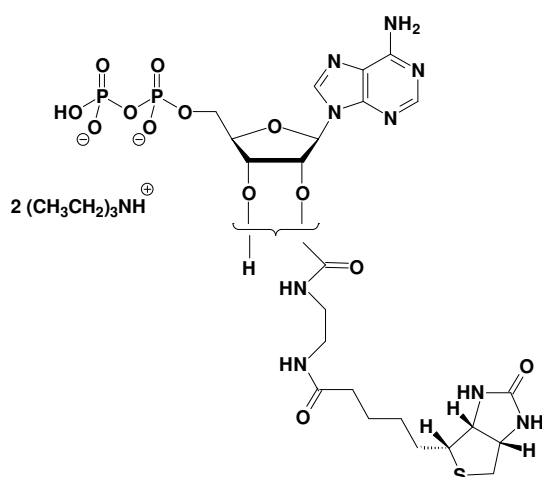


EDA-ADP-Biotin

2'/3'-O-(2-Aminoethyl-carbamoyl)-Adenosine-5'-diphosphate - Biotin

Cat. No.	Amount
NU-802-BIO	50 μ l / 5 mM



Cat. No.: NU-802-BIO

Molecular Formula: $C_{24}H_{37}N_9O_{13}P_2S$ (Anion)

Molecular Weight: 753.61 (Anion)

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

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!