

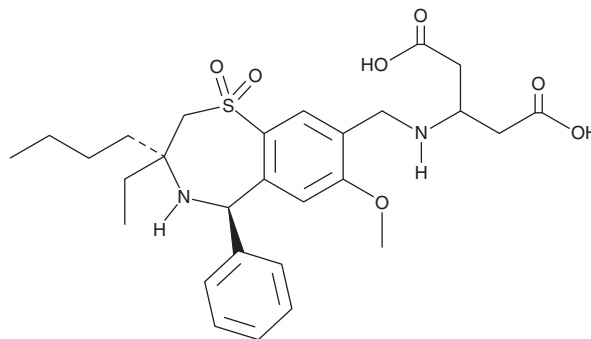
PRODUCT INFORMATION



GSK2330672

Item No. 23843

CAS Registry No.: 1345982-69-5
Formal Name: 3-[[[(3R,5R)-3-butyl-3-ethyl-2,3,4,5-tetrahydro-7-methoxy-1,1-dioxido-5-phenyl-1,4-benzothiazepin-8-yl]methyl]amino]-pentanedioic acid
MF: C₂₈H₃₈N₂O₇S
FW: 546.7
Purity: ≥98%
UV/Vis.: λ_{max}: 206, 238 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

GSK2330672 is supplied as a crystalline solid. A stock solution may be made by dissolving the GSK2330672 in the solvent of choice, which should be purged with an inert gas. GSK2330672 is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of GSK2330672 in these solvents is approximately 2 and 5 mg/ml, respectively.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of GSK2330672 can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of GSK2330672 in PBS (pH 7.2) is approximately 0.2 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

GSK2330672 is an inhibitor of the sodium/bile acid and sulfated solute cotransporter (ASBT; IC₅₀s = 42, 2.1, and 1.9 nM for human, mouse, and rat ASBT, respectively).¹ *In vivo*, GSK2330672 (0.05 mg/kg) stimulates fecal bile acid secretion, reduces hemoglobin A1c (HbA1c) and plasma glucose levels, and increases total GLP-1 (Item No. 24460) and plasma insulin in Zucker diabetic fatty (ZDF) rats.

Reference

1. Wu, Y., Aquino, C.J., Cowan, D.J., *et al.* Discovery of a highly potent, nonabsorbable apical sodium-dependent bile acid transporter inhibitor (GSK2330672) for treatment of type 2 diabetes. *J. Med. Chem.* **56**(12), 5094-5114 (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, 11/30/2022

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