

Puromycin . 2HCl

Protein synthesis inhibitor

Inhibits protein synthesis. Inhibits translation by causing premature release of nascent polypeptide. Induces apoptosis.

Citations: 34

[View Online »](#)

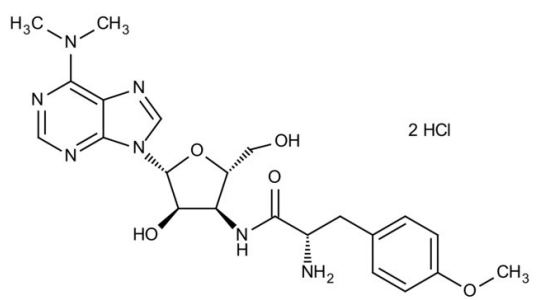
Ordering Information

[Order Online »](#)

BML-GR312-0050	50mg
BML-GR312-0250	250mg

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Stable for at least 1 year after receipt when stored, as supplied, at -20°C. Stock solutions are stable for up to 3 months at -20°C.
Handling	Protect from moisture.
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	3'-[[2-Amino-3-(4-methoxyphenyl)-1-oxopropyl]amino]-3'-deoxy-N,N-dimethyl-adenosine . 2HCl, 3'-(α -Amino-p-methoxy-hydrocinnamamido-3'-deoxy-N,N-dimethyl-adenosine . 2HCl
Appearance	White to off-white solid.
CAS	58-58-2
Formula	$C_{22}H_{29}N_7O_5 \cdot 2HCl$
MI	14: 7943
MW	544.4
Purity	$\geq 98\%$ (HPLC, TLC)
RTECS	AU7355000
Solubility	Soluble in water (50mg/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-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

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

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

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