

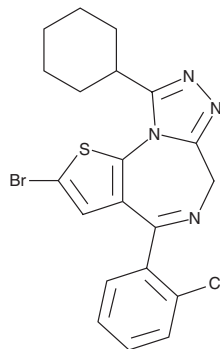
PRODUCT INFORMATION



Ciclotizolam

Item No. 37999

CAS Registry No.: 58765-21-2
Formal Name: 2-bromo-4-(2-chlorophenyl)-9-cyclohexyl-6H-thieno[3,2-f][1,2,4]triazolo[4,3-a][1,4]diazepine
Synonym: WE-973
MF: C₂₀H₁₈BrClN₄S
FW: 461.8
Purity: ≥98%
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥2 years



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

Description

Ciclotizolam (Item No. 37999) is an analytical reference standard categorized as a thienotriazolodiazepine.¹ Ciclotizolam has anticonvulsant and anti-aggressive activities in mice and decreases total sleep time in cats. This product is intended for research and forensic applications.

Reference

1. K.H., W., Kuhn, F.J., Böke-Kuhn, K., *et al.* Pharmacological and neurochemical properties of 1,4-diazepines with two annelated heterocycles ('hetrazepines'). *Eur. J. Pharmacol.* **109(1)**, 19-31 (1985).

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, 03/29/2023

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