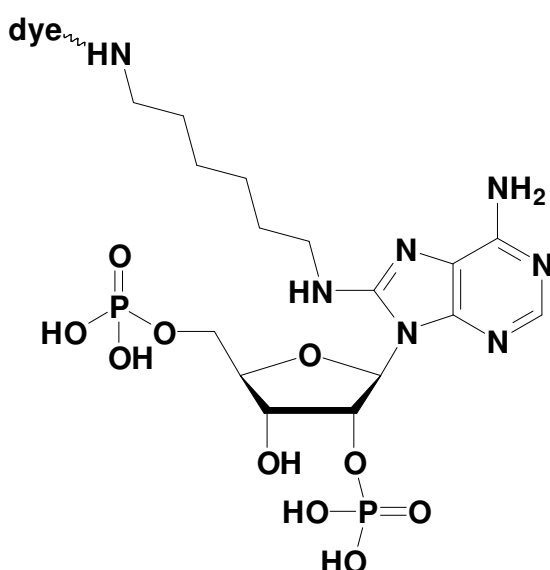


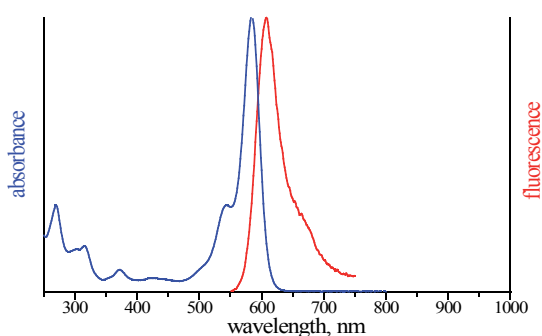
## 8-[(6-Amino)hexyl]-amino-adenosine-2'/5'-bisphosphate - ATTO Rho101

8-[(6-Amino)hexyl]-amino-adenosine-2'/5'-bisphosphate, labeled with ATTO Rho101,  
 Triethylammonium salt

Cat. No.	Amount
NU-812-RHO101	40 $\mu$ l (1 mM)



structural formula of 8-[(6-Amino)hexyl]-amino-adenosine-2',5'-bisphosphate



excitation and emission spectrum of ATTO Rho101

**Cat. No.:** NU-812-RHO101

**Molecular Formula:** C<sub>16</sub>H<sub>29</sub>N<sub>7</sub>O<sub>10</sub>P<sub>2</sub> - ATTO Rho101  
 (free acid)

**Molecular Weight:** 1112,39 (free acid)

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

**Spectroscopic properties:**

$\lambda_{exc}$  586 nm;  $\lambda_{em}$  610 nm;  $\epsilon$  1,2 x 10<sup>5</sup> 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!**