

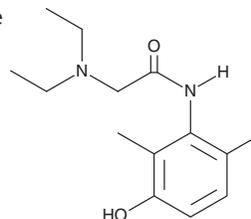
PRODUCT INFORMATION



3'-hydroxy Lidocaine

Item No. 22899

CAS Registry No.: 34604-55-2
Formal Name: 2-(diethylamino)-N-(3-hydroxy-2,6-dimethylphenyl)-acetamide
Synonyms: *m*-hydroxy Lidocaine, *meta*-hydroxy Lidocaine
MF: C₁₄H₂₂N₂O₂
FW: 250.3
Purity: ≥98%
UV/Vis.: λ_{max}: 282 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥4 years



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

Laboratory Procedures

3'-hydroxy Lidocaine is supplied as a crystalline solid. A stock solution may be made by dissolving the 3'-hydroxy lidocaine in the solvent of choice, which should be purged with an inert gas. 3'-hydroxy Lidocaine is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of 3'-hydroxy lidocaine in these solvents is approximately 30 mg/ml.

3'-hydroxy Lidocaine is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, 3'-hydroxy lidocaine should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. 3'-hydroxy Lidocaine has a solubility of approximately 0.11 mg/ml in a 1:8 solution of ethanol:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

3'-hydroxy Lidocaine is an active metabolite of lidocaine formed by the cytochrome P450 (CYP) isoforms CYP1A2 and CYP3A4.^{1,2}

References

1. Wang, J.-S., Backman, J.T., Taavitsainen, P., *et al.* Involvement of CYP1A2 and CYP3A4 in lidocaine N-deethylation and 3-hydroxylation in humans. *Drug Metab. Dispos.* **28(8)**, 959-965 (2000).
2. Hermansson, J., Glaumann, H., Karlén, B., *et al.* Metabolism of lidocaine in human liver *in vitro*. *Acta. Pharmacol. Toxicol. (Copenh)* **47(1)**, 49-52 (1980).

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, 12/02/2022

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