

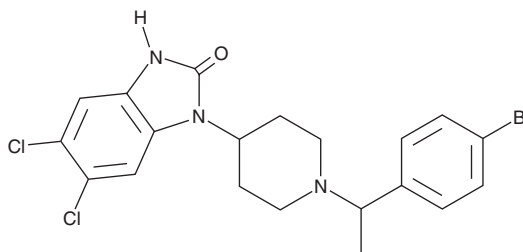
# PRODUCT INFORMATION



## 5,6-dichloro Brorphine

Item No. 40532

**CAS Registry No.:** 2133455-40-8  
**Formal Name:** 1-[1-[1-(4-bromophenyl)ethyl]-4-piperidinyl]-5,6-dichloro-1,3-dihydro-2H-benzimidazol-2-one  
**Synonym:** SR-14968  
**MF:** C<sub>20</sub>H<sub>20</sub>BrCl<sub>2</sub>N<sub>3</sub>O  
**FW:** 469.2  
**Purity:** ≥98%  
**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.

### Description

5,6-dichloro Brorphine (Item No. 40532) is an analytical reference standard categorized as an opioid.<sup>1</sup> 5,6-dichloro Brorphine induces antinociception, antinociceptive tolerance, naloxone-precipitated withdrawal symptoms, and conditioned place preference in mice.<sup>2</sup> This product is intended for research and forensic applications.

### References

1. Stahl, E.L., Schmid, C.L., Acevedo-Canabal, A., *et al.* G protein signaling-biased mu opioid receptor agonists that produce sustained G protein activation are noncompetitive agonists. *Proc. Natl. Acad. Sci. USA* **118(48)**, e2102178118 (2021).
2. Kudla, L., Bugno, R., Podlewska, S., *et al.* Comparison of an addictive potential of  $\mu$ -opioid receptor agonists with G protein bias: Behavioral and molecular modeling studies. *Pharmaceutics* **14(1)**, 55 (2021).

#### 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, 09/10/2025

#### 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