



8-(6-Aminoethyl)-amino-GMP-ATTO-655

8-(6-Aminoethyl)-amino-guanosine-5'-monophosphate, labeled with ATTO 655, Triethylammonium salt

Cat. No.	Amount
NU-829-655	80 µl (1 mM)

For research use only!

Shipping: shipped on blue ice

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: C₄₃H₅₉N₁₀O₁₃PS (free acid)

Molecular Weight: 987.04 g/mol (free acid)

Exact Mass: 986.37 g/mol (free acid)

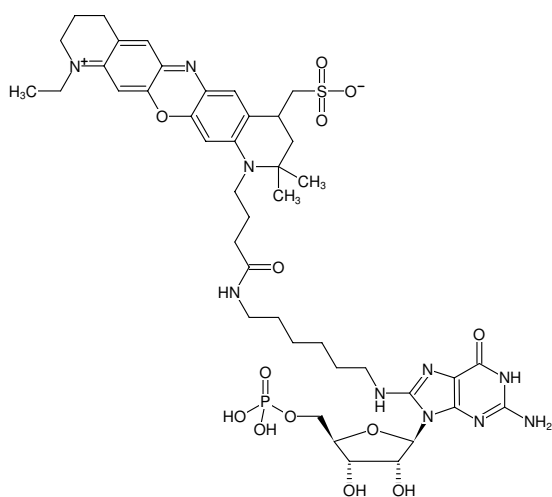
Purity: ≥ 95 % (HPLC)

Form: solution in water

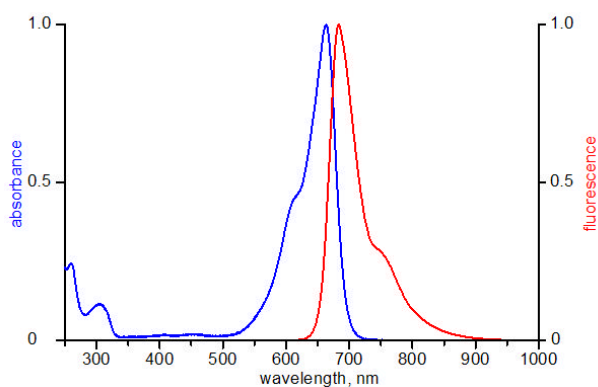
Concentration: 1.0 mM - 1.1 mM

pH: 7.5 ± 0.5

Spectroscopic Properties: λ_{exc} 663 nm, λ_{em} 680 nm, ε 125.0 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)



Structural formula of 8-(6-Aminoethyl)-amino-GMP-ATTO-655



excitation and emission spectrum of ATTO 655