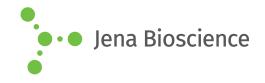
## **DATA SHEET**



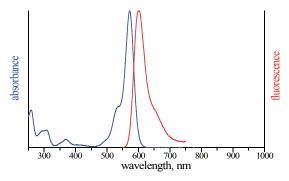


## EDA-m<sup>7</sup>GDP-ATTO-Rho12

2'/3'-O-(2-Aminoethyl-carbamoyl)-7-methyl-guanosine-5'-diphosphate, labeled with ATTO Rho12, Triethylammonium salt

Cat. No.	Amount
NU-827-RHO12	40 μl (1 mM)

Structural formula of EDA-m<sup>7</sup>GDP-ATTO-Rho12



excitation and emission spectrum of ATTO Rho12

## 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>55</sub>H<sub>72</sub>N<sub>10</sub>O<sub>15</sub>P<sub>2</sub> (free acid) **Molecular Weight:** 1175.18 g/mol (free acid)

Exact Mass: 1174.47 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:  $\lambda_{exc}$  576 nm,  $\lambda_{em}$  601 nm,

ε 120.0 L mmol<sup>-1</sup> cm<sup>-1</sup> (Tris-HCl pH 7.5)