

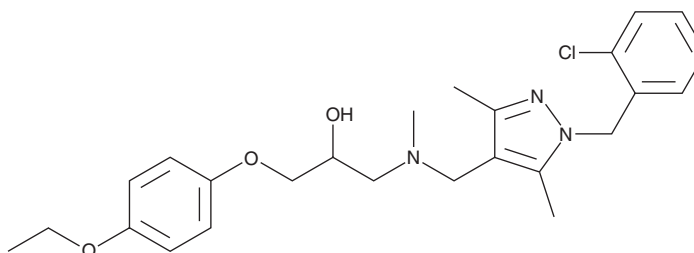
PRODUCT INFORMATION



T-5342126

Item No. 33099

CAS Registry No.: 956507-49-6
Formal Name: 1-[[[1-[(2-chlorophenyl)methyl]-3,5-dimethyl-1H-pyrazol-4-yl]methyl]methylamino]-3-(4-ethoxyphenoxy)-2-propanol
MF: C₂₅H₃₂ClN₃O₃
FW: 458.0
Purity: ≥98%
Supplied as: A solution in methyl acetate
Storage: -20°C
Stability: ≥2 years



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

Description

T-5342126 is a toll-like receptor 4 (TLR4) antagonist.¹ It reduces LPS-induced production of nitric oxide (NO) in RAW 264.7 cells (IC₅₀ = 27.8 μM), as well as decreases LPS-induced IL-8, TNF-α, and IL-6 production in isolated human whole blood (IC₅₀s = 110.5, 315.6, and 318.4 μM, respectively). T-5342126 (82 mg/kg) reduces ethanol intake and the abundance of ionized calcium-binding adapter molecule 1 (Iba1), a marker of microglial activation, in the central nucleus of the amygdala in ethanol-dependent mice.²

References

1. Chavez, S.A., Martinko, A.J., Lau, C., *et al.* Development of β-amino alcohol derivatives that inhibit toll-like receptor 4 mediated inflammatory response as potential antiseptics. *J. Med. Chem.* **54(13)**, 4659-4669 (2011).
2. Bajo, M., Montgomery, S.E., Cates, L.N., *et al.* Evaluation of TLR4 inhibitor, T5342126, in modulation of ethanol-drinking behavior in alcohol-dependent mice. *Alcohol Alcohol.* **51(5)**, 541-548 (2016).

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, 11/01/2023

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