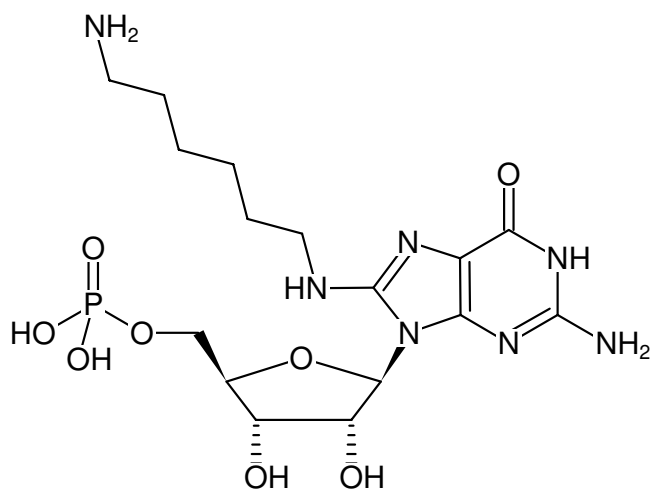




8-(6-Aminoheptyl)-amino-GMP

8-(6-Aminoheptyl)-amino-guanosine-5'-monophosphate, Sodium salt

Cat. No.	Amount
NU-829S	100 μ l (10 mM)
NU-829L	5 x 100 μ l (10 mM)



Structural formula of 8-(6-Aminoheptyl)-amino-GMP

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: C₁₆H₂₈N₇O₈P (free acid)

Molecular Weight: 477.41 g/mol (free acid)

Exact Mass: 477.17 g/mol (free acid)

Purity: \geq 95 % (HPLC)

Form: solution in water

Color: colorless to slightly yellow

Concentration: 10 mM - 11 mM

pH: 7.5 \pm 0.5

Spectroscopic Properties: λ_{max} 256 nm, ϵ 17.0 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)