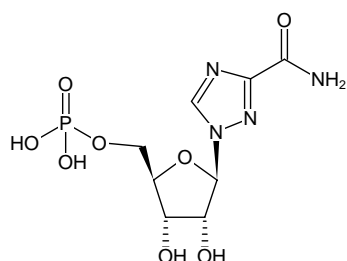




Ribavirin-5'-monophosphate

Sodium salt

Cat. No.	Amount
NU-021-2	2 mg



Structural formula of Ribavirin-5'-monophosphate

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: C₈H₁₃N₄O₈P (free acid)

Molecular Weight: 324.18 g/mol (free acid)

CAS#: 40925-28-8

Purity: ≥ 95 % (HPLC)

Form: white powder

Spectroscopic Properties: λ_{max} 220 nm; ε 7.7 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)

Selected References:

Melendez *et al.* (2009) HPLC-MS/MS method for the intracellular determination of ribavirin monophosphate and ribavirin triphosphate in CEMss cells. *Journal of Pharmaceutical and Biomedical Analysis* **49** (5):1233.

Prosize *et al.* (2002) Crystal Structure of *Tritrichomonas foetus* Inosine Monophosphate Dehydrogenase in Complex with the Inhibitor Ribavirin Monophosphate Reveals a Catalysis-dependent Ion-binding Site. *Journal of Biological Chemistry* **277** (52):50654.