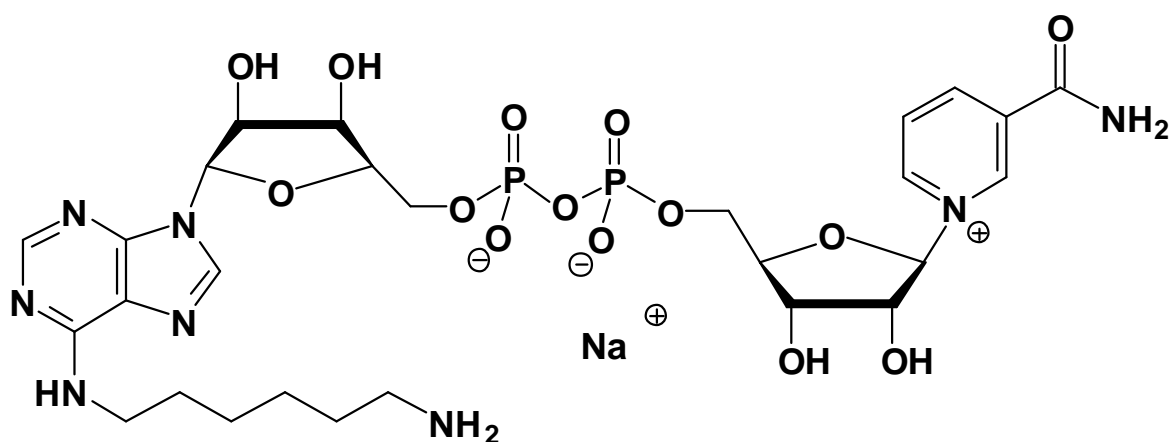


## N<sup>6</sup>-(6-Amino)hexyl-AP<sub>2</sub>(Nic)

Nicotinamide N<sup>6</sup>-(6-Amino)hexyl-adenine dinucleotide, Sodium salt

Cat. No.	Amount
NU-527	300 Units



**Cat. No.:** NU-527

**Molecular Formula:** C<sub>27</sub>H<sub>39</sub>N<sub>8</sub>O<sub>14</sub>P<sub>2</sub> (Anion)

**Molecular Weight:** 761.59 (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