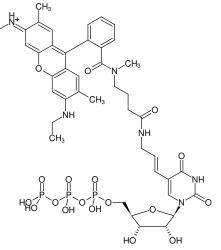




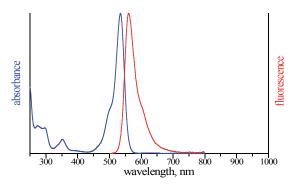
## Aminoallyl-UTP-ATTO-Rho6G

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

| Cat. No.     | Amount       |
|--------------|--------------|
| NU-821-RHO6G | 40 μl (1 mM) |



Structural formula of Aminoallyl-UTP-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:** C<sub>43</sub>H<sub>53</sub>N<sub>6</sub>O<sub>18</sub>P<sub>3</sub> (free acid)

Molecular Weight: 1034.84 g/mol (free acid)

Exact Mass: 1034.26 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:  $\lambda_{exc}$  533 nm,  $\lambda_{em}$  557 nm,  $\epsilon$  115.0 L mmol<sup>-1</sup> cm<sup>-1</sup> (Tris-HCl pH 7.5)

