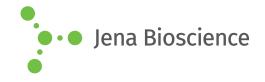
DATA SHEET



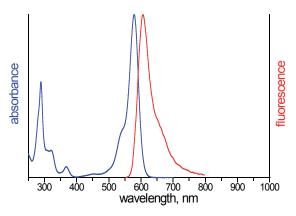


EDA-m⁷GDP-ATTO-Thio12

2'/3'-O-(2-Aminoethyl-carbamoyl)-7-methyl-guanosine-5'-diphosphate, labeled with ATTO Thio12

Cat. No.	Amount
NU-827-THIO12	40 μl (1 mM)

Structural formula of EDA- $\mathrm{m^7GDP}$ -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: $C_{43}H_{52}N_{10}O_{14}P_2S$ (free acid) Molecular Weight: 1026.94 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)