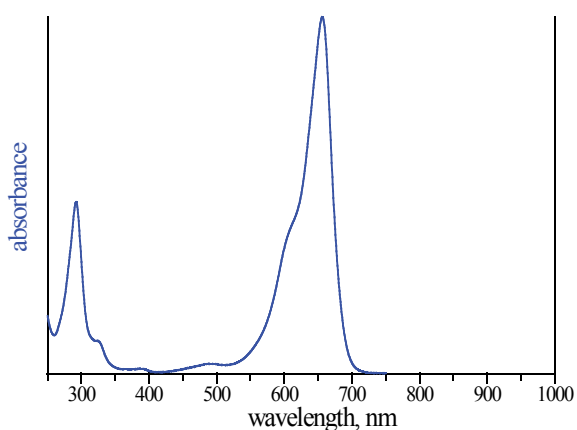
**Concentration:** 1.0 mM - 1.1 mM

**pH:** 7.5 ±0.5

**Spectroscopic Properties:** λ<sub>abs</sub> 668 nm, ε 100.0 L mmol<sup>-1</sup> cm<sup>-1</sup> (Tris-HCl pH 7.5)



Structural formula of 8-(6-Aminoethyl)-amino-GTP-ATTO-MB2



absorption spectrum of ATTO MB2