

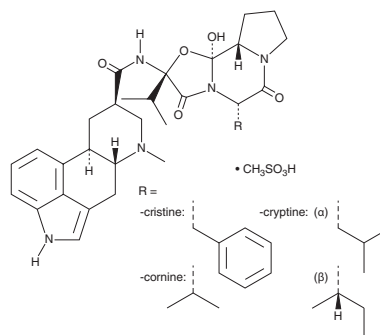
# PRODUCT INFORMATION



## Ergoloid (mesylates)

Item No. 24095

**CAS Registry No.:** 8067-24-1  
**Formal Name:** dihydro-ergotoxine, monomethanesulfonate  
**MF:** C<sub>134</sub>H<sub>184</sub>N<sub>20</sub>O<sub>32</sub>S<sub>4</sub>  
**FW:** 2,715.3  
**Purity:** ≥95%  
**UV/Vis.:** λ<sub>max</sub>: 222 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

Ergoloid (mesylates) is supplied as a solid. A stock solution may be made by dissolving the ergoloid (mesylates) in the solvent of choice. Ergoloid (mesylates) is soluble in organic solvents such as methanol, which should be purged with an inert gas. The solubility of ergoloid (mesylates) in methanol is approximately 1 mg/ml. Ergoloid (mesylates) is slightly soluble in DMSO.

### Description

Ergoloid is a mixture of the ergot alkaloids dihydroergocornine, dihydroergocristine, and dihydroergocryptine.<sup>1</sup> It blocks reflex vasodilation and vasoconstriction induced by norepinephrine (Item No. 16673) in dogs when injected into the kidney at a dose of 0.15 mg. Ergoloid blocks norepinephrine-induced cAMP production in rat brain cortical homogenates *ex vivo* in a dose-dependent manner.<sup>2</sup> It also acts as an erectogenic agent in rats.<sup>3</sup> Formulations containing ergoloid have been used in the treatment of migraine headaches and vascular dementias.

### References

1. Matsushita, S., Ying, L.H., Kaneko, J., *et al.* Pharmacological studies on reflex renal vasodilatation. *Tohoku J. Exp. Med.* **99(1)**, 51-57 (1969).
2. Markstein, R. and Wagner, H. The effect of dihydroergotoxin, phentolamine and pindolol on catecholamine-stimulated adenylyl cyclase in rat cerebral cortex. *FEBS Lett.* **55(1)**, 275-277 (1975).
3. Radosavljevic, M., Pajovic, B., Radunovic, M., *et al.* Influence of dihydroergotoxine, bromocriptine, and ergotamine on penile erection in Wistar rats. *J. Androl.* **33(5)**, 866-871 (2012).

#### 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, 11/15/2022

#### 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