

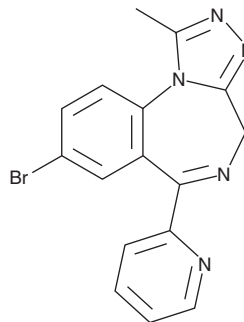
PRODUCT INFORMATION



Pyrazolam

Item No. 14901

CAS Registry No.: 39243-02-2
Formal Name: 8-bromo-1-methyl-6-(2-pyridinyl)-4H-[1,2,4]triazolo[4,3-a][1,4]benzodiazepine
MF: C₁₆H₁₂BrN₅
FW: 354.2
Purity: ≥98%
UV/Vis.: λ_{max}: 222 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

Pyrazolam (Item No. 14901) is an analytical reference standard categorized as a benzodiazepine.¹ Pyrazolam has been abused recreationally and has anxiolytic-like effects in mice.^{1,2} This product is intended for research and forensic applications.

References

1. Moosmann, B., Huppertz, L.M., Hutter, M., *et al.* Detection and identification of the designer benzodiazepine flubromazepam and preliminary data on its metabolism and pharmacokinetics. *J. Mass. Spectrom.* **48(11)**, 1150-1159 (2013).
2. Hester, J.B., Jr. and Von Voigtlander, P. 6-Aryl-4H-s-triazolo[4,3- α][1,4]benzodiazepines. Influence of 1-substitution on pharmacological activity. *J. Med. Chem.* **22(11)**, 1390-1398 (1979).

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, 09/27/2023

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