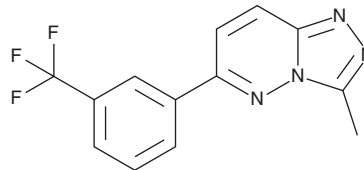


# PRODUCT INFORMATION



CL 218,872  
Item No. 34037

CAS Registry No.: 66548-69-4  
Formal Name: 3-methyl-6-[3-(trifluoromethyl)phenyl]-1,2,4-triazolo[4,3-b]pyridazine  
MF: C<sub>13</sub>H<sub>9</sub>F<sub>3</sub>N<sub>4</sub>  
FW: 278.2  
Purity: ≥95%  
UV/Vis.: λ<sub>max</sub>: 247 nm  
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.

## Laboratory Procedures

CL 218,872 is supplied as a solid. A stock solution may be made by dissolving the CL 218,872 in the solvent of choice, which should be purged with an inert gas. CL 218,872 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of CL 218,872 in these solvents is approximately 1 mg/ml.

## Description

CL 218,872 is an anxiolytic agent.<sup>1</sup> It selectively binds to  $\alpha_1\beta_1\gamma_2$  subunit-containing GABA<sub>A</sub> receptors over  $\alpha_2\beta_1\gamma_2$  or  $\alpha_3\beta_1\gamma_2$  subunit-containing GABA<sub>A</sub> receptors ( $K_{iS}$  = 0.131, 1.82, and 1.53  $\mu$ M, respectively).<sup>2</sup> CL 218,872 increases drinking in the Vogel punished drinking task and inhibits foot-shock-induced fighting, indicating anxiolytic-like activities, in mice in a dose-dependent manner.<sup>1</sup> It also increases the duration of pentobarbitone-induced anesthesia, reduces spontaneous locomotor activity, and inhibits convulsions induced by pentylenetetrazole (PTZ; Item No. 18682), picrotoxin (Item No. 20771), or bicuculline (Item No. 11727) with ED<sub>50</sub> values of 12.8, 28.5, and 35.3 mg/kg, respectively, in mice.<sup>1</sup>

## References

1. 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).
2. Pritchett, D.B., Lüddens, H., and Seeburg, P.H. Type I and type II GABA<sub>A</sub>-benzodiazepine receptors produced in transfected cells. *Science* **245(4924)**, 1389-1392 (1989).

**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, 11/14/2022

## 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