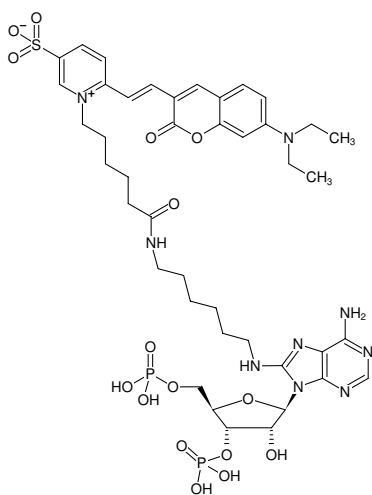




## 8-(6-Aminoethyl)-amino-adenosine-3',5'-bisphosphate-DY-480XL

8-(6-Aminoethyl)-amino-adenosine-3',5'-bisphosphate, labeled with DY 480XL, Triethylammonium salt

Cat. No.	Amount
NU-811-480XL	40 µl (1 mM)



Structural formula of 8-(6-Aminoethyl)-amino-adenosine-3',5'-bisphosphate-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:** C<sub>42</sub>H<sub>57</sub>N<sub>9</sub>O<sub>16</sub>P<sub>2</sub>S (free acid)

**Molecular Weight:** 1037.97 g/mol (free acid)

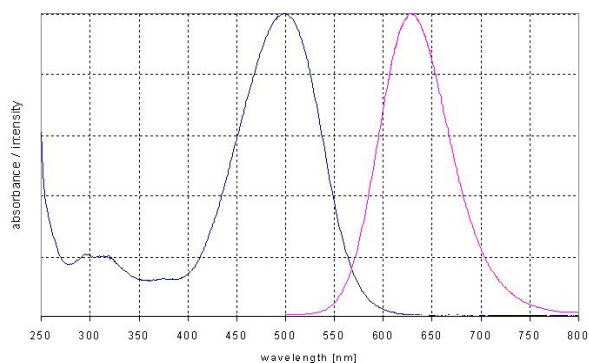
**Purity:** ≥ 95 % (HPLC)

**Form:** clear aqueous solution

**Concentration:** 1.0 mM - 1.1 mM

**pH:** 7.5 ± 0.5

**Spectroscopic Properties:** λ<sub>exc</sub> 500 nm, λ<sub>em</sub> 630 nm, ε 50.0 L mmol<sup>-1</sup> cm<sup>-1</sup> (Tris-HCl pH 7.5)



excitation and emission spectrum of DY 480XL