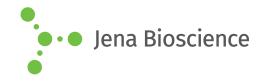
DATA SHEET



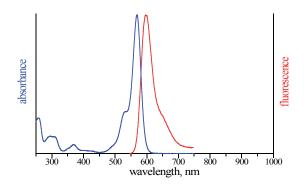


8-(6-Aminohexyl)-amino-cGMP-ATTO-Rho11

8-(6-Aminohexyl)-amino-guanosine-3',5'-cyclic monophosphate, labeled with ATTO Rho11, Triethylammonium salt

Cat. No.	Amount
NU-832-RHO11	80 μl (1 mM)

Structural formula of 8-(6-Aminohexyl)-amino-cGMP-ATTO-Rho11



excitation and emission spectrum of ATTO Rho11

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: $C_{51}H_{63}N_{10}O_{10}P$ (free acid) Molecular Weight: 1007.10 g/mol (free acid)

Exact Mass: 1006.45 g/mol (free acid)

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

Concentration: 1.0 mM - 1.1 mM

Spectroscopic Properties: λ_{exc} 572 nm, λ_{em} 595 nm,

ε 120.0 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)