



8-(6-Aminoethyl)-amino-GMP-ATTO-425

8-(6-Aminoethyl)-amino-guanosine-5'-monophosphate, labeled with ATTO 425, Triethylammonium salt

Cat. No.	Amount
NU-829-425	120 µl (1 mM)

For research use only!

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: C₃₈H₅₃N₈O₁₃P (free acid)

Molecular Weight: 860.85 g/mol (free acid)

Purity: ≥ 95 % (HPLC)

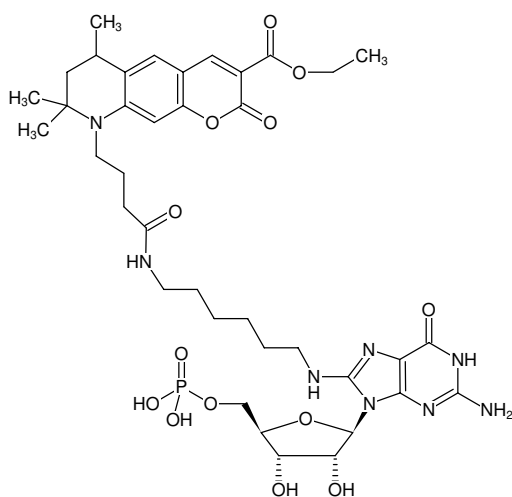
Form: solution in water

Color: green-yellow

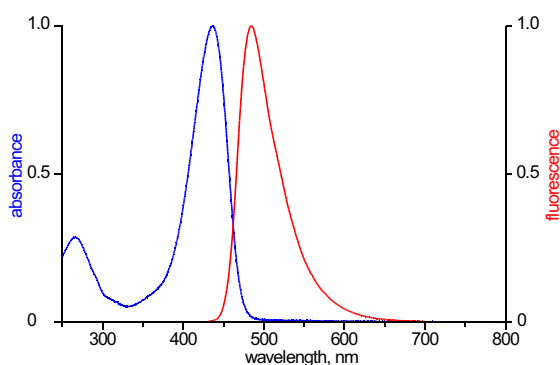
Concentration: 1.0 mM - 1.1 mM

pH: 7.5 ± 0.5

Spectroscopic Properties: λ_{exc} 439 nm, λ_{em} 485 nm, ε 45.0 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)



Structural formula of 8-(6-Aminoethyl)-amino-GMP-ATTO-425



excitation and emission spectrum of ATTO 425