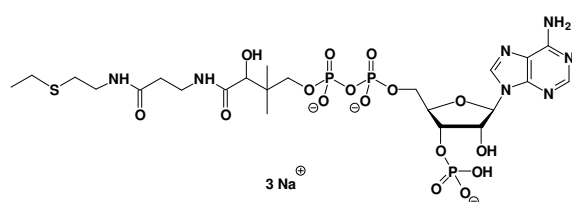


S-Ethyl-CoA

S-Ethyl-Coenzym A, Sodium salt

Cat. No.	Amount
NU-1168S	20 Units
NU-1168L	100 Units



Cat. No.: NU-1168

Molecular Formula: $C_{23}H_{37}N_7O_{16}P_3S$ (Anion)

Molecular Weight: 792.56 (Anion)

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

Spectroscopic properties:

λ_{max} 260 nm; ϵ 16000
(0,1M Tris-HCl 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!