

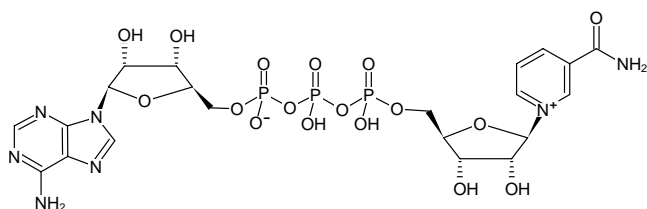
**AP<sub>3</sub>(Nic)**

P<sup>1</sup>-(5'-Adenosyl) P<sup>3</sup>-[5'-(1-β D-Ribofuranosyl-nicotinamide)] triphosphate, Sodium salt

Cat. No.	Amount
NU-514	300 µl (10 mM)

**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.



Structural formula of AP<sub>3</sub>(Nic)

**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>21</sub>H<sub>28</sub>N<sub>7</sub>O<sub>17</sub>P<sub>3</sub> (free acid)

**Molecular Weight:** 743.41 g/mol (free acid)

**Purity:** ≥ 95 % (HPLC)

**Form:** clear aqueous solution

**Concentration:** 10 mM - 11 mM

**pH:** 7.5 ±0.5