

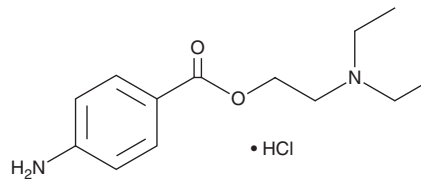
PRODUCT INFORMATION



Procaine (hydrochloride)

Item No. 20093

CAS Registry No.: 51-05-8
Formal Name: 4-amino-benzoic acid, 2-(diethylamino) ethyl ester, monohydrochloride
Synonym: Novocaine
MF: C₁₃H₂₀N₂O₂ • HCl
FW: 272.8
Purity: ≥98%
UV/Vis.: λ_{max}: 225, 299, 484 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

Procaine (hydrochloride) (Item No. 20093) is an analytical reference standard categorized as a local anesthetic that is used as an adulterant. Procaine blocks sodium channels, and formulations containing procaine are used to prevent or relieve pain for short periods, such as during dental procedures.¹ It is commonly found as an adulterant in illicit drugs.² This product is intended for research and forensic applications.

References

1. Epstein, S. and Chilton, N.W. The clinical effectiveness of certain local anesthetics in dental procedures. *Oral Surg. Oral Med. Oral Pathol.* **12(1)**, 93-103 (1959).
2. Cole, C., Jones, L., McVeigh, J., *et al.* CUT: A guide to adulterants, bulking agents and other contaminants found in illicit drugs. Liverpool John Moores University (2010).

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, 06/20/2023

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