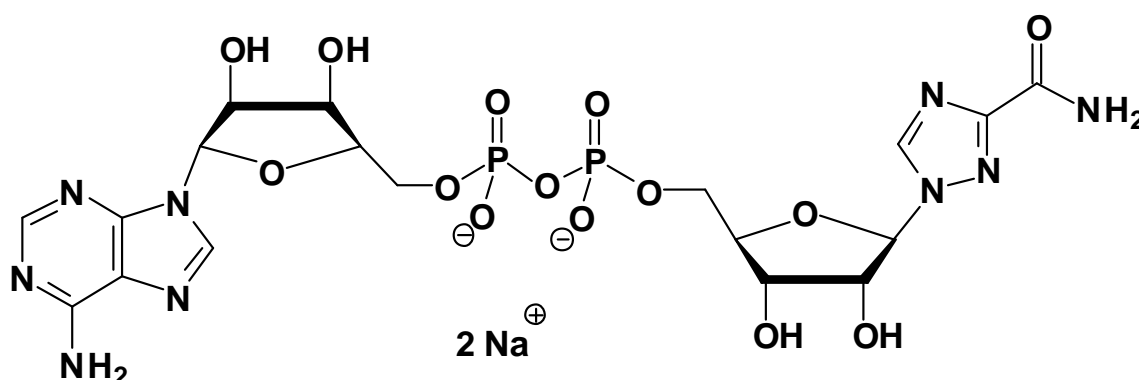


AP₂Ribavirin

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

Cat. No.	Amount
NU-516	300 Units



Cat. No.: NU-516

Molecular Formula: C₁₈H₂₃N₉O₁₄P₂ (Anion)

Molecular Weight: 651.38 (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:

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