

2',5'-

Dideoxyadenosine

Adenylate cyclase inhibitor

A useful, cell-permeable inhibitor of adenylate cyclase ($ID_{50}=250\mu M$).
Inhibits growth and causes cell differentiation ($1.5\mu M$) in K562 cells.

Citations: 6

[View Online »](#)

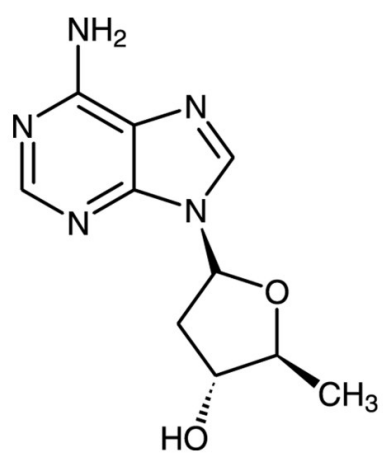
Ordering Information

[Order Online »](#)

BML-CN110-0005	5mg
BML-CN110-0025	25mg

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Solutions are stable up to 3 months when stored at -20°C
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Appearance	White to off-white solid.
CAS	6698-26-6
Couple Target	Adenylate cyclase
Couple Type	Inhibitor
Formula	$C_{10}H_{13}N_5O_2$
MW	235.2
Purity	≥98% (HPLC)
Solubility	Soluble in water (20mg/ml at 50°C) or DMSO (30mg/ml).

Last modified: May 29, 2024



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)