



## γ-(6-Aminohexyl)-ATP-ATTO-Thio12

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

Cat. No.	Amount
NU-833-THIO12	800 μl (0,1 mM)
$H_3C_{-N}F$ $H_3C_{-N}F$ $H_3C_{-N}F$ $H_3C_{-N}F$ $CH_3$	$P_{HN} \rightarrow P_{HN} \rightarrow P$

Structural formula of  $\gamma$ -(6-Aminohexyl)-ATP-ATTO-Thio12

## 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:  $C_{45}H_{58}N_9O_{15}P_3S$  (free acid)

Molecular Weight: 1089.98 g/mol (free acid)

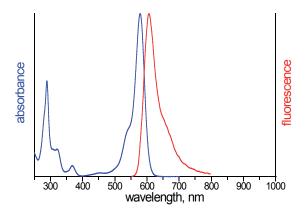
**Purity:** ≥ 95 % (HPLC)

Form: solution in water

Concentration: 0.10 mM - 0.11 mM

**pH:** 7.5 ±0.5

Spectroscopic Properties:  $\lambda_{exc}$  582 nm,  $\lambda_{em}$  607 nm,  $\epsilon$  110.0 L mmol<sup>-1</sup> cm<sup>-1</sup> (Tris-HCl pH 7.5)



excitation and emission spectrum of ATTO Thio12

