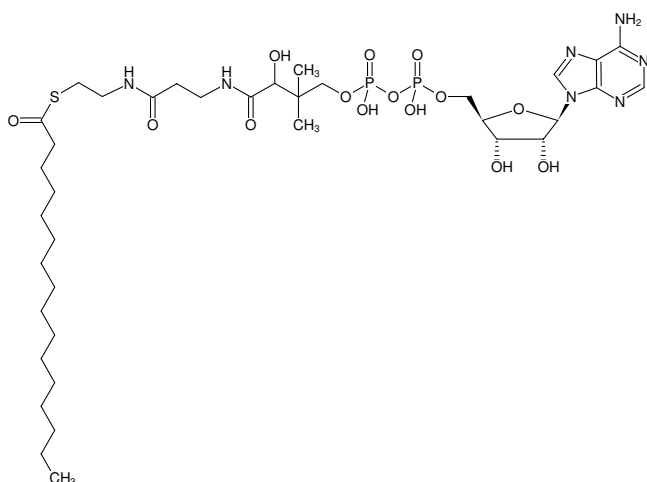




3'-Dephospho-Palmitoyl-CoA

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

Cat. No.	Amount
NU-1180S	20 µl (10 mM)
NU-1180L	5 x 20 µl (10 mM)



Structural formula of 3'-Dephospho-Palmitoyl-CoA

For general laboratory use.

Shipping: shipped on gel packs

Storage Conditions: store at -20 °C

Short term exposure (up to 1 week cumulative) to ambient temperature possible.

Shelf Life: 12 months after date of delivery

Molecular Formula: C₃₇H₆₅N₇O₁₄P₂S (free acid)

Molecular Weight: 925.96 g/mol (free acid)

Exact Mass: 925.38 g/mol (free acid)

Purity: ≥ 95 % (HPLC)

Form: solution in water

Color: colorless to slightly yellow

Concentration: 10 mM - 11 mM

pH: 7.5 ±0.5

Spectroscopic Properties: λ_{max} 258 nm, ε 10.5 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)