

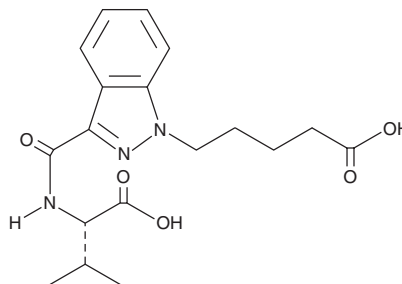
PRODUCT INFORMATION



5-fluoro AMB metabolite 6

Item No. 37058

CAS Registry No.: 1890250-15-3
Formal Name: N-[[1-(4-carboxybutyl)-1H-indazol-3-yl]carbonyl]-L-valine
Synonym: AMB metabolite 6
MF: C₁₈H₂₃N₃O₅
FW: 361.4
Purity: ≥98%
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

5-fluoro AMB metabolite 6 (Item No. 37058) is an analytical reference standard categorized as a synthetic cannabinoid metabolite.¹ 5-fluoro AMB metabolite 6 is a metabolite of 5-fluoro AMB (Item No. 15489) and AMB (Item No. 15488). This product is intended for research and forensic applications.

Reference

1. Andersson, M., Diao, X., Wohlfarth, A., *et al.* Metabolic profiling of new synthetic cannabinoids AMB and 5F-AMB by human hepatocyte and liver microsome incubations and high-resolution mass spectrometry. *Rapid Commun. Mass Spectrom.* **30**(8), 1067-1078 (2016).

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, 07/07/2023

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