PRODUCT INFORMATION



CYM 5478

Item No. 29024

CAS Registry No.: 870762-83-7

Formal Name: 1-[2-[2,5-dimethyl-1-

> (phenylmethyl)-1H-pyrrol-3-yl]-2-oxoethyl]-5-(trifluoromethyl)-

2(1H)-pyridinone

MF: $C_{21}H_{19}F_3N_2O_2$

FW: 388.4 **Purity:** ≥98% UV/Vis.: λ_{max} : 297 nm Supplied as: A crystalline solid

-20°C Storage: Stability: ≥2 years

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

Laboratory Procedures

CYM 5478 is supplied as a crystalline solid. A stock solution may be made by dissolving the CYM 5478 in the solvent of choice, which should be purged with an inert gas. CYM 5478 is soluble in the organic solvent DMSO.

Description

CYM 5478 is a sphingosine-1-phosphate receptor 2 (S1P₂) agonist (EC₅₀ = 0.78 μ M in a reporter assay).¹ It is selective for S1P₂ over S1P₁ and S1P₃₋₅ in a TGF α -shedding assay (EC₅₀s = 119, 1,690, 1,950, >10,000, and >10,000 nM, respectively). CYM 5478 (0.1-10 μM) reduces serum starvation-induced decreases in C6 rat glioma cell viability. It also reduces accumulation of reactive oxygen species (ROS) and apoptosis induced by cisplatin (Item No. 13119) in C6 cells when used at a concentration of 10 μM.

Reference

1. Chen, Y.L., Obach, R.S., Braselton, J., et al. 2-Aryloxy-4-alkylaminopyridines: Discovery of novel corticotropin-releasing factor 1 antagonists. J. Med. Chem. 51(5), 1385-1392 (2008).

WARNING
THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

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.

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, 12/4/2019

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