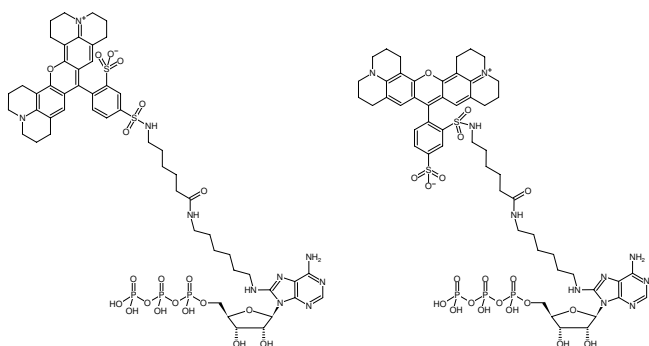




### 8-(6-Aminoethyl)-amino-ATP-Texas Red

8-(6-Aminoethyl)-amino-adenosine-5'-triphosphate, labeled with Texas Red, Triethylammonium salt

Cat. No.	Amount
NU-807-TXR	160 $\mu$ l (1 mM)



Structural formula of 8-(6-Aminoethyl)-amino-ATP-Texas Red

**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>53</sub>H<sub>69</sub>N<sub>10</sub>O<sub>20</sub>P<sub>3</sub>S<sub>2</sub> (free acid)

**Molecular Weight:** 1323.22 g/mol (free acid)

**Purity:**  $\geq$  95 % (HPLC)

**Form:** red-violet solution in water

**Concentration:** 1.0 mM - 1.1 mM

**pH:** 7.5  $\pm$  0.5

**Spectroscopic Properties:**  $\lambda_{exc}$  588 nm,  $\lambda_{em}$  609 nm,  $\epsilon$  80.0 L mmol<sup>-1</sup> cm<sup>-1</sup> (Tris-HCl pH 7.5)