

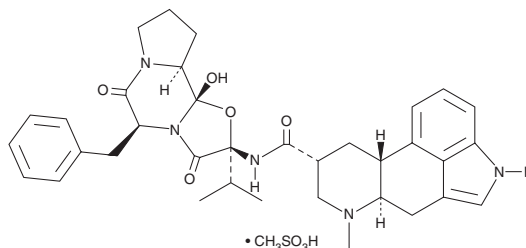
PRODUCT INFORMATION



Dihydroergocristine (mesylate)

Item No. 29492

CAS Registry No.: 24730-10-7
Formal Name: 9,10 α -dihydro-12'-hydroxy-2'-(1-methylethyl)-5' α -(phenylmethyl)-ergotaman-3',6',18-trione, monomethanesulfonate
MF: C₃₅H₄₁N₅O₅ • CH₃SO₃H
FW: 707.8
Purity: \geq 95%
Supplied as: A solid
Storage: -20°C
Stability: \geq 4 years



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

Laboratory Procedures

Dihydroergocristine (mesylate) is supplied as a solid. A stock solution may be made by dissolving the dihydroergocristine (mesylate) in the solvent of choice, which should be purged with an inert gas. Dihydroergocristine (mesylate) is slightly soluble in chloroform and methanol.

Description

Dihydroergocristine is an alkaloid with diverse biological activities.¹⁻⁴ It inhibits contraction of isolated rat vas deferens induced by norepinephrine ($pA_2 = 7.88$) but not calcium.¹ Dihydroergocristine (0.1-100 μ g/kg, i.v.) reduces blood pressure in a rat model of norepinephrine-induced hypertension and increases blood pressure in hypotensive rats.² It reduces aqueous humor production and intraocular pressure in a rabbit model of α -chymotrypsin-induced ocular hypertension.³ Dihydroergocristine also reduces hypermotility induced by ethanol in mice.⁴

References

1. Lucchelli, A., Zonta, F., Grana, E., *et al.* Dualism in the α -adrenoceptor blocking action of dihydroergocristine. *Pharmacol. Res. Commun.* **13(3)**, 231-240 (1981).
2. Morpurgo, C., Faini, D., and Falcone, A. Effects of phentolamine, dihydroergocristine and isoxsuprine on the blood pressure and heart rate in normotensive, hypotensive and hypertensive rats. *Naunyn Schmiedebergs Arch. Pharmacol.* **290(4)**, 335-346 (1975).
3. Melena, J., Santafe, J., and Segarra, J. The effect of topical dihydroergocristine on the intraocular pressure in alpha-chymotrypsin-induced ocular hypertensive rabbits. *Methods Find. Exp. Clin. Pharmacol.* **20(10)**, 861-867 (1998).
4. Boada, J.N., Fera, M., Serrano, M., *et al.* Inhibitory effect of dihydroergocristine, phenoxybenzamine and propranolol on the hypermotility induced by low doses of ethyl alcohol in mice. *Arch. Farmacol. Toxicol.* **5(2)**, 117-122 (1979).

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, 12/09/2022

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