

PRODUCT INFORMATION



Articaine (hydrochloride)

Item No. 44973

CAS Registry No.: 23964-57-0

Formal Name: 4-methyl-3-[[1-oxo-2-(propylamino) propyl]amino]-2-thiophenecarboxylic acid, methyl ester, monohydrochloride

Synonyms: Carticaine, HOE 045, HOE 40045

MF: $C_{13}H_{20}N_2O_3S \cdot HCl$

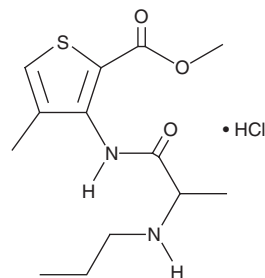
FW: 320.8

Purity: $\geq 98\%$

Supplied as: A solid

Storage: $-20^{\circ}C$

Stability: ≥ 4 years



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

Description

Articaine (hydrochloride) (Item No. 44973) is an analytical reference standard categorized as an anesthetic.¹ Articaine has been found as an adulterant in cocaine.² This product is intended for analytical forensic applications. This product is also available as a general research tool (Item No. 17883).

References

1. Weaver, J.M. Articaine, a new local anesthetic for American dentists: Will it supercede lidocaine? *Anesth. Prog.* **46**(4), 111-112 (1999).
2. Pawlik, E., Mahler, H., Hartung, B., et al. Drug-related death: Adulterants from cocaine preparations in lung tissue and blood. *Forensic Sci. Int.* **249**, 294-303 (2015).

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, 09/01/2026

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