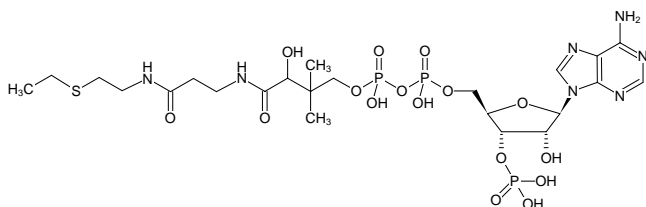


**S-Ethyl-CoA**

S-Ethyl-Coenzyme A, Sodium salt

Cat. No.	Amount
NU-1168S	20 µl (10 mM)
NU-1168L	5 x 20 µl (10 mM)



Structural formula of S-Ethyl-CoA

For research use only!**Shipping:** shipped on blue ice**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:** C₂₃H₄₀N₇O₁₆P₃S (free acid)**Molecular Weight:** 795.58 g/mol (free acid)**Purity:** ≥ 95 % (HPLC)**Form:** clear aqueous solution**Concentration:** 10 mM - 11 mM**pH:** 7.5 ± 0.5**Spectroscopic Properties:** λ_{max} 260 nm; ε 16.0 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)