

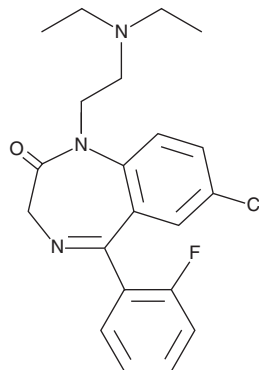
PRODUCT INFORMATION



Flurazepam

Item No. 41590

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
Synonym: Ro 56901/3
MF: C₂₁H₂₃ClFN₃O
FW: 387.9
Purity: ≥98%
Supplied as: A 25 mg/ml solution in methanol
Storage: -20°C
Stability: ≥4 years



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

Description

Flurazepam (Item No. 41590) is an analytical reference standard categorized as a benzodiazepine.¹ Flurazepam has sedative effects in mice.² Formulations containing flurazepam have been used in the treatment of insomnia. Flurazepam is regulated as a Schedule IV compound in the United States. This product is intended for research and forensic applications.

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. Marley, R.J., Freund, R.K., and Wehner, J.M. Differential response to flurazepam in long-sleep and short-sleep mice. *Pharmacol. Biochem. Behav.* **31**(2), 453-458 (1988).

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, 11/21/2025

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