

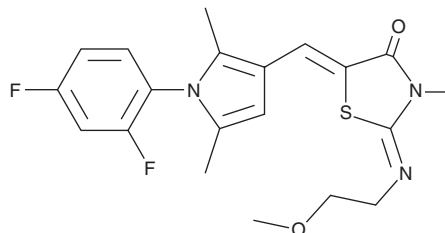
PRODUCT INFORMATION



CYM 50308

Item No. 14667

CAS Registry No.: 1345858-76-5
Formal Name: (2Z,5Z)-5-[[1-(2,4-difluorophenyl)-2,5-dimethyl-1H-pyrrol-3-yl]methylene]-2-[(2-methoxyethyl)imino]-3-methyl-4-thiazolidinone
Synonyms: CID-49835928, ML-248
MF: C₂₀H₂₁F₂N₃O₂S
FW: 405.5
Purity: ≥98%
UV/Vis.: λ_{max}: 363 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 50308 is supplied as a crystalline solid. A stock solution may be made by dissolving the CYM 50308 in the solvent of choice, which should be purged with an inert gas. CYM 50308 is soluble in the organic solvent DMSO at a concentration of approximately 1 mg/ml.

Description

CYM 50308 is a novel agonist of S1P₄ (EC₅₀ = 56 nM) that displays 37-fold selectivity against S1P₅ (EC₅₀ = 2.1 μM) and no appreciable activity over the S1P₁₋₃ subtypes at concentrations up to 25 μM.¹

Reference

1. Urbano, M., Guerrero, M., Velaparthi, S., *et al.* Discovery, synthesis and SAR analysis of novel selective small molecule S1P₄-R agonists based on a (2Z,5Z)-5-((pyrrol-3-yl)methylene)-3-alkyl-2-(alkylimino)thiazolidin-4-one chemotype. *Bioorg. Med. Chem. Lett.* **21(22)**, 6739-6745 (2011).

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, 11/10/2022

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