

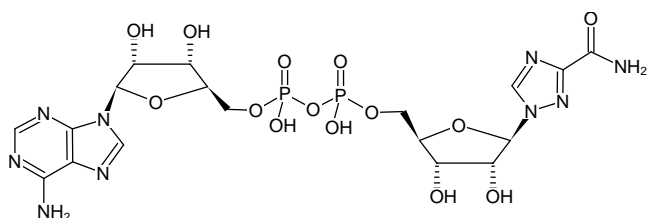
**AP₂Ribavirin**

1,2,4-Triazole-3-carboxamide adenine dinucleotide, Sodium salt

Cat. No.	Amount
NU-516	300 µl (10 mM)

Selected References:

Gebeyehu *et al.* (1985) Ribavirin, tiazofurin, and selenazofurin: mononucleotides and nicotinamide adenine dinucleotide analogues. Synthesis, structure and interactions with IMP dehydrogenase. *J. Med. Chem.* **28**: 99.

Structural formula of AP₂Ribavirin**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:** C₁₈H₂₅N₉O₁₄P₂ (free acid)**Molecular Weight:** 653.39 g/mol (free acid)**Purity:** ≥ 95 % (HPLC)**Form:** clear aqueous solution**Concentration:** 10 mM - 11 mM**pH:** 7.5 ±0.5