DATA SHEET



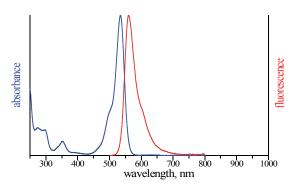


Aminoallyl-dUTP-ATTO-Rho6G

5-(3-Aminoallyl)-2'-deoxyuridine-5'-triphosphate, labeled with ATTO Rho6G, Triethylammonium salt

| Cat. No. | Amount |
|--------------|--------------|
| NU-803-RHO6G | 40 μl (1 mM) |

Structural formula of Aminoallyl-dUTP-ATTO-Rho6G



excitation and emission spectrum of ATTO Rho6G

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:** C₄₃H₅₃N₆O₁₇P₃ (free acid) **Molecular Weight:** 1018.84 g/mol (free acid)

Exact Mass: 1018.27 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_{exc} 533 nm, λ_{em} 557 nm,

 ϵ 115.0 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)