

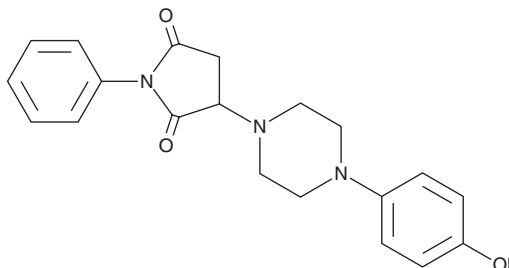
PRODUCT INFORMATION



Malic Enzyme Inhibitor ME1

Item No. 44383

CAS Registry No.: 522649-59-8
Formal Name: 3-[4-(4-hydroxyphenyl)-1-piperazinyl]-1-phenyl-2,5-pyrrolidinedione
MF: C₂₀H₂₁N₃O₃
FW: 351.4
Purity: ≥98%
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

Malic enzyme inhibitor ME1 is supplied as a solid. A stock solution may be made by dissolving the malic enzyme inhibitor ME1 in the solvent of choice, which should be purged with an inert gas. Malic enzyme inhibitor ME1 is soluble (≥10 mg/ml) in DMSO.

Description

Malic enzyme inhibitor ME1 is an inhibitor of malic enzyme (ME; IC₅₀s = 0.22, 0.2, and 0.15 μM for ME1, ME2, and ME3, respectively).¹

Reference

1. Sheth, G., Shah, S.R., Sengupta, P., *et al.* In the quest for potent and selective malic enzyme 3 inhibitors for the treatment of pancreatic ductal adenocarcinoma. *ACS Med. Chem. Lett.* **14**(1), 41-50 (2022).

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, 10/06/2025

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