

3'-N₃-Ado

Reactive adenosine analogue

3'-N₃-Ado is an analogue of adenosine where the 3'-hydroxy group has been replaced by a reactive azido group which allows for reaction with suitable ligands.

BLG-A243-25 (5 x 5 µmol pack size) is not sold in the U.S. or Canada. Please [contact us](#) for available options.

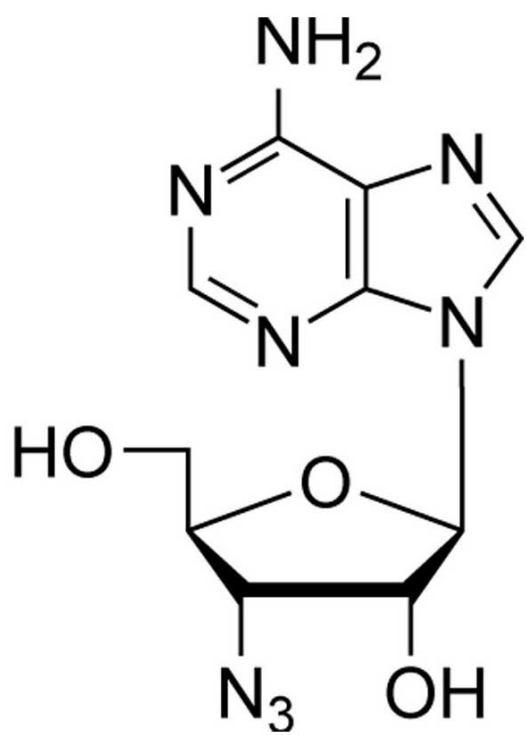
Ordering Information

[Order Online »](#)

BLG-A243-05	5µmol
BLG-A243-25	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 3'-Azido-3'-deoxyadenosine

CAS 58699-62-0

Formula $C_{10}H_{12}N_8O_3$

MW 292.3

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