

Ribavirin

Broad-spectrum antiviral nucleoside analog

Ribavirin is a synthetic guanosine analog with broad-spectrum antiviral activity. It is metabolized intracellularly to its active triphosphate form, which inhibits viral RNA-dependent RNA polymerase and induces lethal mutagenesis in viral genomes. Ribavirin also inhibits inosine monophosphate dehydrogenase (IMPDH), reducing GTP synthesis and viral replication. It has demonstrated EC_{50} values ranging from 2-95 $\mu\text{g/mL}$ across various RNA and DNA viruses and 109.5 μM against SARS-CoV-2 in Vero cells.

Key features and applications include:

- **Broad Antiviral Spectrum:** Active against both RNA and DNA viruses, including RSV, HCV, HIV, influenza, and coronaviruses.
- **Multiple Mechanisms of Action:** Inhibits viral polymerases, depletes GTP pools via IMPDH inhibition, and induces error catastrophe in viral genomes.
- **Versatile Administration:** Used in oral, intravenous, and aerosolized forms in clinical and research settings.
- **Combination Therapy:** Commonly used with interferon- α or direct-acting antivirals (DAAs) in hepatitis C treatment.

Research Applications:

- Antiviral drug screening
- Viral replication and mutagenesis studies
- Host-virus interaction models
- Immunomodulatory and anti-inflammatory research

Relevant disease states include:

- **Hepatitis C Virus (HCV):** Historically used in combination therapy to enhance viral clearance.
- **Respiratory Syncytial Virus (RSV):** Administered via aerosol in severe pediatric infections.

- **Viral Hemorrhagic Fevers:** Investigated for use in Lassa fever and other high-mortality viral infections.
- **COVID-19:** Studied for its potential to inhibit SARS-CoV-2 replication.
- **Cancer Research:** Shown to inhibit tumor growth and metastasis in soft tissue sarcomas by downregulating PRMT1 and PRMT5.

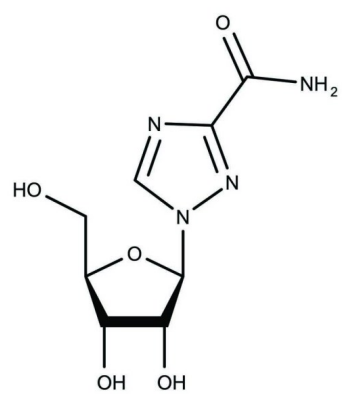
Ordering Information

Order Online »

ENZ-CHM342-0250	250mg
-----------------	-------

Manuals, SDS & CofA

View Online »



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended.
Short Term Storage	-20°C
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name	Ribasphere, 1-[(2R,3R,4S,5R)-3,4-dihydroxy-5-(hydroxymethyl)oxolan-2-yl]-1,2,4-triazole-3-carboxamide
Appearance	White solid.
CAS	36791-04-5
Formula	$C_8H_{12}N_4O_5$
Identity	Determined by NMR.
MW	244.21
Purity	≥98% (TLC)
Solubility	Soluble in DMSO (up to 25 mg/mL) or in water (up to 25 mg/mL).



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
[info-
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
[info-
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)