PRODUCT INFORMATION



Adinazolam

Item No. 18205

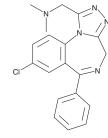
CAS Registry No.: 37115-32-5

Formal Name: 8-chloro-N,N-dimethyl-6-phenyl-

4H-[1,2,4]triazolo[4,3-a][1,4]

benzodiazepine-1-methanamine

MF: $C_{19}H_{18}CIN_5$ 351.8 FW: ≥98% **Purity:** Supplied as: A neat solid -20°C Storage: Stability: ≥4 years



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

Description

Adinazolam (Item No. 18205) is an analytical reference material categorized as a benzodiazepine. 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.

Reference

1. Fraser, A.D., Isner, A.F., and Bryan, W. Urinary screening for adinazolam and its major metabolites by the Emit d.a.u. and FPIA benzodiazepine assays with confirmation by HPLC. J. Anal. Toxicol. 17(7), 427-431 (1993).

WARNING
THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

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.

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, 04/06/2020

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