

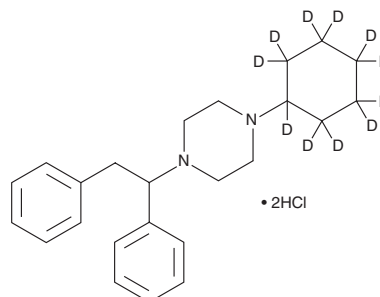
PRODUCT INFORMATION



MT-45-d₁₁ (hydrochloride) (exempt preparation)

Item No. 36876

CAS Registry No.: 2747914-21-0
Formal Name: 1-(cyclohexyl-d₁₁)-4-(1,2-diphenylethyl) piperazine, dihydrochloride
Synonym: l-C6-d₁₁
MF: C₂₄H₂₁D₁₁N₂ • 2HCl
FW: 432.5
Purity: ≥98%
Supplied as: A 100 µg/ml solution in methanol
Storage: -20°C
Stability: ≥3 years



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

Description

MT-45-d₁₁ (hydrochloride) (exempt preparation) (Item No. 36876) is an analytical reference standard intended for use as an internal standard for the quantification of MT-45 (Item Nos. 19737 | 14082) by GC- or LC-MS. MT-45 is a novel synthetic opioid that is structurally unrelated to most opioids but has analgesic activity similar to morphine (Item Nos. ISO60147 | 15464) in animal models. MT-45-d₁₁ is regulated as a Schedule I compound in the United States. MT-45-d₁₁ (hydrochloride) (exempt preparation) (Item No. 36876) is provided as a DEA exempt preparation. This product is intended for research and forensic applications.

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, 01/10/2025

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