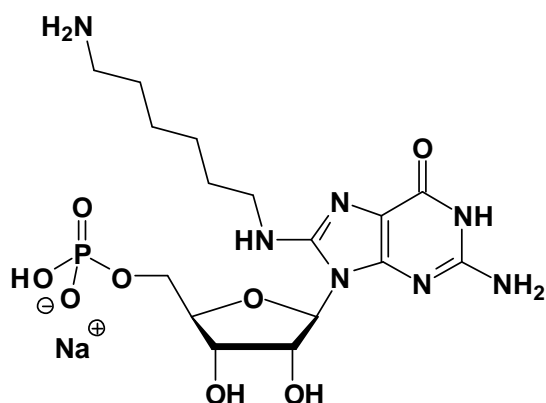


8-[(6-Amino)hexyl]-amino-GMP

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

Cat. No.	Amount
NU-829S	100 Units
NU-829L	500 Units



Cat. No.: NU-829

Molecular Formula: C₁₆H₂₇N₇O₈P (Anion)

Molecular Weight: 476.4 (Anion)

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

Spectroscopic properties: λ_{max} 256 nm ; ε 17000

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 12 months after date of delivery.

For research use only!

1 unit = 1 μl of a 10 mM solution