

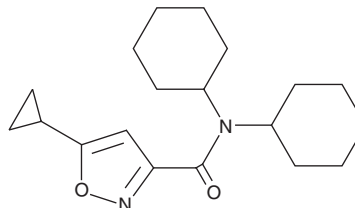
PRODUCT INFORMATION



CYM 5541

Item No. 15190

CAS Registry No.: 945128-26-7
Formal Name: N,N-dicyclohexyl-5-cyclopropyl-3-isoxazolecarboxamide
Synonyms: CID-17253208, ML-249
MF: C₁₉H₂₈N₂O₂
FW: 316.4
Purity: ≥98%
UV/Vis.: λ_{max}: 232 nm
Supplied as: A crystalline 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

CYM 5541 is supplied as a crystalline solid. A stock solution may be made by dissolving the CYM 5541 in the solvent of choice, which should be purged with an inert gas. CYM 5541 is soluble in the organic solvent DMSO at a concentration of approximately 5 mg/ml.

Description

The sphingosine 1-phosphate (S1P) receptor S1P₃ has roles in vasodilation, epithelial barrier function, cardioprotection, fibrosis, immunity, and inflammation.^{1,2} CYM 5541 is a selective, allosteric agonist of S1P₃ (EC₅₀ = 72-132 nM).^{3,4} It less effectively activates S1P₁ (EC₅₀ = 28-38 μM) and is ineffective at S1P₂, S1P₄, and S1P₅ as well as a panel of other receptors, ion channels and transporters.³

References

1. Choi, J.W., Lee, C.W., and Chun, J. Biological roles of lysophospholipid receptors revealed by genetic null mice: An update. *Biochim. Biophys. Acta* **1781**, 531-539 (2008).
2. Rosen, H., Stevens, R.C., Hanson, M., et al. Sphingosine-1-phosphate and its receptors: Structure, signaling, and influence. *Annu. Rev. Biochem.* **82**, 637-662 (2013).
3. Guerrero, M., Poddutoori, R., Pinacho-Crisostomo, F., et al. Probe development efforts for an allosteric agonist of the sphingosine 1-phosphate receptor 3 (S1P₃), in Probe Reports from the NIH Molecular Libraries Program,¹ (2011).
4. Jo, E., Bhatarai, B., Repetto, E., et al. Novel selective allosteric and bitopic ligands for the S1P3 receptor. *ACS Chem. Biol.* **7(12)**, 1975-1983 (2012).

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, 11/04/2022

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