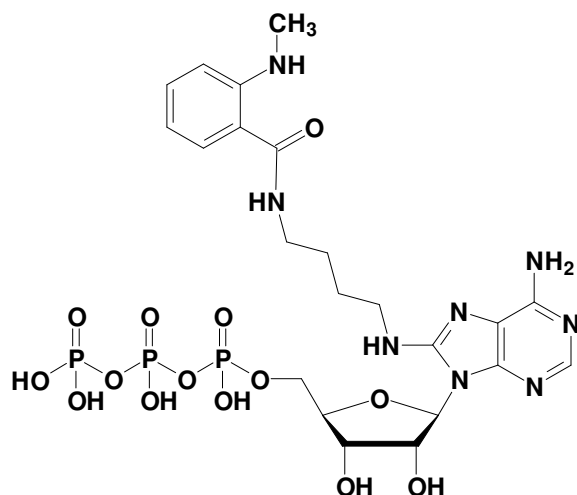


8-[(4-Amino)butyl]-amino-ATP-MNT (MABA-ATP) 8-[(4-Amino)butyl]-amino-adenosine-5'-triphosphate, labeled with MANT, Triethylammonium salt

Cat. No.	Amount
NU-806-MNT	50 μ l / 5 mM



structural formula of 8-[(4-Amino)butyl]-amino-adenosine-5'-triphosphate-MNT

Cat. No.: NU-806-MNT

Molecular Formula: C₂₂H₃₃N₈O₁₄P₃ (free acid)

Molecular Weight: 726.47 (free acid)

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

Spectroscopic properties:

λ_{exc} 335 nm; λ_{em} 440 nm; ϵ_{280} 18000 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!