

# Caffeine

CNS stimulant

CNS stimulant. Blocks adenosine A1 and A2A receptors. cAMP phosphodiesterase inhibitor. Interferes with the uptake and storage of  $\text{Ca}^{2+}$  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 »](#)

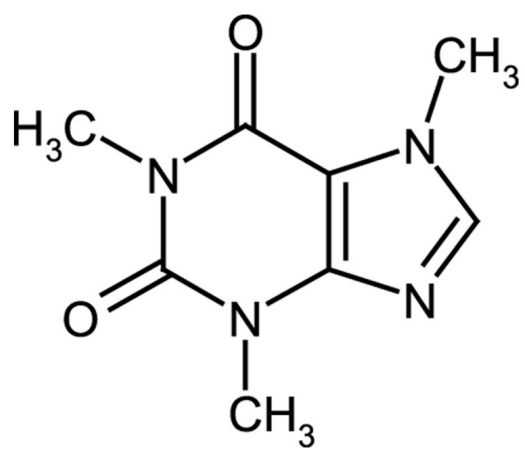
## Ordering Information

[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.

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-be@enzolifesciences.com)

Belgium, The Netherlands  
& Luxembourg  
Phone: +32 3 466 0420  
[info-  
be@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)