

## 8-[(6-Amino)hexyl]-amino-GTP - Texas Red

8-[(6-Amino)hexyl]-amino-guanosine-5'-triphosphate, labeled with Texas Red,  
 Triethylammonium salt

Cat. No.	Amount
NU-830-TXR	160 $\mu$ l (1 mM)

**Cat. No.:** NU-830-TXR

**Molecular Formula:** C<sub>53</sub>H<sub>69</sub>N<sub>10</sub>O<sub>21</sub>P<sub>3</sub>S<sub>2</sub> (free acid)

**Molecular Weight:** 1339,22 (free acid)

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

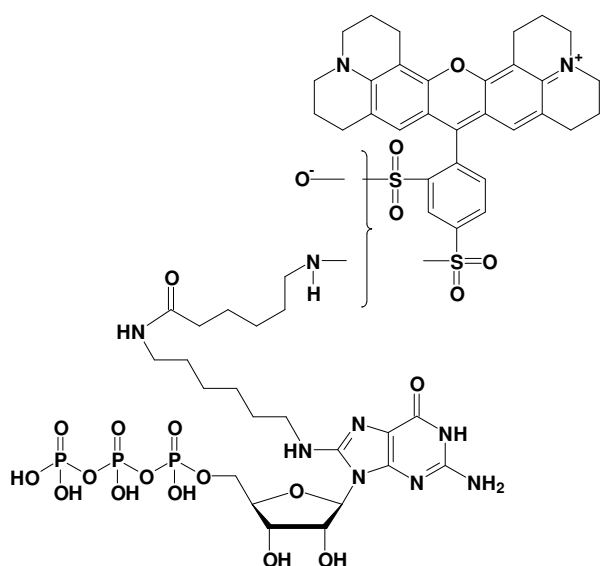
**Spectroscopic properties:**

$\lambda_{exc}$  588 nm;  $\lambda_{em}$  609 nm;  $\epsilon$  80.000 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!**



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