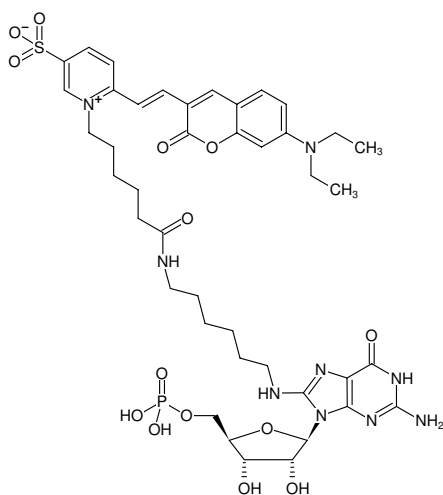




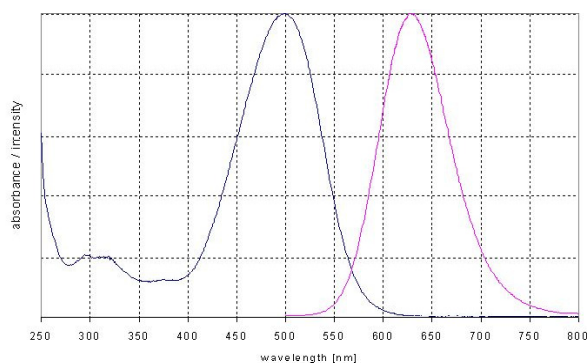
8-(6-Aminoethyl)-amino-GMP-DY-480XL

8-(6-Aminoethyl)-amino-guanosine-5'-monophosphate, labeled with DY 480XL, Triethylammonium salt

Cat. No.	Amount
NU-829-480XL	80 µl (1 mM)



Structural formula of 8-(6-Aminoethyl)-amino-GMP-DY-480XL



excitation and emission spectrum of DY 480XL

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: C₄₂H₅₆N₉O₁₄PS (free acid)

Molecular Weight: 973.99 g/mol (free acid)

Purity: ≥ 95 % (HPLC)

Form: clear aqueous solution

Concentration: 1.0 mM - 1.1 mM

pH: 7.5 ± 0.5

Spectroscopic Properties: λ_{exc} 500 nm, λ_{em} 630 nm, ε 50.0 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)