



### 8-(6-Aminoethyl)-amino-GMP-Cy5

8-(6-Aminoethyl)-amino-guanosine-5'-monophosphate, labeled with Cy5, Triethylammonium salt

Cat. No.	Amount
NU-829-CY5	160 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:** C<sub>48</sub>H<sub>64</sub>N<sub>9</sub>O<sub>15</sub>PS<sub>2</sub> (free acid)

**Molecular Weight:** 1102.18 g/mol (free acid)

**Purity:** ≥ 95 % (HPLC)

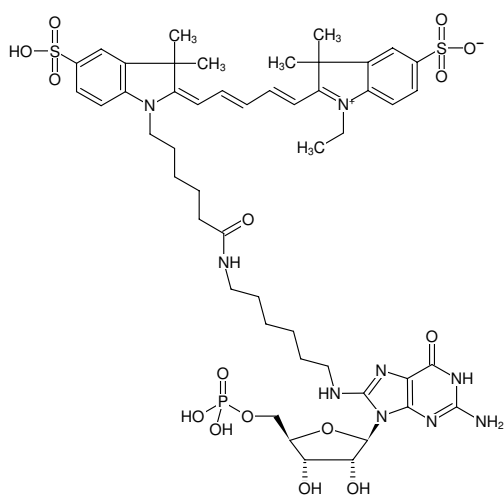
**Form:** solution in water

**Color:** blue

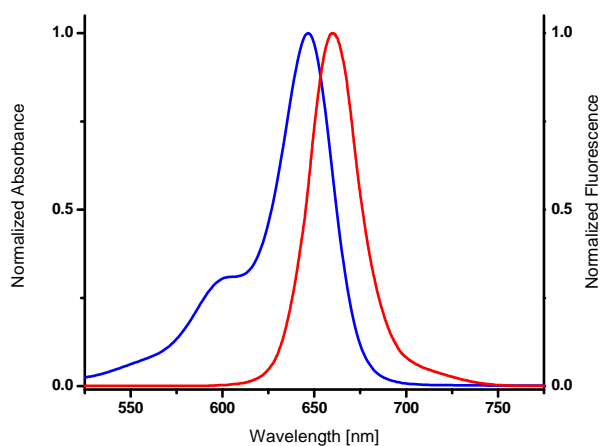
**Concentration:** 1.0 mM - 1.1 mM

**pH:** 7.5 ± 0.5

**Spectroscopic Properties:** λ<sub>exc</sub> 649 nm, λ<sub>em</sub> 670 nm, ε 250.0 L mmol<sup>-1</sup> cm<sup>-1</sup> (Tris-HCl pH 7.5)



Structural formula of 8-(6-Aminoethyl)-amino-GMP-Cy5



excitation and emission spectrum of Cy5