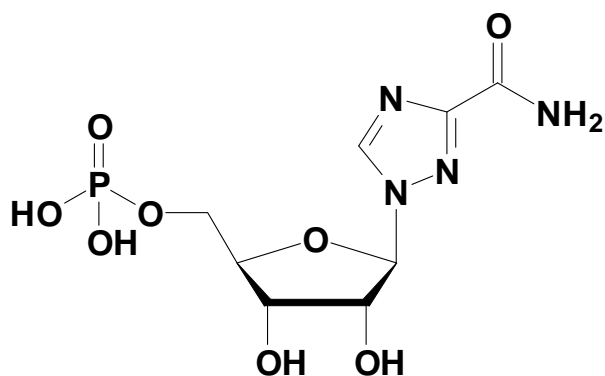


Ribavirin-5'-monophosphate Sodium salt

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



CAS number: 40925-28-8

Molecular Formula: C₈H₁₃N₄O₈P (free acid)

Molecular Weight: 324.18 g/mol (free acid)

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.

Purity: >95% (HPLC)

Appearance: white solid

Spectroscopic properties:

λ_{abs} 220 nm; ϵ 7700 cm⁻¹M⁻¹ (Tris-HCl, pH 7.5)

For research use only!

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 Trichomonas 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.