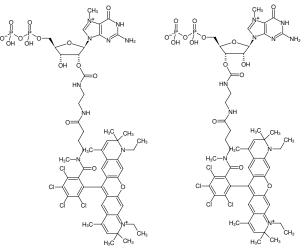




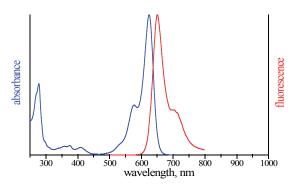
## EDA-m<sup>7</sup>GDP-ATTO-Rho14

2'/3'-O-(2-Aminoethyl-carbamoyl)-7-methyl-guanosine-5'-diphosphate, labeled with ATTO Rho14, Triethylammonium salt

Cat. No.	Amount
NU-827-RHO14	40 μl (1 mM)



Structural formula of EDA-m<sup>7</sup>GDP-ATTO-Rho14



excitation and emission spectrum of ATTO Rho14

## 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:  $C_{55}H_{64}Cl_4N_{10}O_{15}P_2$  (free acid)

Molecular Weight: 1308.92 g/mol (free acid)

Exact Mass: 1306.28 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:  $\lambda_{exc}$  626 nm,  $\lambda_{em}$  646 nm,  $\epsilon$  140.0 L mmol<sup>-1</sup> cm<sup>-1</sup> (Tris-HCl pH 7.5)

