

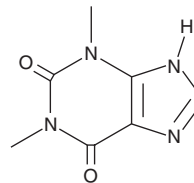
PRODUCT INFORMATION



Theophylline

Item No. 32993

CAS Registry No.: 58-55-9
Formal Name: 3,9-dihydro-1,3-dimethyl-1H-purine-2,6-dione
Synonyms: 1,3-Dimethylxanthine, NSC 2066
MF: C₇H₈N₄O₂
FW: 180.2
Purity: ≥98%
UV/Vis.: λ_{max}: 271 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥5 years



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

Description

Theophylline (Item No. 32993) is an analytical reference standard categorized as a caffeine metabolite.¹ Theophylline has been found as an adulterant in seized heroin samples.² It has also been used as a marker of authentic human urine specimens.¹ This product is intended for use in analytical forensic applications. This product is also available as a general research tool (Item No. 23760).

References

1. Goggin, M.M., Tann, C.-M., Miller, A., *et al.* Catching fakes: New markers of urine sample validity and invalidity. *J. Anal. Toxicol.* **41(2)**, 121-126 (2016).
2. Furst, R.T. The re-engineering of heroin: An emerging heroin “cutting” trend in New York City. *Addict. Res. Theory* **8(4)**, 357-379 (2000).

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, 06/22/2023

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