

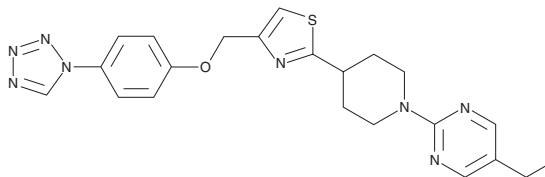
PRODUCT INFORMATION



MBX-2982

Item No. 22206

CAS Registry No.: 1037792-44-1
Formal Name: 5-ethyl-2-[4-[4-[[4-(1H-tetrazol-1-yl)phenoxy]methyl]-2-thiazolyl]-1-piperidinyl]-pyrimidine
MF: $C_{22}H_{24}N_8OS$
FW: 448.5
Purity: $\geq 98\%$
UV/Vis.: λ_{max} : 250 nm
Supplied as: A crystalline solid
Storage: $-20^{\circ}C$
Stability: ≥ 4 years



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

Laboratory Procedures

MBX-2982 is supplied as a crystalline solid. A stock solution may be made by dissolving the MBX-2982 in the solvent of choice, which should be purged with an inert gas. MBX-2982 is soluble in the organic solvent chloroform at a concentration of approximately 30 mg/ml.

Description

MBX-2982 is an agonist of GPR119.¹ It reduces nuclear and total protein levels of sterol regulatory element binding protein 1 (SREBP-1) in HepG2 cells and rat primary hepatocytes under high-glucose and -insulin conditions and increases phosphorylation of the inhibitory form, SREBP-1c. MBX-2982 (10 mg/kg) inhibits hepatic lipid accumulation in wild-type, but not GPR119 knockout, mice fed a high-fat diet. It also increases plasma levels of glucagon-like peptide 1 (GLP-1; Item No. 24460) in mice when administered at a dose of 10 mg/kg prior to, and to a greater extent following, glucose administration.² MBX-2982 increases glucokinase activity in an enzyme assay with an EC_{50} value of $45.11 \mu M$.³

References

1. 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).
2. Lan, H., Lin, H.V., Wang, C.F., *et al.* Agonists at GPR119 mediate secretion of GLP-1 from mouse enteroendocrine cells through glucose-independent pathways. *Br. J. Pharmacol.* **165**(8), 2799-2807 (2012).
3. Spasov, A.A., Kosolapov, V.A., Babkov, D.A., *et al.* Effect of GRP119 receptor agonist, compound MBX-2982, on activity of human glucokinase. *Bull. Exp. Biol. Med.* **163**(5), 695-698 (2017).

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, 12/16/2022

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