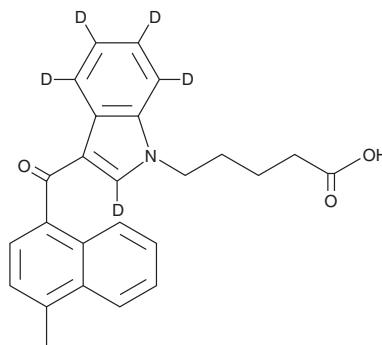


# PRODUCT INFORMATION



## MAM2201 N-pentanoic acid metabolite-d<sub>5</sub> Item No. 14366

**CAS Registry No.:** 2748469-43-2  
**Formal Name:** 5-(3-(4-methyl-1-naphthoyl)-1H-indol-1-yl)-2,4,5,6,7-d<sub>5</sub>pentanoic acid  
**Synonym:** JWH 122 N-pentanoic acid metabolite-d<sub>5</sub>  
**MF:** C<sub>25</sub>H<sub>18</sub>D<sub>5</sub>NO<sub>3</sub>  
**FW:** 390.5  
**Chemical Purity:** ≥98% (MAM2201 N-pentanoic acid metabolite)  
**Deuterium Incorporation:** ≥99% deuterated forms (d<sub>1</sub>- d<sub>5</sub>); ≤1% d<sub>0</sub>  
**UV/Vis.:** λ<sub>max</sub>: 221, 311 nm  
**Supplied as:** A solution in acetonitrile  
**Storage:** -20°C  
**Stability:** ≥4 years



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

### Description

MAM2201 N-pentanoic acid metabolite-d<sub>5</sub> (Item No. 14366) is intended for use as an internal standard for the quantification of MAM2201 N-pentanoic acid metabolite (Item No. 11779) by GC- or LC-MS. The accuracy of the sample weight in this vial is between 5% over and 2% under the amount shown on the vial. If better precision is required, the deuterated standard should be quantitated against a more precisely weighed unlabeled standard by constructing a standard curve of peak intensity ratios (deuterated versus unlabeled).

MAM2201 (Item No. 9001219) is a synthetic cannabinoid (CB) that is closely related, structurally, to AM2201 (Item No. 10707) and JWH 122 (Item No. 10591), two compounds which display high affinities for both CB receptors. MAM2201 N-pentanoic acid metabolite is a potential phase 1 metabolite of MAM2201 or JWH 122, based on the known metabolism of similar compounds. The physiological and toxicological properties of this compound have not been evaluated. This product is intended for forensic and research 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, 06/16/2023

#### 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