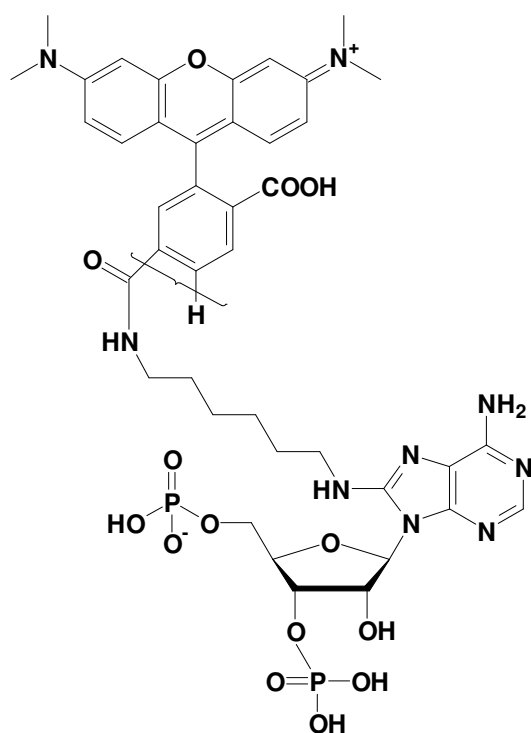


## 8-[(6-Amino)hexyl]-amino-adenosine-3'/5'-bisphosphate - 5/6-TAMRA 8-[(6-Amino)hexyl]-amino-adenosine-3'/5'-bisphosphate, labeled with 5/6-TAMRA, Triethylammonium salt

Cat. No.	Amount
NU-811-TAM	160 $\mu$ l (1 mM)



structural formula of 8-[(6-Amino)hexyl]-amino-adenosine-3',5'-bisphosphate - 5/6-TAMRA

**Cat. No.:** NU-811-TAM

**Molecular Formula:** C<sub>41</sub>H<sub>49</sub>N<sub>9</sub>O<sub>14</sub>P<sub>2</sub> (free acid)

**Molecular Weight:** 953.83 (free acid)

**Purity:** > 95%, clear aqueous solution, pH 7.5

**Spectroscopic properties:**

$\lambda_{exc}$  545 nm;  $\lambda_{em}$  575 nm;  $\epsilon$  90,000 cm<sup>-1</sup> M<sup>-1</sup>

**Storage conditions:**

Short term exposure (up to 1 week cumulative) to ambient temperature possible. Long term storage at < -20°C. If stored as recommended, Jena Bioscience guarantees optimal performance of this product for 12 months after date of delivery.

**For research use only!**