



### 8-(6-Aminoethyl)-amino-GMP-5/6-TAMRA

8-(6-Aminoethyl)-amino-guanosine-5'-monophosphate, labeled with 5/6-TAMRA

Cat. No.	Amount
NU-829-TAM	160 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:** C<sub>41</sub>H<sub>48</sub>N<sub>9</sub>O<sub>16</sub>P (free acid)

**Molecular Weight:** 889.85 g/mol (free acid)

**Purity:** ≥ 95 % (HPLC)

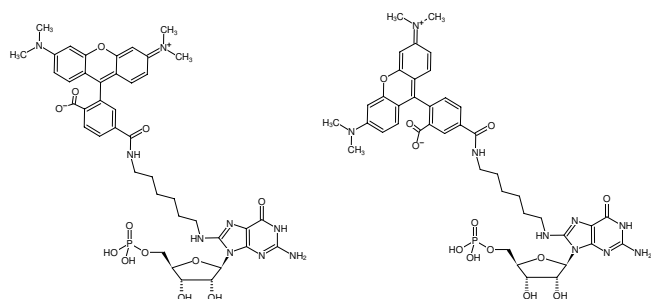
**Form:** solution in water

**Color:** pink to red

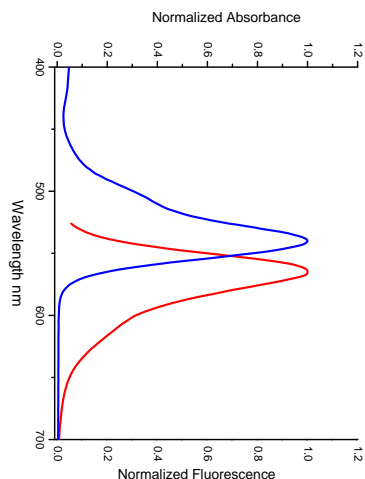
**Concentration:** 1.0 mM - 1.1 mM

**pH:** 7.5 ± 0.5

**Spectroscopic Properties:** λ<sub>exc</sub> 545 nm, λ<sub>em</sub> 575 nm, ε 90.0 L mmol<sup>-1</sup> cm<sup>-1</sup> (Tris-HCl pH 7.5)



Structural formula of 8-(6-Aminoethyl)-amino-GMP-5/6-TAMRA



excitation and emission spectrum of 5/6-TAMRA