

# PRODUCT INFORMATION



## O-Desmethyl Mebeverine Acid

Item No. 35523

CAS Registry No.: 586357-02-0

Formal Name: 4-[ethyl[2-(4-hydroxyphenyl)-1-methylethyl]amino]-butanoic acid

MF:  $C_{15}H_{23}NO_3$

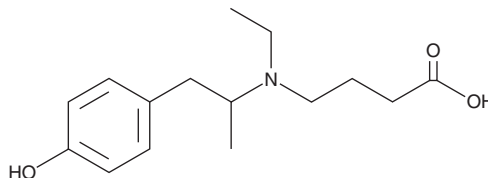
FW: 265.4

Purity:  $\geq 95\%$

Supplied as: A solid

Storage:  $-20^{\circ}\text{C}$

Stability:  $\geq 4$  years



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

### Laboratory Procedures

O-Desmethyl mebeverine acid is supplied as a solid. A stock solution may be made by dissolving the O-desmethyl mebeverine acid in the solvent of choice, which should be purged with an inert gas. O-Desmethyl mebeverine acid is slightly soluble in DMSO and is slightly soluble with sonication in methanol.

### Description

O-Desmethyl mebeverine acid is a metabolite of the antispasmodic agent mebeverine.<sup>1</sup> It is formed from mebeverine by hydrolysis and subsequent oxidation and demethylation.

### Reference

1. Moskaleva, N.E., Baranov, P.A., Mesonzhnik, N.V., *et al.* HPLC-MS/MS method for the simultaneous quantification of desmethyimebeverine 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