

PRODUCT INFORMATION

para-hydroxy 2-methyl AP-237 (hydrochloride)

Item No. 34456

Formal Name: (E)-1-(4-(3-(4-hydroxyphenyl)allyl)-2-methylpiperazin-1-yl)butan-1-one, monohydrochloride

Synonyms: 4'-hydroxy 2-methyl AP-237, *p*-hydroxy 2-methyl AP-237

MF: C₁₈H₂₆N₂O₂ • HCl

FW: 338.9

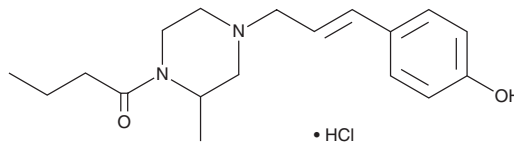
Purity: ≥98%

UV/Vis.: λ_{max}: 272 nm

Supplied as: A crystalline solid

Storage: -20°C

Stability: ≥5 years



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

Description

para-hydroxy 2-methyl AP-237 (hydrochloride) (Item No. 34456) is an analytical reference standard that is structurally similar to known opioids. *para*-hydroxy 2-methyl AP-237 is a potential metabolite of 2-methyl AP-237 (Item No. 26485) based on the published metabolism of AP-237 (Item No. 26484).¹ At the time *para*-hydroxy 2-methyl AP-237 (hydrochloride) (Item No. 34456) was made available for purchase, specific metabolism data had not been published. We suspect that this compound could be the first eluting metabolite of 2-methyl AP-237 (Item No. 26485) based on data presented in our 2-methyl AP-237 NPS Metabolism Monograph (Issue 1). Contact us if updated information on this molecule is now available. This product is intended for research and forensic applications.

Reference

1. Baba, S. and Morishita, S.-I. Studies on drug metabolism by use of isotopes. XVI. Species differences in metabolism of 1-butyryl-4-cinnamylpiperazine hydrochloride. *Chem. Pharm. Bull. (Tokyo)* **23(9)**, 1949-1954 (1975).

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, 06/23/2023

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