

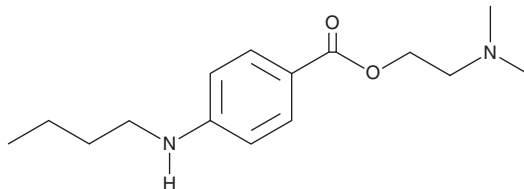
PRODUCT INFORMATION



Tetracaine

Item No. 30059

CAS Registry No.: 94-24-6
Formal Name: 4-(butylamino)-benzoic acid,
2-(dimethylamino)ethyl ester
Synonym: Tetracain
MF: C₁₅H₂₄N₂O₂
FW: 264.4
Purity: ≥98%
UV/Vis.: λ_{max}: 226, 309 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

Tetracaine (Item No. 30059) is an analytical reference standard categorized as an anesthetic.¹ Tetracaine has been found as an adulterant in products marketed as male sexual function enhancers.² This product is intended for research and forensic applications.

References

1. Ireland, P.E., Ferguson, J.K.W., and Stark, E.J. The clinical and experimental comparison of cocaine and pontocaine as topical anesthetics in otolaryngological practice. *Trans. Am. Laryngol. Rhinol. Otol. Soc.* **57(55th Meeting)**, 438-450 (1951).
2. Lee, J.H., Cho, S.-H., Kim, J.Y., *et al.* Determination and quantification of nine adulterant local anaesthetics in illegal treatments for male premature ejaculation by GC-FID and GC-MS. *Int. J. Pharm. Pharm. Sci.* **8(3)**, 135-140 (2016).

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, 07/28/2023

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