

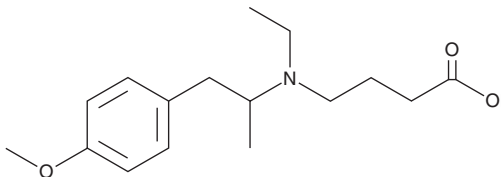
PRODUCT INFORMATION



Mebeverine Acid

Item No. 35498

CAS Registry No.: 475203-77-1
Formal Name: 4-[ethyl[2-(4-methoxyphenyl)-1-methylethyl]amino]-butanoic acid
MF: C₁₆H₂₅NO₃
FW: 279.4
Purity: ≥98%
UV/Vis.: λ_{max}: 226 nm
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years



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

Laboratory Procedures

Mebeverine acid is supplied as a solid. A stock solution may be made by dissolving the mebeverine acid in the solvent of choice, which should be purged with an inert gas. Mebeverine acid is soluble in DMSO.

Description

Mebeverine acid is a metabolite of the antispasmodic agent mebeverine.¹

Reference

1. Moskaleva, N.E., Baranov, P.A., Mesonzhnik, N.V., *et al.* HPLC-MS/MS method for the simultaneous quantification of desmethylmebeverine acid, mebeverine acid and mebeverine alcohol in human plasma along with its application to a pharmacokinetics study. *J. Pharm. Biomed. Anal.* **138**, 118-125 (2017).

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, 12/12/2022

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