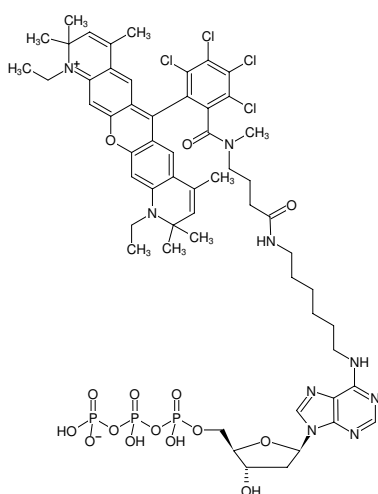




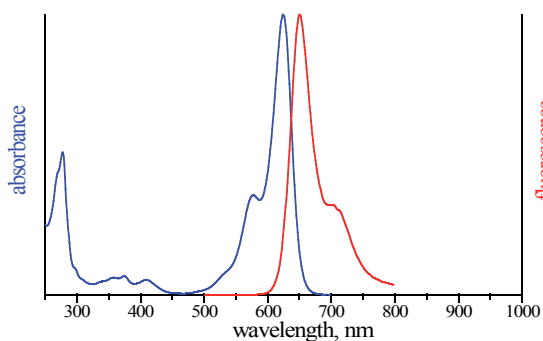
## N<sup>6</sup>-(6-Aminohexyl)-dATP-ATTO-Rho14

N<sup>6</sup>-(6-Aminohexyl)-2'-deoxyadenosine-5'-triphosphate, labeled with ATTO Rho14, Triethylammonium salt

| Cat. No.     | Amount       |
|--------------|--------------|
| NU-835-RHO14 | 40 µl (1 mM) |



Structural formula of N<sup>6</sup>-(6-Aminohexyl)-dATP-ATTO-Rho14



excitation and emission spectrum of ATTO Rho14

**For general laboratory use.**

**Please centrifuge briefly before opening (volume ≤2 ml).**

**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>70</sub>Cl<sub>4</sub>N<sub>9</sub>O<sub>15</sub>P<sub>3</sub> (free acid)

**Molecular Weight:** 1355.96 g/mol (free acid)

**Exact Mass:** 1353.30 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> 626 nm, λ<sub>em</sub> 646 nm, ε 140.0 L mmol<sup>-1</sup> cm<sup>-1</sup> (Tris-HCl pH 7.5)