

Caffeine

CNS stimulant

CNS stimulant. Blocks adenosine A1 and A2A receptors. cAMP phosphodiesterase inhibitor. Interferes with the uptake and storage of Ca^{2+} by the sarcoplasmic reticulum in skeletal muscle. Prevents apoptosis and cell cycle effects induced by various chemicals. Inhibits cellular DNA repair mechanisms. Anti-inflammatory.

Citations: 15

[View Online »](#)

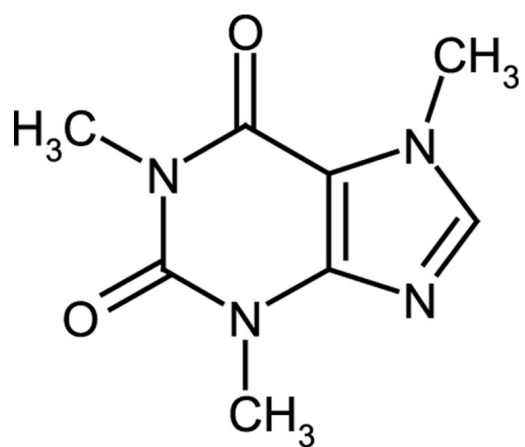
Ordering Information

[Order Online »](#)

ALX-550-322-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 Ambient

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name 1,3,7-Trimethylxanthine, 3,7-Dihydro-1,3,7-trimethyl-1H-purine-2,6-dione

Appearance White to colorless powder.

CAS 58-08-2

Couple Target Adenosine receptor, PDE

Couple Type Blocker, Inhibitor

Formula $C_8H_{10}N_4O_2$

Identity Determined by IR.

MI 14: 1636

MW 194.2

Purity ≥99% (HPLC)

Solubility Soluble in chloroform, 100% alcohol or hot water.

Source Found in tea leaves, coffee beans, cocoa beans, maté leaves, guarana paste and kola nuts.



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)