

PRODUCT INFORMATION



U-47700 (hydrochloride)

Item No. 21078

Formal Name: *trans*-3,4-dichloro-N-[2-(dimethylamino)cyclohexyl]-N-methyl-benzamide, monohydrochloride

MF: C₁₆H₂₂Cl₂N₂O • HCl

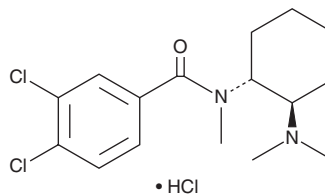
FW: 365.7

Purity: ≥98%

Supplied as: A neat solid

Storage: -20°C

Stability: ≥2 years



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

Description

U-47700 (hydrochloride) (Item No. 21078) is an analytical reference material categorized as a utopioid.¹ U-47700 has been associated with overdoses, one of which was reversed with naloxone, and several reported fatalities.^{2,3} U-47700 is regulated as a Schedule I compound in the United States. This product is intended for forensic and research 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.

References

1. Vasudevan, L., Vandeputte, M.M., Deventer, M., *et al.* Assessment of structure-activity relationships and biased agonism at the mu opioid receptor of novel synthetic opioids using a novel, stable bio-assay platform. *Biochem. Pharmacol.* **177**, 113910 (2020).
2. Schneir, A., Metushi, I.G., Sloane, C., *et al.* Near death from a novel synthetic opioid labeled U-47700: Emergence of a new opioid class. *Clin. Toxicol. (Phila)* **55(1)**, 51-54 (2017).
3. Dziadosz, M., Klintschar, M., and Teske, J. Postmortem concentration distribution in fatal cases involving the synthetic opioid U-47700. *Int. J. Legal Med.* **131(6)**, 1555-1556 (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, 10/12/2021

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