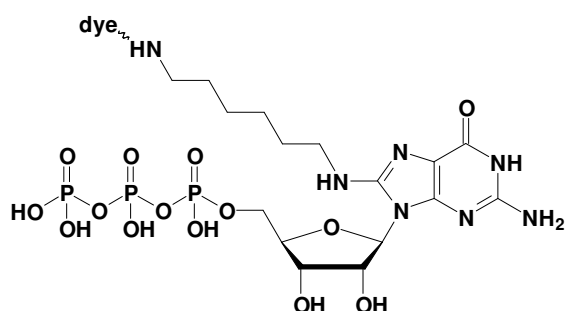


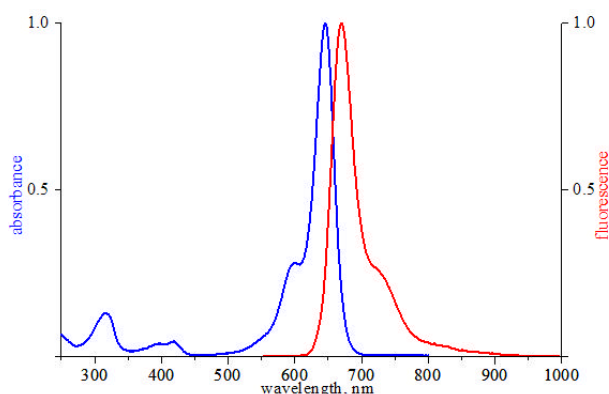
## 8-[(6-Amino)hexyl]-amino-GTP - ATTO 647N

8-[(6-Amino)hexyl]-amino-guanosine-5'-triphosphate, labeled with ATTO 647N,  
 Triethylammonium salt

Cat. No.	Amount
NU-830-647N	80 $\mu$ l (1 mM)



structural formula of 8-[(6-Amino)hexyl]-amino-GTP



excitation and emission spectrum of ATTO 647N

**Cat. No.:** NU-830-647N

**Molecular Formula:**  $C_{16}H_{30}N_7O_{14}P_3$  - ATTO 647N  
 (free acid)

**Molecular Weight:** 1264.37 (free acid)

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

**Spectroscopic properties:**

$\lambda_{exc}$  644 nm;  $\lambda_{em}$  669 nm;  $\epsilon$  150,000  $cm^{-1} M^{-1}$

**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!**