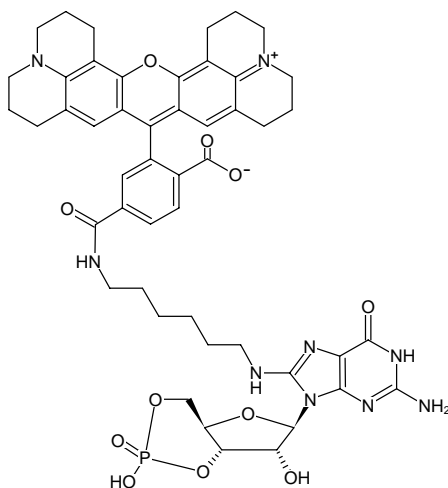




## 8-[(6-Amino)hexyl]-amino-cGMP - 6-ROX

8-[(6-Amino)hexyl]-amino-guanosine-3',5'-cyclic monophosphate, labeled with 6-ROX

Cat. No.	Amount
NU-832-ROX	160 µl (1 mM)



Structural formula of 8-[(6-Amino)hexyl]-amino-cGMP - 6-ROX

**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>49</sub>H<sub>54</sub>N<sub>9</sub>O<sub>11</sub>P (free acid)

**Molecular Weight:** 975.99 g/mol (free acid)

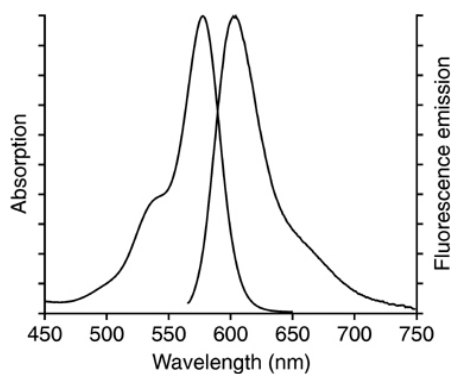
**Purity:** ≥ 95 % (HPLC)

**Form:** clear aqueous solution

**Concentration:** 1 mM - 1.1 mM

**pH:** 7.5 ± 0.5

**Spectroscopic Properties:** λ<sub>exc</sub> 575 nm; λ<sub>em</sub> 600 nm; ε 82.0 L mmol<sup>-1</sup> cm<sup>-1</sup> (Tris-HCl pH 7.5)



excitation and emission spectrum of 6-ROX