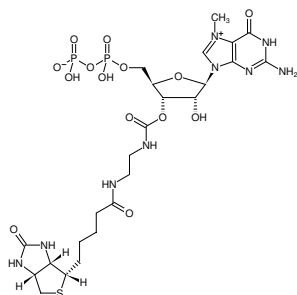
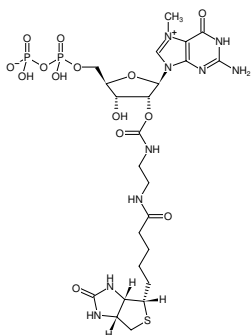


**EDA-m⁷GDP-Biotin**

2'/3'-O-(2-Aminoethyl-carbamoyl)-N7-methyl-guanosine-5'-diphosphate-Biotin, Triethylammonium salt

Cat. No.	Amount
NU-827-BIO	50 µl (5 mM)

Structural formula of EDA-m⁷GDP-Biotin**For research use only!****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:** C₂₄H₃₇N₉O₁₄P₂S (free acid)**Molecular Weight:** 769.62 g/mol (free acid)**Exact Mass:** 769.17 g/mol (free acid)**Purity:** ≥ 95 % (HPLC)**Form:** solution in water**Color:** colorless to slightly yellow**Concentration:** 5.0 mM - 5.5 mM**pH:** 7.5 ±0.5**Spectroscopic Properties:** λ_{max} 258/280 nm, ε 9.8/8.0 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)