

PRODUCT INFORMATION

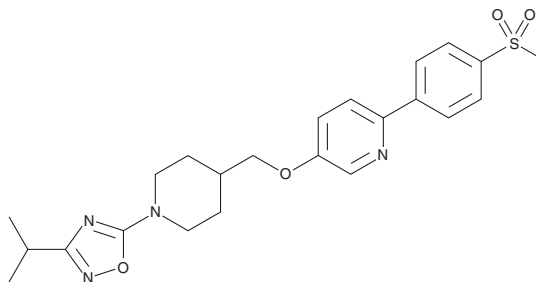


GSK1292263

Item No. 26181

CAS Registry No.: 1032823-75-8
Formal Name: 5-[[1-[3-(1-methylethyl)-1,2,4-oxadiazol-5-yl]-4-piperidinyl]methoxy]-2-[4-(methylsulfonyl)phenyl]-pyridine

MF: C₂₃H₂₈N₄O₄S
FW: 456.6
Purity: ≥98%
UV/Vis.: λ_{max}: 224, 276, 296 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

GSK1292263 is supplied as a crystalline solid. A stock solution may be made by dissolving the GSK1292263 in the solvent of choice. GSK1292263 is soluble in the organic solvent DMSO, which should be purged with an inert gas, at a concentration of approximately 30 mg/ml.

Description

GSK1292263 is an agonist of G protein-coupled receptor 119 (GPR119; EC₅₀ = 43.9 nM in CHO-K1 cells expressing human receptors).¹ It inhibits increases in sterol regulatory element binding protein-1 (SREBP-1) levels induced by the LXR agonist T0901317 and increases activation of AMPK in HepG2 cells in a concentration-dependent manner.² GSK1292263 also inhibits breast cancer resistant protein (BCRP) and organic anion transporting polypeptide 1B1 (OATP1B1) by 52 and 40%, respectively, when used at a concentration of 3 μM.³

References

1. Park, Y.H., Choi, H.H., Lee, D.H., *et al.* YH18421, a novel GPR119 agonist exerts sustained glucose lowering and weight loss in diabetic mouse model. *Arch. Pharm. Res.* **40(6)**, 772-782 (2017).
2. Yang, J.W., Kim, H.S., Im, J.H., *et al.* GPR119: A promising target for nonalcoholic fatty liver disease. *FASEB J.* **30(1)**, 324-335 (2016).
3. Polli, J.W., Hussey, E., Bush, M., *et al.* Evaluation of drug interactions of GSK1292263 (a GPR119 agonist) with statins: From *in vitro* data to clinical study design. *Xenobiotica* **43(6)**, 498-508 (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, 11/03/2022

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