

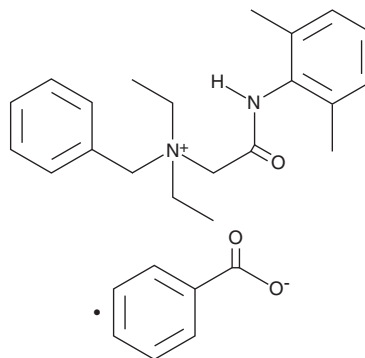
PRODUCT INFORMATION



Denatonium (benzoate salt)

Item No. 26350

CAS Registry No.: 3734-33-6
Formal Name: N-[2-[(2,6-dimethylphenyl)amino]-2-oxoethyl]-N,N-diethylbenzenemethanaminium, monobenzoate
Synonyms: Lidocaine benzyl benzoate, NSC 157658, WIN 16,568
MF: $C_{21}H_{29}N_2O \cdot C_7H_5O_2$
FW: 446.6
Purity: $\geq 98\%$
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

Denatonium (benzoate salt) (Item No. 26350) is an analytical reference standard categorized as a bittering agent.¹ Formulations containing denatonium have been used as bittering agents to reduce ingestion of toxic substances by children. This product is intended for research and forensic applications.

Reference

1. Hansen, S.R., Janssen, C., and Beasley, V.R. Denatonium benzoate as a deterrent to ingestion of toxic substances: Toxicity and efficacy. *Vet. Hum. Toxicol.* **35(3)**, 234-236 (1993).

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, 07/07/2023

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