

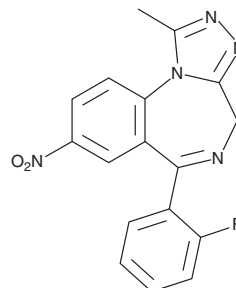
PRODUCT INFORMATION



Flunitrazolam

Item No. 23108

CAS Registry No.: 2243815-18-9
Formal Name: 6-(2-fluorophenyl)-1-methyl-8-nitro-4H-[1,2,4]triazolo[4,3-a][1,4]benzodiazepine
MF: C₁₇H₁₂FN₅O₂
FW: 337.3
Purity: ≥98%
UV/Vis.: λ_{max}: 227, 247 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

Flunitrazolam (Item No. 23108) is an analytical reference standard categorized as a benzodiazepine.¹ Flunitrazolam has been found in biological samples from intoxication cases.² This product is intended for research and forensic applications.

References

1. Ameline, A., Richeval, C., Gaulier, J.M., *et al.* Characterization of flunitrazolam, a new designer benzodiazepine, in oral fluid after a controlled single administration. *J. Anal. Toxicol.* (2018).
2. Helander, A., Bäckberg, M., Signell, P., *et al.* Intoxications involving acrylfentanyl and other novel designer fentanyls - results from the Swedish STRIDA project. *Clin. Toxicol. (Phila)* **55(6)**, 589-599 (2017).

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, 07/05/2023

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