

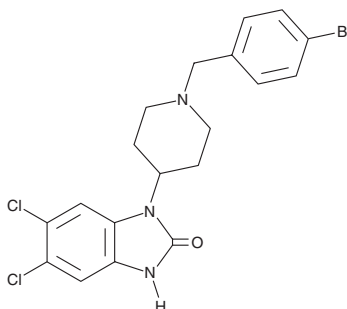
PRODUCT INFORMATION



5,6-dichloro Desmethylbromphine

Item No. 46168

CAS Registry No.: 2134602-44-9
Formal Name: 1-[1-[(4-bromophenyl)methyl]-4-piperidinyl]-5,6-dichloro-1,3-dihydro-2H-benzimidazol-2-one
Synonym: SR 15099
MF: C₁₉H₁₈BrCl₂N₃O
FW: 455.2
Purity: ≥98%
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years



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

Description

5,6-dichloro Desmethylbromphine (Item No. 46168) is an analytical reference standard categorized as an opioid.¹ 5,6-dichloro Desmethylbromphine increases latency to withdrawal in the hot plate and warm water tail-flick withdrawal assays and induces respiratory depression to a lesser extent than morphine (Item Nos. ISO60147 | 15464 | 25955) and fentanyl (Item Nos. ISO60197 | 14719 | 23580). This product is intended for research and forensic applications.

Reference

1. Schmid, C.L., Kennedy, N.M., Ross, N.C., *et al.* Bias factor and therapeutic window correlate to predict safer opioid analgesics. *Cell* **171**(5), 1165-1175 (2017).

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, 07/07/2026

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