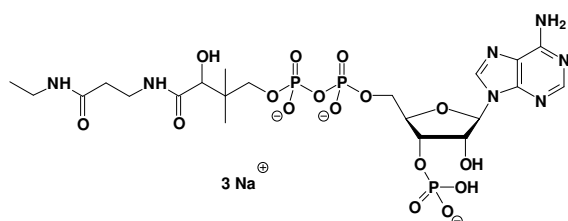


## Desulpho-CoA

Desulpho-Coenzym A, Sodium salt

Cat. No.	Amount
NU-1162S	20 Units
NU-1162L	100 Units



**Cat. No.:** NU-1162

**Molecular Formula:** C<sub>21</sub>H<sub>33</sub>N<sub>7</sub>O<sub>16</sub>P<sub>3</sub> (Anion)

**Molecular Weight:** 732.45 (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 µl of a 10 mM solution

**For research use only!**