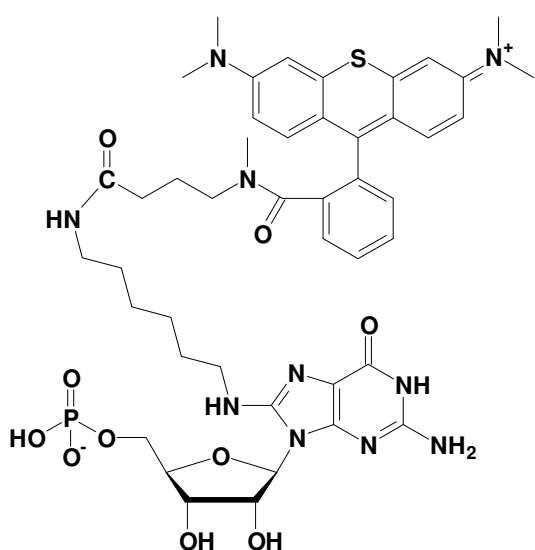


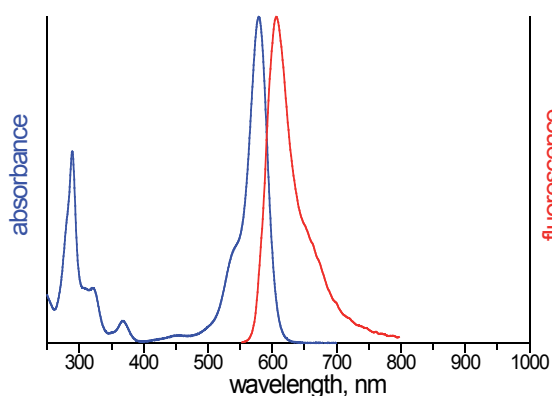
## 8-[(6-Amino)hexyl]-amino-GMP - ATTO Thio12

### 8-[(6-Amino)hexyl]-amino-guanosine-5'-monophosphate, labeled with ATTO Thio12

Cat. No.	Amount
NU-829-THIO12	80 $\mu$ l (1 mM)



structural formula of 8-[(6-Amino)hexyl]-amino-GMP - ATTO Thio12



excitation and emission spectrum of ATTO Thio12

**Cat. No.:** NU-829-THIO12

**Molecular Formula:** C<sub>45</sub>H<sub>57</sub>N<sub>10</sub>O<sub>10</sub>PS (free acid)

**Molecular Weight:** 961,04 (free acid)

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

**Spectroscopic properties:**

$\lambda_{exc}$  579 nm;  $\lambda_{em}$  609 nm;  $\epsilon$  1,1 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!**