

PRODUCT INFORMATION



XLR11 N-(4-hydroxypentyl) metabolite-d₅ Item No. 14377

Formal Name: (1-(5-fluoro-4-hydroxypentyl)-1H-indol-3-yl)
(2,2,3,3-tetramethylcyclopropyl)
methanone-2,4,5,6,7-d₅

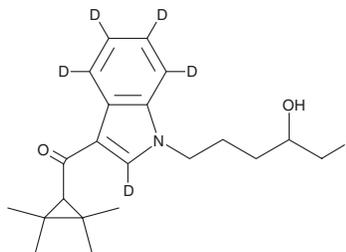
MF: C₂₁H₂₃D₅FNO₂

FW: 350.5

Purity: ≥98%

Supplied as: A neat oil

Storage: -20°C



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

Description

XLR11 N-(4-hydroxypentyl) metabolite-d₅ RM contains five deuterium atoms at the 2, 4, 5, 6, and 7 positions. It is intended for use as an internal standard for the quantification of XLR11 N-(4-hydroxypentyl) metabolite by GC- or LC-mass spectrometry. XLR11 (Item No. 11565) is a synthetic cannabinoid (CB) featuring a tetramethylcyclopropyl group, which reportedly confers selectivity for the peripheral CB₂ receptor over the central CB₁ receptor.¹ XLR11 also has an N-(5-fluoropentyl) chain, which increases binding to both CB receptors.² XLR11 N-(4-hydroxypentyl) metabolite is an expected phase 1 metabolite of XLR11, based on the known metabolism of similar compounds.^{3,4} The physiological properties of this compound have not been evaluated. 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.

References

1. Frost, J.M., Dart, M.J., Tietje, K.R., *et al.* Indol-3-ylcycloalkyl ketones: Effects of N1 substituted indole side chain variations on CB₂ cannabinoid receptor activity. *J. Med. Chem.* **53**, 295-315 (2010).
2. Makriyannis, A. and Deng, H. Cannabimimetic indole derivatives. WO 01/28557 A1 (2001), 1-25, PCT/US00/28832.
3. Zhang, Q., Ma, P., Cole, R.B., *et al.* Identification of *in vitro* metabolites of JWH-015, an aminoalkylindole agonist for the peripheral cannabinoid receptor (CB₂) by HPLC-MS/MS. *Anal. Bioanal. Chem.* **386**, 1345-1355 (2006).
4. Wintermeyer, A., Möller, I., Thevis, M., *et al.* *In vitro* phase I metabolism of the synthetic cannabimimetic JWH-018. *Anal. Bioanal. Chem.* **398**, 2141-53 (2010).

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, 08/12/2019

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