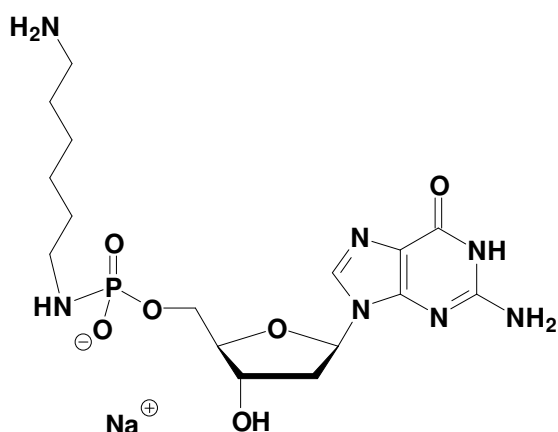


α -[(6-Aminohexyl)-imido]-dGMP

α -[(6-Aminohexyl)-imido]-2'-Deoxy-Guanosine-5'-monophosphate, Sodium salt

Cat. No.	Amount
NU-845S	5 mg
NU-845L	25 mg



Cat. No.: NU-845

Molecular Formula: C₁₆H₂₇N₇O₆P (Anion)

Molecular Weight: 444.40 (Anion)

Purity: > 95%, lyophilized powder

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!