

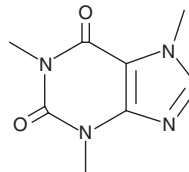
PRODUCT INFORMATION



Caffeine

Item No. 22975

CAS Registry No.: 58-08-2
Formal Name: 3,7-dihydro-1,3,7-trimethyl-1H-purine-2,6-dione
Synonyms: Methyltheobromine, NSC 5036
MF: C₈H₁₀N₄O₂
FW: 194.2
Purity: ≥98%
Supplied as: A neat solid
Storage: -20°C
Stability: ≥4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Description

Caffeine (Item No. 22975) is an analytical reference material categorized as a stimulant.¹ It is found in kratom extract that has been used to overcome opioid withdrawal.² This product is intended for analytical forensic applications. This product is also available as a general research tool (Item No. 14118).

This product is qualified as a Reference Material that has been manufactured and tested to ISO/IEC 17025 and ISO 17034 international standards.

References

1. Judelson, D.A., Preston, A.G., Miller, D.L., *et al.* Effects of theobromine and caffeine on mood and vigilance. *J. Clin. Psychopharmacol.* **33**(4), 499-506 (2013).
2. Chittrakarn, S., Penjamras, P., and Keawpradub, N. Quantitative analysis of mitragynine, codeine, caffeine, chlorpheniramine and phenylephrine in a kratom (*Mitragyna speciosa* Korth.) cocktail using high-performance liquid chromatography. *Forensic Sci. Int.* **217**(1-3), 81-83 (2011).

WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the [complete](#) Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 04/15/2019

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM