

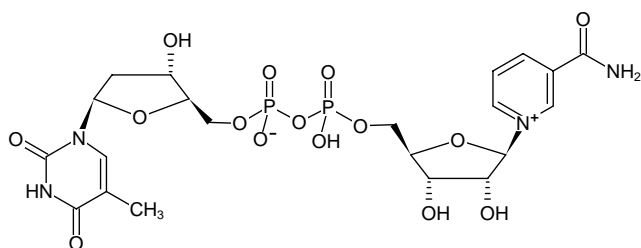
**TP₂(Nic)**

Nicotinamide thymine dinucleotide, Sodium salt

Cat. No.	Amount
NU-518	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 TP₂(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:** C₂₁H₂₈N₄O₁₅P₂ (free acid)**Molecular Weight:** 638.41 g/mol (free acid)**Purity:** ≥ 95 % (HPLC)**Form:** clear aqueous solution**Concentration:** 10 mM - 11 mM**pH:** 7.5 ±0.5