

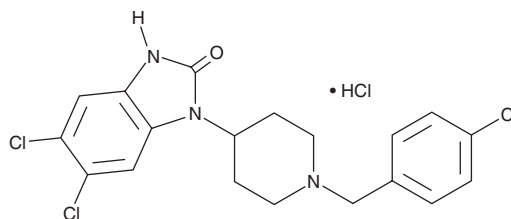
PRODUCT INFORMATION



SR-17018 (hydrochloride)

Item No. 45831

Formal Name: 5,6-dichloro-1-[1-[(4-chlorophenyl)methyl]-4-piperidinyl]-1,3-dihydro-2H-benzimidazol-2-one, monohydrochloride
Synonym: 5,6-dichloro Desmethylchlorphine
MF: $C_{19}H_{18}Cl_3N_3O \cdot HCl$
FW: 447.2
Purity: $\geq 98\%$
Supplied as: A solid
Storage: $-20^{\circ}C$
Stability: ≥ 4 years



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

Description

SR-17018 (hydrochloride) (Item No. 45831) is an analytical reference standard categorized as an opioid.¹ SR-17018 increases latency to withdrawal in the hot plate and warm water tail-flick assays without inducing respiratory depression in mice, as well as induces place preference in a conditioned place preference test in mice.^{1,2} It also induces antinociceptive tolerance after repeated administration in mice.² This product is intended for research and forensic applications.

References

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).
2. Kudla, L., Bugno, R., Podlewska, S., *et al.* Comparison of an addictive potential of μ -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, 06/09/2026

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