



## $\gamma$ -(6-Aminoethyl)-GTP-ATTO-Rho12

$\gamma$ -(6-Aminoethyl)-guanosine-5'-triphosphate, labeled with ATTO Rho12, Triethylammonium salt

Cat. No.	Amount
NU-834-RHO12	80 $\mu$ l (1 mM)

**For general laboratory use.**

**Shipping:** shipped on gel packs

**Storage Conditions:** store at -20 °C

Short term exposure (up to 1 week cumulative) to ambient temperature possible.

**Shelf Life:** 12 months after date of delivery

**Molecular Formula:** C<sub>57</sub>H<sub>78</sub>N<sub>9</sub>O<sub>17</sub>P<sub>3</sub> (free acid)

**Molecular Weight:** 1254.22 g/mol (free acid)

**Exact Mass:** 1253.47 g/mol (free acid)

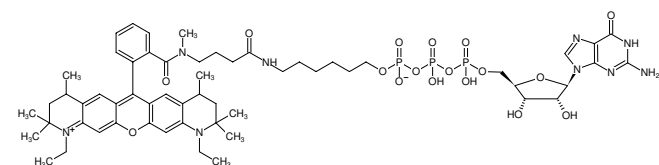
**Purity:**  $\geq$  95 % (HPLC)

**Form:** solution in water

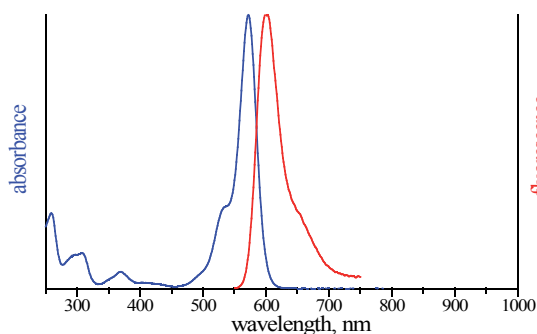
**Concentration:** 1.0 mM - 1.1 mM

**pH:** 7.5  $\pm$  0.5

**Spectroscopic Properties:**  $\lambda_{exc}$  576 nm,  $\lambda_{em}$  601 nm,  $\epsilon$  120.0 L mmol<sup>-1</sup> cm<sup>-1</sup> (Tris-HCl pH 7.5)



Structural formula of  $\gamma$ -(6-Aminoethyl)-GTP-ATTO-Rho12



excitation and emission spectrum of ATTO Rho12