



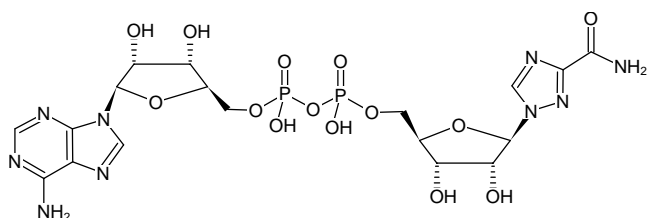
## AP<sub>2</sub>Ribavirin

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.

**Shelf Life:** 12 months after date of delivery

**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

**Concentration:** 10 mM - 11 mM

**pH:** 7.5 ± 0.5