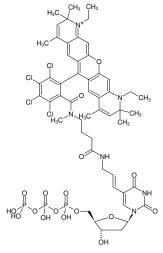




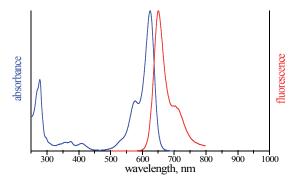
## Aminoallyl-dUTP-ATTO-Rho14

5-(3-Aminoallyl)-2'-deoxyuridine-5'-triphosphate, labeled with ATTO Rho14, Triethylammonium salt

Cat. No.	Amount
NU-803-RHO14	40 μl (1 mM)



Structural formula of Aminoallyl-dUTP-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<sub>53</sub>H<sub>61</sub>Cl<sub>4</sub>N<sub>6</sub>O<sub>17</sub>P<sub>3</sub> (free acid)

Molecular Weight: 1288.82 g/mol (free acid)

Exact Mass: 1286.21 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)

