

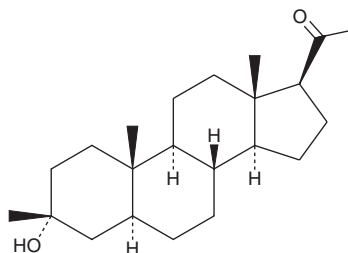
PRODUCT INFORMATION



Ganaxolone

Item No. 20919

CAS Registry No.: 38398-32-2
Formal Name: (3 α ,5 α)-3-hydroxy-3-methyl-pregnan-20-one
Synonym: CCD 1042
MF: C₂₂H₃₆O₂
FW: 332.5
Purity: \geq 98%
UV/Vis.: λ_{max} : 282 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: \geq 5 years



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

Description

Ganaxolone (Item No. 20919) is an analytical reference standard categorized as a GABA_A receptor positive allosteric modulator.¹ Ganaxolone has anticonvulsant activity in mice. Ganaxolone is regulated as a Schedule V compound in the United States. This product is intended for research and forensic applications.

Reference

1. Carter, R.B., Wood, P.L., Wieland, S., *et al.* Characterization of the anticonvulsant properties of ganaxolone (CCD 1042; 3 α -hydroxy-3 β -methyl-5 α -pregnan-20-one), a selective, high-affinity, steroid modulator of the γ -aminobutyric acid_A receptor. *J. Pharmacol. Exp. Ther.* **280**(3), 1284-1295 (1997).

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, 07/24/2023

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