



## EDA-m<sup>7</sup>GTP-ATTO-Rho6G

2'/3'-O-(2-Aminoethyl-carbamoyl)-7-methyl-guanosine-5'-triphosphate, labeled with ATTO Rho6G, Triethylammonium salt

Cat. No.	Amount
NU-824-RHO6G	40 µl (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>45</sub>H<sub>57</sub>N<sub>10</sub>O<sub>18</sub>P<sub>3</sub> (free acid)

**Molecular Weight:** 1118.92 g/mol (free acid)

**Exact Mass:** 1118.31 g/mol (free acid)

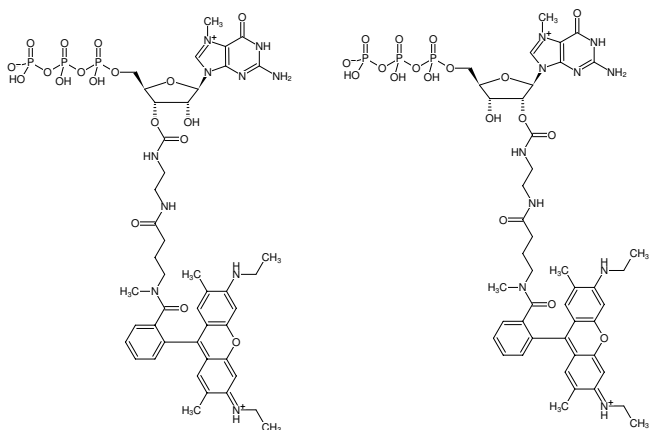
**Purity:** ≥ 95 % (HPLC)

**Form:** solution in water

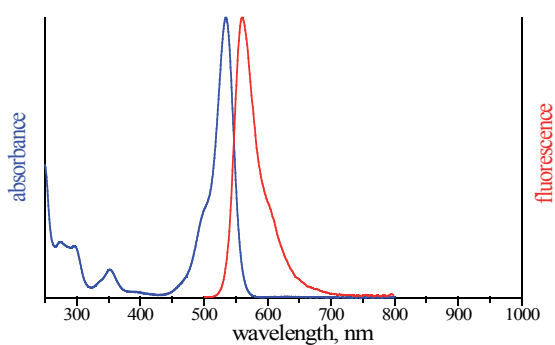
**Concentration:** 1.0 mM - 1.1 mM

**pH:** 7.5 ± 0.5

**Spectroscopic Properties:** λ<sub>exc</sub> 533 nm, λ<sub>em</sub> 557 nm, ε 115.0 L mmol<sup>-1</sup> cm<sup>-1</sup> (Tris-HCl pH 7.5)



Structural formula of EDA-m<sup>7</sup>GTP-ATTO-Rho6G



excitation and emission spectrum of ATTO Rho6G