

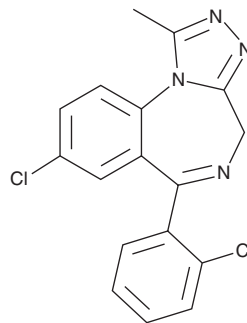
PRODUCT INFORMATION



Triazolam

Item No. 22577

CAS Registry No.: 28911-01-5
Formal Name: 8-chloro-6-(2-chlorophenyl)-1-methyl-4H-[1,2,4]triazolo[4,3-a][1,4]benzodiazepine
Synonym: U-33030
MF: C₁₇H₁₂Cl₂N₄
FW: 343.2
Purity: ≥98%
UV/Vis.: λ_{max}: 225 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥5 years



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

Description

Triazolam (Item No. 22577) is an analytical reference standard categorized as a benzodiazepine. It substitutes for cocaine (Item Nos. ISO60176 | 16186 | 22165) in baboons.¹ Formulations containing triazolam have been used in the treatment of insomnia. Triazolam is regulated as a Schedule IV compound in the United States. This product is intended for research and forensic applications.

Reference

1. Griffiths, R.R., Lamb, R.J., Sannerud, C.A., *et al.* Self-injection of barbiturates, benzodiazepines and other sedative-anxiolytics in baboons. *Psychopharmacology (Berl.)* **103(2)**, 154-161 (1991).

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, 01/04/2024

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