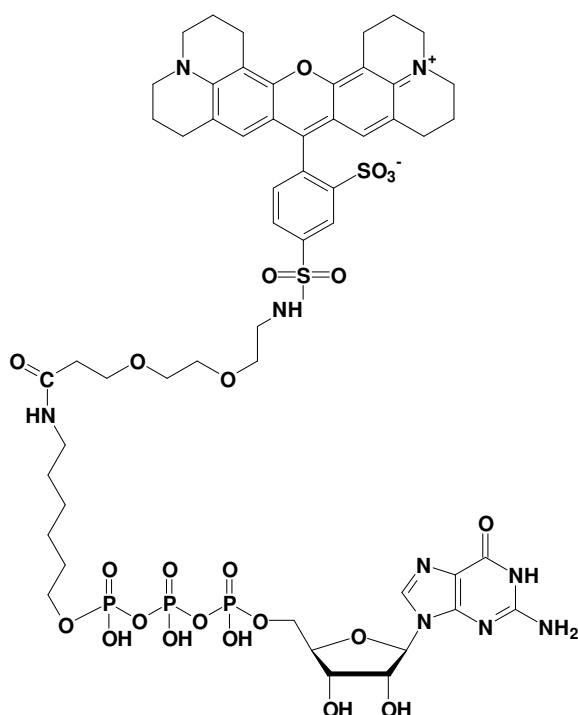


γ -[6-Aminoethyl]-GTP - Texas Red

γ -[6-Aminoethyl]-guanosine-5'-triphosphate, labeled with Texas Red,
 Triethylammonium salt

Cat. No.	Amount
NU-834-TXR	160 μ l (1 mM)



structural formula of γ -[6-Aminoethyl]-GTP - Texas Red

Cat. No.: NU-834-TXR

Molecular Formula: C₅₄H₇₀N₉O₂₃P₃S₂ (free acid)

Molecular Weight: 1370,23 (free acid)

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

Spectroscopic properties:

λ_{exc} 588 nm; λ_{em} 609 nm; ϵ 80.000 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!