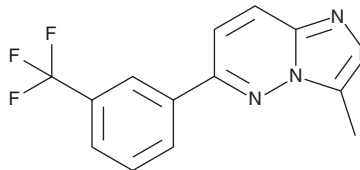


PRODUCT INFORMATION



CL 218,872
Item No. 40450

CAS Registry No.: 66548-69-4
Formal Name: 3-methyl-6-[3-(trifluoromethyl)phenyl]-
1,2,4-triazolo[4,3-b]pyridazine
MF: C₁₃H₉F₃N₄
FW: 278.2
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

CL 218,872 (Item No. 40450) is an analytical reference standard categorized as a triazolopyridazine.¹ CL 218,872 has anxiolytic and sedative effects in mice.² This product is intended for analytical forensic applications. This product is also available as a general research tool (Item No. 34037).

References

1. Yamamura, H.I., Mimaki, T., Yamamura, S.H., et al. [³H]CL 218,872, a novel triazolopyridazine which labels the benzodiazepine receptors in rat brain. *Eur. J. Pharmacol.* **77(4)**, 351-354 (1982).
2. Oakley, N.R., Jones, B.J., and Straughan, D.W. The benzodiazepine receptor ligand CL218,872 has both anxiolytic and sedative properties in rodents. *Neuropharmacology* **23(7A)**, 797-802 (1984).

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, 02/02/2024

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