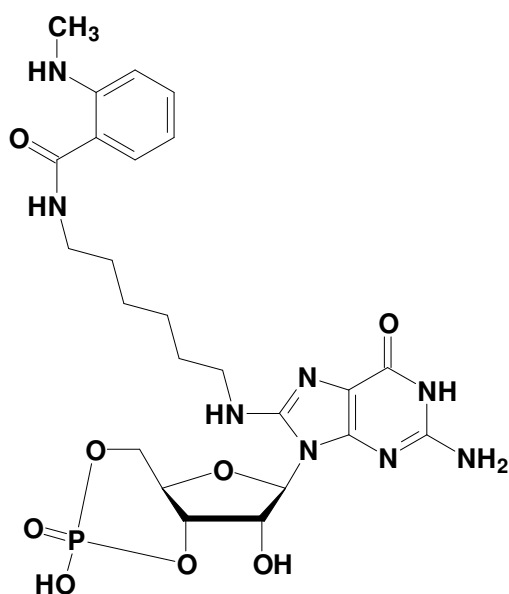


8-[(6-Amino)hexyl]-amino-cGMP - MANT

8-[(6-Amino)hexyl]-amino-guanosine-3',5'-cyclic monophosphate, labeled with MANT,
 Triethylammonium salt

Cat. No.	Amount
NU-832-MNT	200 μ l (5 mM)



Structural formula of 8-[(6-Amino)hexyl]-amino-cGMP - MANT

Cat. No.: NU-832-MNT

Molecular Formula: C₂₄H₃₃N₈O₈P (free acid)

Molecular Weight: 592.54 (free acid)

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

Spectroscopic properties:

λ_{exc} 335 nm; λ_{em} 440 nm; ϵ 2,700 cm⁻¹ M⁻¹

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!