

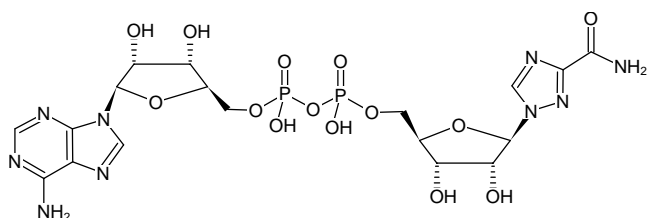
**AP<sub>2</sub>Ribavirin**

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 AP<sub>2</sub>Ribavirin**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. If stored as recommended, Jena Bioscience guarantees optimal performance of this product for 12 months after date of delivery.

**Shelf Life:** 12 months**Molecular Formula:** C<sub>18</sub>H<sub>25</sub>N<sub>9</sub>O<sub>14</sub>P<sub>2</sub> (free acid)**Molecular Weight:** 653.39 g/mol (free acid)**Purity:** ≥ 95 % (HPLC)**Form:** clear aqueous solution, pH 7.5 ±0.5**Concentration:** 10 mM - 11 mM**pH:** 7.5 ±0.5