

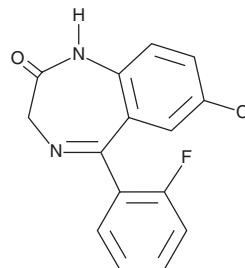
PRODUCT INFORMATION



Desalkylflurazepam

Item No. 18484

CAS Registry No.: 2886-65-9
Formal Name: 7-chloro-5-(2-fluorophenyl)-1,3-dihydro-2H-1,4-benzodiazepin-2-one
Synonyms: N-Desalkylflurazepam, N-Desalkyl-2-oxoquazepam, nor-Flurazepam, Norfludiazepam, Ro 5-3367
MF: C₁₅H₁₀ClFN₂O
FW: 288.7
Purity: ≥98%
Supplied as: A neat solid
Storage: -20°C
Stability: ≥2 years



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

Description

Desalkylflurazepam (Item No. 18484) is an analytical reference material that is structurally categorized as a benzodiazepine. It is an active metabolite of several benzodiazepines, including flurazepam (Item No. 18160), flutoprazepam, fludiazepam, midazolam, and quazepam. Desalkylflurazepam inhibits L-type voltage-gated calcium channels (Ca_v; IC₅₀s = 55 and 37 μM for Ca_v1.2 and 1.3, respectively) by positively modulating GABA_A receptors.¹ It can be detected in urine, serum, and meconium by LC-MS/MS.² 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. Earl, D.E. and Tietz, E.I. Inhibition of recombinant L-type voltage-gated calcium channels by positive allosteric modulators of GABA_A receptors. *J. Pharmacol. Exp. Ther.* **337**(1), 301-311 (2011).
2. Marin, S.J., Coles, R., Merrell, M., *et al.* Quantitation of benzodiazepines in urine, serum, plasma, and meconium by LC-MS-MS. *J. Anal. Toxicol.* **32**(7), 491-498 (2008).

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, 05/01/2020

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