

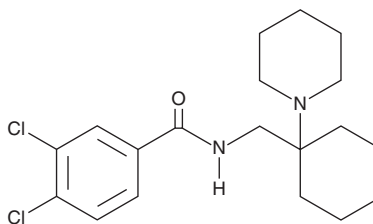
PRODUCT INFORMATION



AH 7959

Item No. 19366

CAS Registry No.: 763023-14-9
Formal Name: 3,4-dichloro-N-[[1-(1-piperidiny) cyclohexyl]methyl]-benzamide
MF: C₁₉H₂₆Cl₂N₂O
FW: 369.3
Purity: ≥98%
UV/Vis.: λ_{max}: 239 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

AH 7959 (Item No. 19366) is an analytical reference standard that is structurally categorized as an opioid. Unlike the related compound AH 7921 (Item Nos. 12036 | 19732), AH 7959 is ineffective as an analgesic in mice, by either the phenylquinone or hot plate test (ED_{50s} > 100 mg/kg for both tests).¹ This product is intended for research and forensic applications.

Reference

1. Brittain, R. T., Kellett, D. N., Neat, M. L., *et al.* Proceedings: Anti-nociceptive effects in N-substituted cyclohexylmethylbenzamides. *Br. J. Pharmacol.* **49(1)**, 158P-159P (1973).

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, 06/19/2023

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