

Pyrrolidinedithiocarbamic acid . ammonium salt

Inhibitor of NF- κ B activation

Inhibitor of NF- κ B activation (induced by either PMA or TNF- α) in various cell types.

Citations: 14

[View Online »](#)

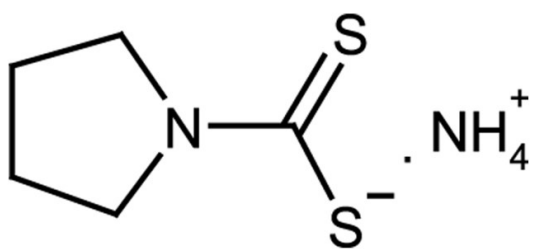
Ordering Information

[Order Online »](#)

ALX-400-002-G005	5g
------------------	----

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability As indicated on product label or CoA when stored as recommended.

Long Term Storage +4°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Ammoniumpyrrolidinedithiocarbamate,
Ammoniumtetramethylenedithiocarbamate

Appearance White to-off white solid.

CAS 5108-96-3

Couple Target NF-kappaB

Couple Type Inhibitor

Formula $C_5H_8NS_2 \cdot NH_4$

MW 146.3 . 18.0

Purity ≥98% (HPLC)

Solubility Soluble in water.

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-
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)