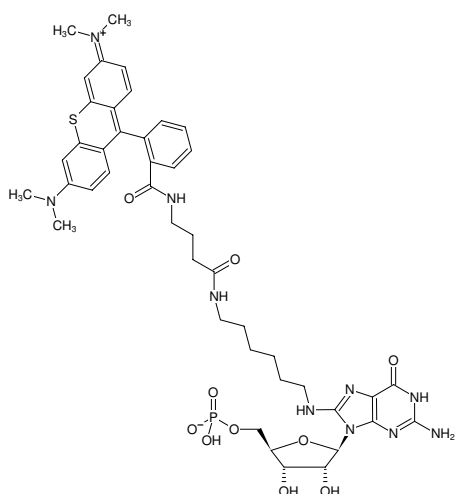




8-(6-Aminoethyl)-amino-GMP-ATTO-Thio12

8-(6-Aminoethyl)-amino-guanosine-5'-monophosphate, labeled with ATTO Thio12

Cat. No.	Amount
NU-829-THIO12	80 μ l (1 mM)



Structural formula of 8-(6-Aminoethyl)-amino-GMP-ATTO-Thio12

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: C₄₅H₅₇N₁₀O₁₀PS (free acid)

Molecular Weight: 961.04 g/mol (free acid)

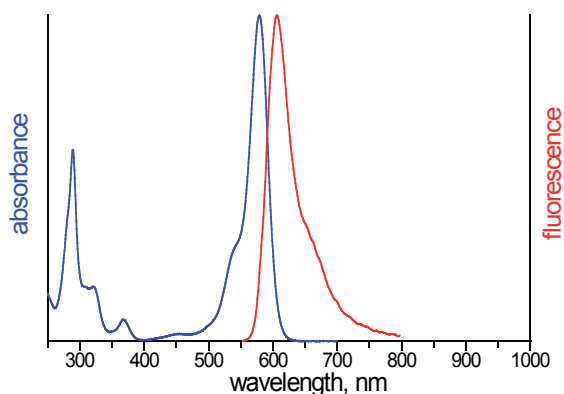
Purity: \geq 95 % (HPLC)

Form: solution in water

Concentration: 1.0 mM - 1.1 mM

pH: 7.5 \pm 0.5

Spectroscopic Properties: λ_{exc} 582 nm, λ_{em} 607 nm, ϵ 110.0 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)



excitation and emission spectrum of ATTO Thio12