

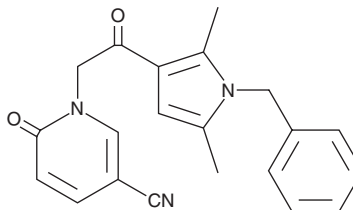
PRODUCT INFORMATION



CYM 5520

Item No. 17638

CAS Registry No.: 1449747-00-5
Formal Name: 1-[2-[2,5-dimethyl-1-(phenylmethyl)-1H-pyrrol-3-yl]-2-oxoethyl]-1,6-dihydro-6-oxo-3-pyridinecarbonitrile
MF: C₂₁H₁₉N₃O₂
FW: 345.4
Purity: ≥95%
UV/Vis.: λ_{max}: 254, 301 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥2 years



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

Laboratory Procedures

CYM 5520 is supplied as a crystalline solid. A stock solution may be made by dissolving the CYM 5520 in the solvent of choice, which should be purged with an inert gas. CYM 5520 is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of CYM 5520 in these solvents is approximately 30 mg/ml.

Description

Sphingosine-1-phosphate (S1P; Item No. 62570) is an extracellular lipid mediator whose major effects are mediated through five distinct G protein-coupled receptors, S1P₁/EDG-1, S1P₂/EDG-5, S1P₃/EDG-3, S1P₄/EDG-6, and S1P₅/EDG-8. S1P₂ has been implicated in multiple biological functions, including Rho activation, inhibition of Rac and cell migration, and in “feed forward” autocrine signaling in NF-κB survival signaling of tumor cells.¹ CYM 5520 is a noncompetitive allosteric agonist of S1P₂ (EC₅₀ = 0.48 μM) that does not display agonist activity towards S1P₁, S1P₃, S1P₄, or S1P₅.¹

Reference

1. Satsu, H., Schaeffer, M.T., Guerrero, M., *et al.* A sphingosine 1-phosphate receptor 2 selective allosteric agonist. *Bioorg. Med. Chem.* **21(17)**, 5373-5382 (2013).

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, 02/21/2023

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