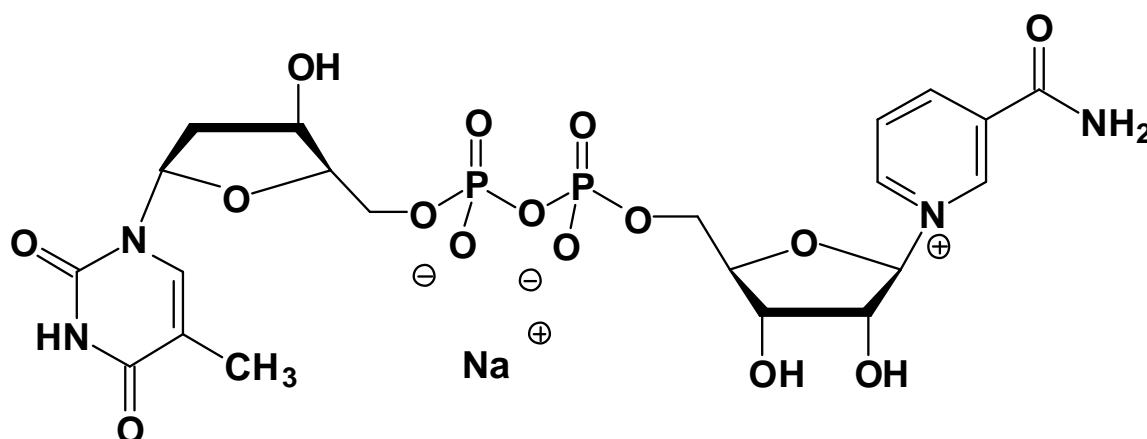


## TP<sub>2</sub>(Nic)

Nicotinamide thymine dinucleotide, Sodium salt

Cat. No.	Amount
NU-518	300 Units



**Cat. No.:** NU-518

**Molecular Formula:** C<sub>21</sub>H<sub>27</sub>N<sub>4</sub>O<sub>15</sub>P<sub>2</sub> (Anion)

**Molecular Weight:** 637.41 (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.

**For research use only!**

1 unit = 1 µl of a 10 mM solution

**Selected References:**

Fawcett *et al.* (1962) Preparation and properties of some nicotinamide adenine dinucleotide analogues with pentose and purine modifications. *J. of Biological Chem.* **237**(5): 1709