

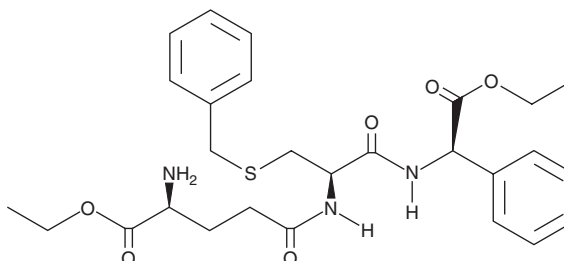
PRODUCT INFORMATION



Ezatiostat

Item No. 16248

CAS Registry No.: 168682-53-9
Formal Name: (2R)-L-γ-glutamyl-S-(phenylmethyl)-L-cysteinyl-2-phenyl-glycine, 1,3 diethyl ester
Synonyms: Terrapin 199, TLK199
MF: C₂₇H₃₅N₃O₆S
FW: 529.7
Purity: ≥98%
Supplied as: A crystalline 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

Ezatiostat is supplied as a crystalline solid. A stock solution may be made by dissolving the ezatiostat in the solvent of choice, which should be purged with an inert gas. Ezatiostat is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of ezatiostat in these solvents is approximately 20 mg/ml.

Ezatiostat is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, ezatiostat should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. Ezatiostat has a solubility of approximately 0.5 mg/ml in a 1:1 solution of ethanol:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Ezatiostat is an inhibitor of glutathione S-transferase pi 1-1 (GSTP1-1; $K_i = 0.4 \mu\text{M}$).¹ It is selective for GSTP1-1 over GST α and GST μ family GSTs (K_i s = 20-75 μM). Ezatiostat also inhibits the protein-protein interaction between GSTP1-1 and JNK1. *In vivo*, ezatiostat (75 mg/kg) increases the number of circulating white blood cells in *Gstp1* wild-type, but not *Gstp1*^{-/-}, mice.

Reference

1. Yount, J.S., Zhang, M.M., and Hang, H.C. Visualization and identification of fatty acylated proteins using chemical reporters. *Curr. Protoc. Chem. Biol.* **3**(2), 65-79 (2011).

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, 10/05/2022

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