

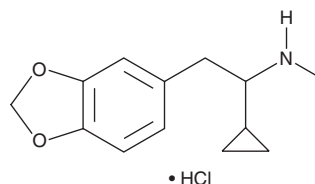
PRODUCT INFORMATION



UWA-101 (hydrochloride)

Item No. 15433

CAS Registry No.: 1431520-52-3
Formal Name: α -cyclopropyl-N-methyl-1,3-benzodioxole-5-ethanamine, monohydrochloride
MF: C₁₃H₁₇NO₂ • HCl
FW: 255.7
Purity: ≥98%
Supplied as: A neat solid
Storage: -20°C
Stability: ≥5 years



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

Description

UWA-101 (hydrochloride) (Item No. 15433) is an analytical reference material that is structurally categorized as an amphetamine. It inhibits both serotonin and dopamine reuptake (K_i s = 467 and 894 nM, respectively), but has no effect on the reuptake of noradrenaline (K_i > 50,000 nM).¹ UWA-101 enhances L-DOPA anti-parkinsonian action without exacerbating dyskinesia.² This product is intended for research and forensic 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. Johnston, T.H., Millar, Z., Huot, P., *et al.* A novel MDMA analogue, UWA-101, that lacks psychoactivity and cytotoxicity, enhances L-DOPA benefit in parkinsonian primates. *FASEB J.* **26(5)**, 2154-2163 (2012).
2. Huot, P., Johnston, T.H., Gandy, M.N., *et al.* The monoamine re-uptake inhibitor UWA-101 improves motor fluctuations in the MPTP-lesioned common marmoset. *PLoS One* **7(9)**, (2012).

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, 08/26/2019

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