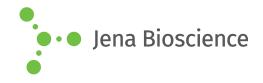
## **DATA SHEET**



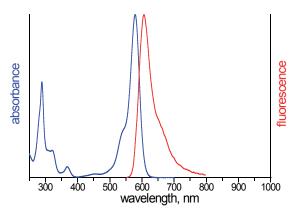


## N<sup>6</sup>-(6-Aminohexyl)-ATP-ATTO-Thio12

N<sup>6</sup>-(6-Aminohexyl)-adenosine-5'-triphosphate, labeled with ATTO Thio12, Triethylammonium salt

Cat. No.	Amount
NU-805-THIO12	80 μl (1 mM)

Structural formula of N<sup>6</sup>-(6-Aminohexyl)-ATP-ATTO-Thio12



excitation and emission spectrum of 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: 1.0 mM - 1.1 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)