

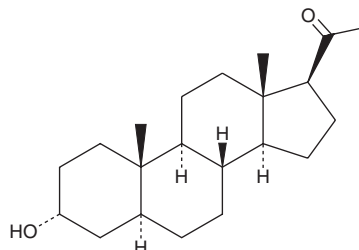
PRODUCT INFORMATION



Allopregnanolone

Item No. 16930

CAS Registry No.: 516-54-1
Formal Name: 3 α -hydroxy-5 α -pregnan-20-one
Synonyms: ALLO, Brexanolone,
3 α ,5 α -tetrahydro Progesterone,
3 α ,5 α -THP
MF: C₂₁H₃₄O₂
FW: 318.5
Purity: \geq 95%
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

Allopregnanolone (Item No. 16930) is an analytical reference standard categorized as a GABA_A and GABA_C receptor positive allosteric modulator.^{1,2} Allopregnanolone has anxiolytic-like and anticonvulsant activities in rodents.^{3,4} Formulations containing allopregnanolone have been used in the treatment of postpartum depression. Allopregnanolone is regulated as a Schedule IV compound in the United States. This product is intended for research and forensic applications.

References

1. Alvarez, L.D., Pecci, A., and Estrin, D.A. In search of GABA_A receptor's neurosteroid binding sites. *J. Med. Chem.* **62(11)**, 5250-5260 (2019).
2. Johnston, G.A.R., Chebib, M., Hanrahan, J.R., *et al.* Neurochemicals for the investigation of GABA_C receptors. *Neurochem. Res.* **35(12)**, 1970-1977 (2010).
3. Fernández-Guasti, A. and Picazo, O. Flumazenil blocks the anxiolytic action of allopregnanolone. *Eur. J. Pharmacol.* **281(1)**, 113-115 (1995).
4. Reddy, D.S., Castaneda, D.C., O'Malley, B.W., *et al.* Anticonvulsant activity of progesterone and neurosteroids in progesterone receptor knockout mice. *J. Pharmacol. Exp. Ther.* **10(1)**, 230-239 (2004).

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, 07/05/2023

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