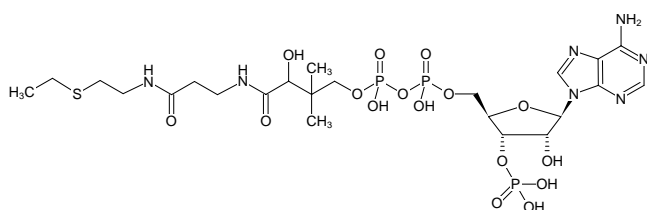


**S-Ethyl-CoA**

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:** C<sub>23</sub>H<sub>40</sub>N<sub>7</sub>O<sub>16</sub>P<sub>3</sub>S (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<sub>max</sub> 260 nm; ε 16.0 L mmol<sup>-1</sup> cm<sup>-1</sup> (Tris-HCl pH 7.5)