

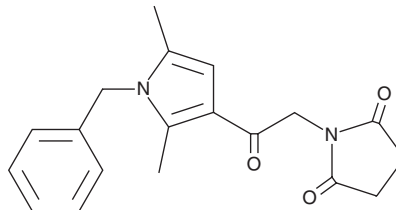
PRODUCT INFORMATION



ML-031

Item No. 15171

CAS Registry No.: 852230-33-2
Formal Name: 1-[2-[2,5-dimethyl-1-(phenylmethyl)-1H-pyrrol-3-yl]-2-oxoethyl]-2,5-pyrrolidinedione
Synonym: CID-2113511
MF: C₁₉H₂₀N₂O₃
FW: 324.4
Purity: ≥95%
UV/Vis.: λ_{max}: 251, 293 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

ML-031 is supplied as a crystalline solid. A stock solution may be made by dissolving the ML-031 in the solvent of choice, which should be purged with an inert gas. ML-031 is soluble in organic solvents such as ethanol and DMSO. The solubility of ML-031 in these solvents is approximately 5 mg/ml.

Description

Sphingosine-1-phosphate (S1P) is a bioactive lipid that exhibits a broad spectrum of biological activities including cell proliferation, survival, migration, cytoskeletal organization, and morphogenesis. It exerts its activity by binding to five distinct G protein-coupled receptors, S1P₁/EDG-1, S1P₂/EDG-5, S1P₃/EDG-3, S1P₄/EDG-6, and S1P₅/EDG-8. ML-031 is one of the first identified agonists of S1P₂ demonstrating an EC₅₀ value of 1 μM in a S1P reporter assay.^{24537}

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, 10/20/2022

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