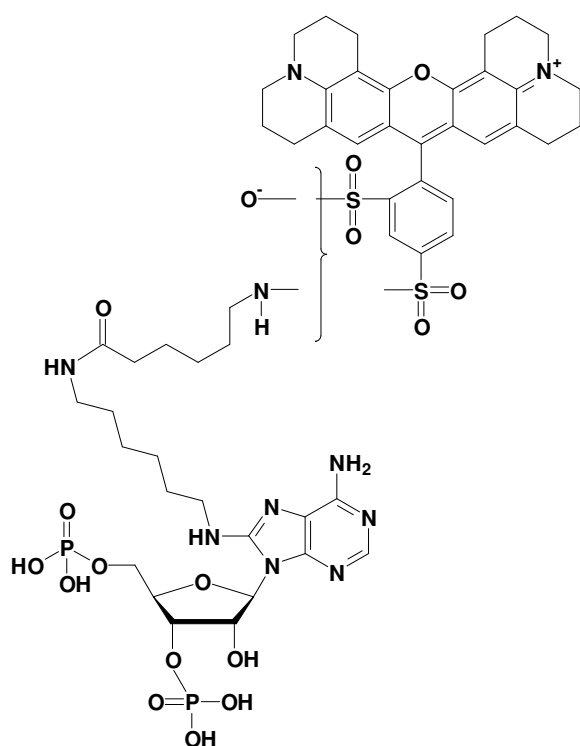


8-[(6-Amino)hexyl]-amino-adenosine-3'/5'-bisphosphate - Texas Red

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

Cat. No.	Amount
NU-811-TXR	160 μ l (1 mM)



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

Cat. No.: NU-811-TXR

Molecular Formula: C₅₃H₆₈N₁₀O₁₇P₂S₂ (free acid)

Molecular Weight: 1243,24 (free acid)

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

Spectroscopic properties:

λ_{exc} 588 nm; λ_{em} 609 nm; ϵ 80.000 cm⁻¹ M⁻¹

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!