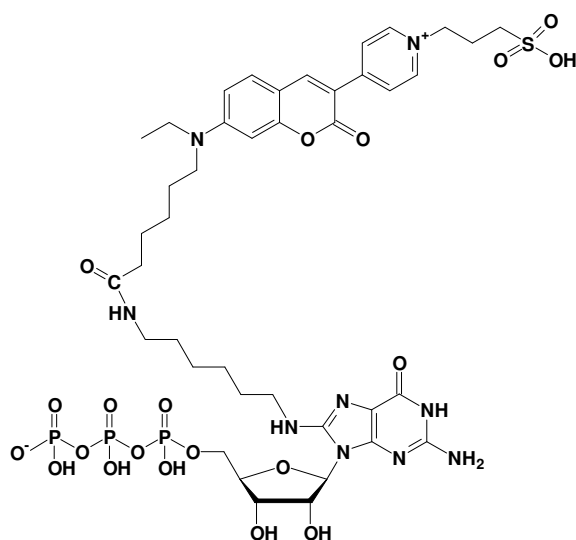


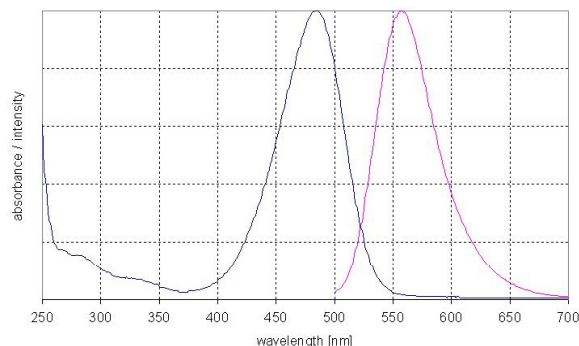
## 8-[(6-Amino)hexyl]-amino-GTP - DY 485XL

8-[(6-Amino)hexyl]-amino-guanosine-5'-triphosphate, labeled with DY 485XL,  
 Triethylammonium salt

Cat. No.	Amount
NU-830-485XL	80 $\mu$ l (1 mM)



structural formula of 8-[(6-Amino)hexyl]-amino-GTP - DY 485XL



excitation and emission spectrum of DY 485XL

**Cat. No.:** NU-830-485XL

**Molecular Formula:** C<sub>41</sub>H<sub>58</sub>N<sub>9</sub>O<sub>20</sub>P<sub>3</sub>S (free acid)

**Molecular Weight:** 1121,94 (free acid)

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

**Spectroscopic properties:**

$\lambda_{exc}$  485 nm;  $\lambda_{em}$  560 nm;  $\epsilon$  50,000 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!**