

Potent competitive inhibitor of inosine monophosphate (IMP) dehydrogenase, viral RNA polymerase and viral mRNA guanylyltransferase, causes mutagenesis and reduction of viral infectivity.

*BLG-R001-25 (5 x 5 μ mol pack size) is not sold in the U.S. or Canada.
Please [contact us](#) for available options.*

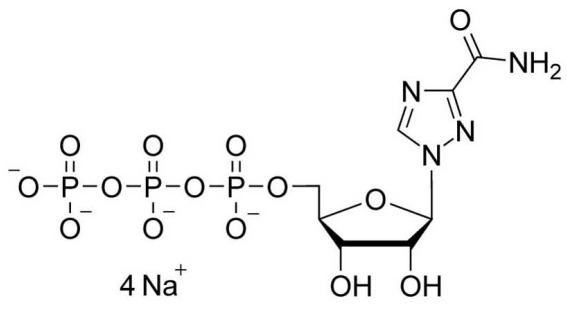
Ordering Information

[Order Online »](#)

| | |
|--------------------|---------------|
| BLG-R001-05 | 5 μ mol |
| BLG-R001-25 | 5x5 μ mol |

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Long Term Storage -80°C

Shipping Dry Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name 1-β-D-Ribofuranosyl-1, 2, 4-triazole-3-carboxamide-5'-O-triphosphate . tetrasodium salt, Ribavirin-5'-O-triphosphate

Appearance 10 mM aqueous solution

CAS 63142-71-2

Couple Type Modified nucleotides

Formula $C_8H_{15}N_4O_{14}P_3$ (free acid)

MW 484.2

Purity > 95% HPLC



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)