



8-(6-Aminoethyl)-amino-cGMP-ATTO-590

8-(6-Aminoethyl)-amino-guanosine-3',5'-cyclic monophosphate, labeled with ATTO 590

Cat. No.	Amount
NU-832-590	120 µl (1 mM)

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: C₅₃H₆₂N₉O₁₁P (free acid)

Molecular Weight: 1032.10 g/mol (free acid)

Purity: ≥ 95 % (HPLC)

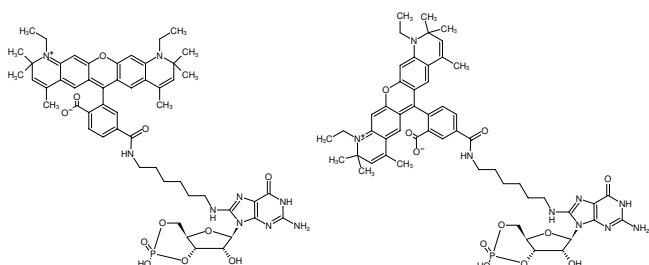
Form: solution in water

Color: red-violet

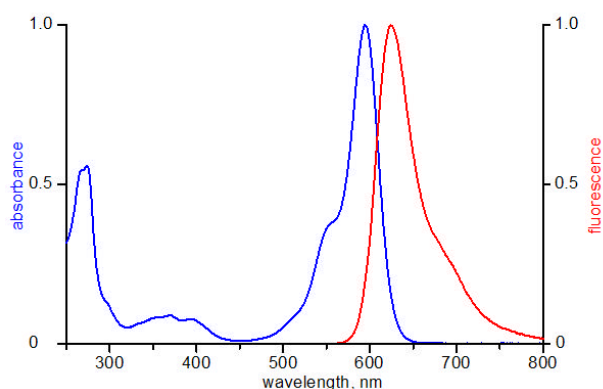
Concentration: 1.0 mM - 1.1 mM

pH: 7.5 ± 0.5

Spectroscopic Properties: λ_{exc} 593 nm, λ_{em} 622 nm, ε 120.0 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)



Structural formula of 8-(6-Aminoethyl)-amino-cGMP-ATTO-590



excitation and emission spectrum of ATTO 590