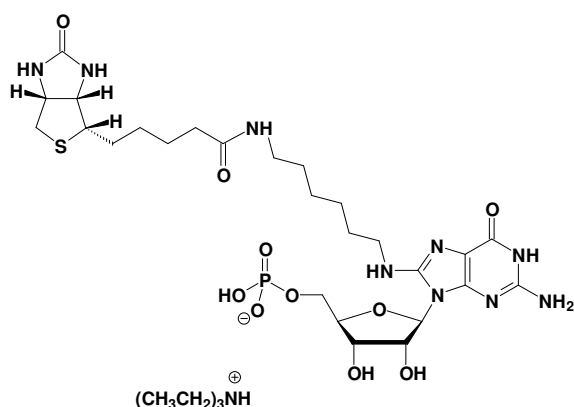


8-[(6-Amino)hexyl]-amino-GMP - Biotin

8-[(6-Amino)hexyl]-amino-guanosine-5'-monophosphate - Biotin, Triethylammonium salt

Cat. No.	Amount
NU-829-BIO	50 μ l / 5 mM



Cat. No.: NU-829-BIO

Molecular Formula: C₂₆H₄₁N₉O₁₀PS (Anion)

Molecular Weight: 702.69 (Anion)

Purity: > 95%, clear aqueous solution, pH 7.5

Storage conditions:

Short term exposure (up to 1 week cumulative) to ambient temperature possible. Long term storage at < -20°C. If stored as recommended, Jena Bioscience guarantees optimal performance of this product for 6 months after date of delivery.

For research use only!