

5'-N₃-Guo

Reactive guanosine analogue

5'-N₃-Guo is an analogue of guanosine, modified in its 5'-position by a reactive azido group. The compound is offered as component of the Ko-Libri library of rare nucleosides and nucleobases (Cat. No. K 001) as well.

*BLG-A079-25 (5 x 5 µmol pack size) is not sold in the U.S. or Canada.
Please [contact us](#) for available options.*

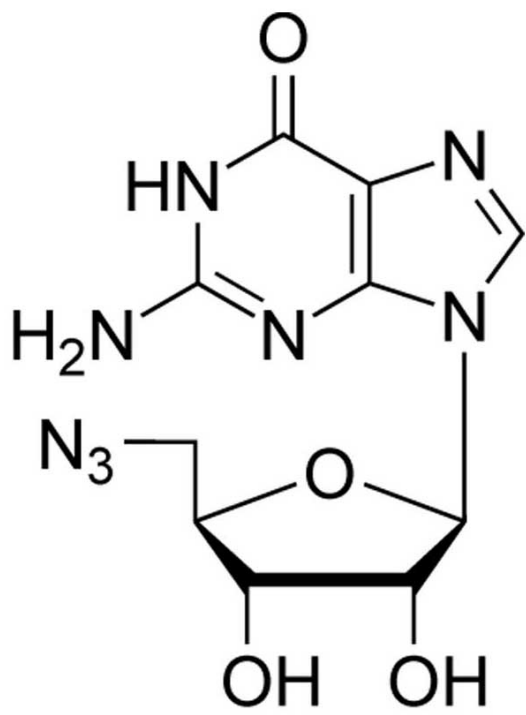
Ordering Information

[Order Online »](#)

BLG-A079-05	5µmol
BLG-A079-25	5x5µmol

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Long Term Storage -20°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name 5'-Azido-5'-deoxyguanosine

CAS 42204-44-4

Formula $C_{10}H_{12}N_8O_4$

MW 308.3

Purity > 95% HPLC

Quantity 5µmol are equivalent to ~1.5 mg



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

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

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

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

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