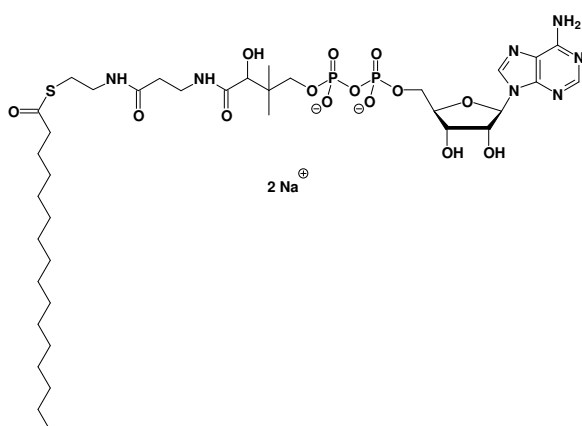


3'-Dephospho-Palmitoyl-CoA

3'-Dephospho-Palmitoyl-Coenzym A, Sodium salt

Cat. No.	Amount
NU-1180S	20 Units
NU-1180L	100 Units



Cat. No.: NU-1180

Molecular Formula: $C_{37}H_{63}N_7O_{14}P_2S$ (Anion)

Molecular Weight: 923.95 (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.

* 1 unit = 1 μ l of a 10 mM solution

For research use only!