

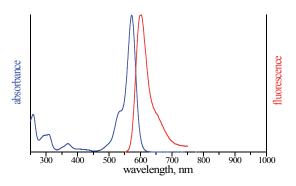


## 8-(6-Aminohexyl)-amino-GTP-ATTO-Rho12

8-(6-Aminohexyl)-amino-guanosine-5'-triphosphate, labeled with ATTO Rho12, Triethylammonium salt

Cat. No.	Amount
NU-830-RHO12	80 μl (1 mM)

Structural formula of 8-(6-Aminohexyl)-amino-GTP-ATTO-Rho12



excitation and emission spectrum of ATTO Rho12

## 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>79</sub>N<sub>10</sub>O<sub>17</sub>P<sub>3</sub> (free acid) **Molecular Weight:** 1269.23 g/mol (free acid)

Exact Mass: 1268.48 g/mol (free acid)

**Purity:** ≥ 95 % (HPLC) **Form:** solution in water

Concentration: 1.0 mM - 1.1 mM

**pH:** 7.5 ±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)