

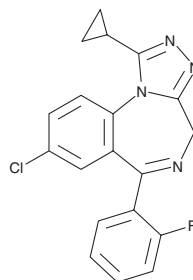
PRODUCT INFORMATION



Cyclopropylflualprazolam

Item No. 45621

CAS Registry No.: 2101840-38-2
Formal Name: 8-chloro-1-cyclopropyl-6-(2-fluorophenyl)-H-[1,2,4]triazolo[4,3-a][1,4]benzodiazepine
MF: C₁₉H₁₄ClFN₄
FW: 352.8
Purity: ≥98%
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years



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

Description

Cyclopropylflualprazolam (Item No. 45621) is an analytical reference standard categorized as a benzodiazepine.¹ Cyclopropylflualprazolam has been found in pressed candies marketed as sleep aids. This product is intended for research and forensic applications.

Reference

1. Liu, X., Zhao, L., Dong, P., et al. Isolation and identification of the designer benzodiazepine cyclopropylflualprazolam in pressed candies marketed for sleep improvement. *J. Pharm. Biomed. Anal.* 275, 117459 (2026).

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, 05/05/2026

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