PRODUCT INFORMATION



Bromadol

Item No. 15608

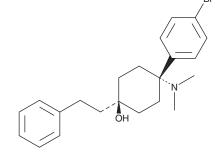
CAS Registry No.: 77239-98-6

Formal Name: trans-4-(4-bromophenyl)-

> 4-(dimethylamino)-1-(2phenylethyl)-cyclohexanol

Synonyms: trans-BDPC, trans-Bromadol

 $C_{22}H_{28}BrNO$ MF: FW: 402.4 **Purity:** ≥98% Supplied as: A neat solid Storage: -20°C Stability: ≥5 years



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

Description

Bromadol (Item No. 15608) is an analytical reference material categorized as an opioid. This product is intended for research and forensic applications.

This product is qualified as a Reference Material that has been manufactured and tested to ISO/IEC 17025 and ISO 17034 international standards for reference materials.

Reference

1. Von Voigtlander, P.F., Lednicer, D., Lewis, R.A., et al. 4-Aryl-4-aminocyclohexanone deivatives: A chemically novel series of analgesics including opioid antagonists and extremely potent agonists. Endogenous and exogenous opiate agonists and antagonists Pergamon Press, 17-21 (1979).

WARNING
THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

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

subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information Buyer agrees to purchase the mater can be found on our website.

Copyright Cayman Chemical Company, 08/26/2019

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