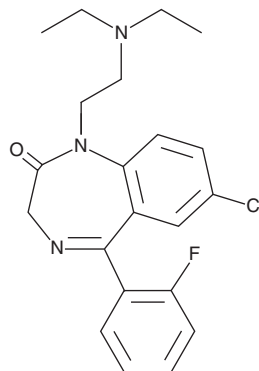


PRODUCT INFORMATION

Flurazepam

Item No. 16190

CAS Registry No.: 17617-23-1
Formal Name: 7-chloro-1-[2-(diethylamino)ethyl]-5-(2-fluorophenyl)-1,3-dihydro-2H-1,4-benzodiazepin-2-one
Synonyms: Dalmadorm, Dalmane, Ro 56901/3
MF: C₂₁H₂₃ClFN₃O
FW: 387.9
Purity: ≥98%
Supplied as: A neat oil
Storage: -20°C
Stability: ≥2 years



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

Description

Flurazepam (Item No. 16190) is an analytical reference material that is categorized as a benzodiazepine. It acts as a partial agonist of the benzodiazepine site on GABA_A receptors, potentiating the action of GABA with an EC₅₀ value of 930 nM in neuronal chick spinal cord cultures.¹ Through this signaling mechanism, flurazepam produces anxiolytic and sedative properties.² Formulations containing flurazepam have been used in the treatment of insomnia but also carry a potential for misuse.^{2,3} Flurazepam is regulated as a Schedule IV compound in the United States. This product is intended for research and forensic applications.

This product is qualified as a Reference Material that has been manufactured and tested to ISO/IEC 17025 and ISO 17034 international standards.

References

1. Chan, C.Y. and Farb, D.H. Modulation of neurotransmitter action: Control of the γ -aminobutyric acid response through the benzodiazepine receptor. *J. Neurosci.* **5**(9), 2365-2373 (1985).
2. Bertisch, S.M., Herzig, S.J., Winkelman, J.W., et al. National use of prescription medications for insomnia: NHANES 1999-2010. *Sleep* **37**(2), 343-349 (2014).
3. Griffiths, R.R. and Johnson, M.W. Relative abuse liability of hypnotic drugs: A conceptual framework and algorithm for differentiating among compounds. *J. Clin. Psychiatry* **66**(Suppl 9), 31-41 (2005).

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, 07/10/2019

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