

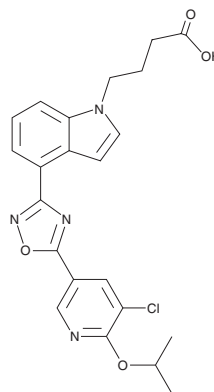
PRODUCT INFORMATION



GSK2018682

Item No. 31437

CAS Registry No.: 1034688-30-6
Formal Name: 4-[5-[5-chloro-6-(1-methylethoxy)-3-pyridinyl]-1,2,4-oxadiazol-3-yl]-1H-indole-1-butanoic acid
MF: C₂₂H₂₁ClN₄O₄
FW: 440.9
Purity: ≥95%
UV/Vis.: λ_{max}: 225, 272, 302 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

GSK2018682 is supplied as a crystalline solid. A stock solution may be made by dissolving the GSK2018682 in the solvent of choice, which should be purged with an inert gas. GSK2018682 is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of GSK2018682 in these solvents is approximately 30 mg/ml.

GSK2018682 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, GSK2018682 should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. GSK2018682 has a solubility of approximately 0.25 mg/ml in a 1:3 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

GSK2018682 is an agonist of sphingosine-1-phosphate receptor 1 (S1P₁) and S1P₅.¹ It is selective for S1P₁ and S1P₅ over S1P₂, S1P₃, and S1P₄ receptors. GSK2018682 is efficacious in a mouse model of experimental autoimmune encephalomyelitis (EAE).

Reference

1. Xu, J., Gray, F., Henderson, A., *et al.* Safety, pharmacokinetics, pharmacodynamics, and bioavailability of GSK2018682, a sphingosine-1-phosphate receptor modulator, in healthy volunteers. *Clin. Pharmacol. Drug Dev.* 3(3), 170-178 (2014).

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, 10/27/2022

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