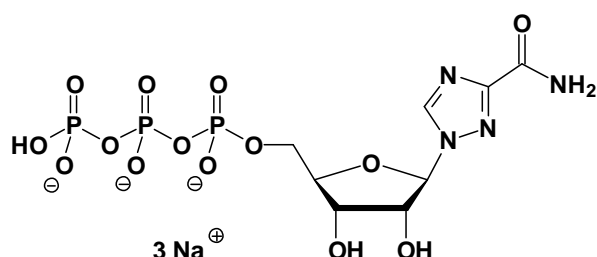


Ribavirin-5'-triphosphate

1-β-D-Ribofuranosyl-1,2,4-triazole-3-carboxamide-5'-triphosphate, Sodium salt

| Cat. No. | Amount |
|----------|-----------|
| NU-1105S | 20 Units |
| NU-1105L | 100 Units |



Cat. No.: NU-1105

Molecular Formula: C₈H₁₂N₄O₁₄P₃ (Anion)

Molecular Weight: 481.12 (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:

Maag *et al.* (2001) Hepatitis C virus RNA-dependent RNA polymerase (NS5B) as a mediator of the antiviral activity of ribavirin. *J. Biol. Chem.* **276** (49):46094.

Lanford *et al.* (2001) Ribavirin induces error-prone replication of GB virus B in primary tamarin hepatocytes. *J. Virol.* **75** (17):8074.

Crotty *et al.* (2000) The broad-spectrum antiviral ribonucleoside ribavirin is an RNA virus mutagen. *Nat. Med.* **6** (12):1375.

Ulrich *et al.* (2000) Towards the engineering of an orthogonal protein kinase/nucleotide triphosphate pair. *Tetrahedron* **56** (48):9495.

Hong *et al.* (1999) Direct antiviral activity of ribavirin: Hepatitis C virus NS5B polymerase incorporates ribavirin triphosphate into nascent RNA products. *Hepatology* **30** (4):773 Part 2 Suppl. S.

Jarvis *et al.* (1998) Ribavirin uptake by human erythrocytes and the involvement of nitrobenzylthioinosine-sensitive (es)- nucleoside transporters. *Brit. J. Pharmacol.* **123** (8):1587.

Hosoya *et al.* (1993) Comparative inhibitory effects of various nucleoside and nonnucleoside analogs on replication of influenza-virus type-A and type-B in-vitro and in-vivo. *J. Infect. Dis.* **168** (3):641.

Wray *et al.* (1985) Effect of ribavirin triphosphate on primer generation and elongation during influenza-virus transcription in-vitro. *Antivir. Res.* **5** (1):39.

Eriksson *et al.* (1977) Inhibition of influenza-virus ribonucleic-acid polymerase by ribavirin triphosphate. *Antimicrob. Agents Ch.* **11** (6):946.