



γ-(6-Aminohexyl)-GTP-ATTO-Thio12

 γ -(6-Aminohexyl)-guanosine-5'-triphosphate, labeled with ATTO Thio12, Triethylammonium salt

Cat. No.	Amount
NU-834-THIO12	80 μl (1 mM)
$H_{0}C_{-N}$	

Structural formula of γ -(6-Aminohexyl)-GTP-ATTO-Thio12



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: C45H58N9O16P3S (free acid)

Molecular Weight: 1105.98 g/mol (free acid)

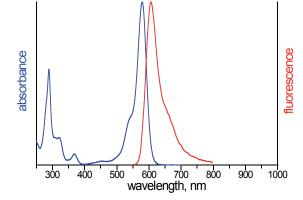
Purity: ≥ 95 % (HPLC)

Form: solution in water

Concentration: 1.0 mM - 1.1 mM

pH: 7.5 ±0.5

Spectroscopic Properties: λ_{exc} 582 nm, λ_{em} 607 nm, ϵ 110.0 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)



excitation and emission spectrum of ATTO Thio12

