

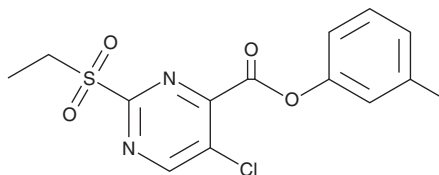
PRODUCT INFORMATION



SBI-115

Item No. 39168

CAS Registry No.: 882366-16-7
Formal Name: 5-chloro-2-(ethylsulfonyl)-
4-pyrimidinecarboxylic acid
3-methylphenyl ester
MF: C₁₄H₁₃ClN₂O₄S
FW: 340.8
Purity: ≥98%
Supplied as: A 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

SBI-115 is supplied as a solid. A stock solution may be made by dissolving the SBI-115 in the solvent of choice, which should be purged with an inert gas. SBI-115 is soluble in acetonitrile.

Description

SBI-115 is an antagonist of G protein-coupled bile acid receptor 1 (GPBAR1), also known as TGR5.^{49821} It inhibits tauro lithocholic acid-induced increases in cAMP levels in, and proliferation of, cholangiocytes isolated from patients with autosomal dominant polycystic kidney disease when used at concentrations of 100 and 200 μM. SBI-115 (10 μM) decreases the viability of, and protein levels of GPBAR1 in, PANC-1 pancreatic cancer cells.^{49822} *In vivo*, SBI-115 (80 mg/kg) prevents *B. pseudocatenuatum*-induced decreases in serum levels of IFN-γ and IL-17A in a mouse model of collagen-induced arthritis.^{49823}

References

1. Masyuk, T.V., Masyuk, A.I., Pisarello, M.L., *et al.* TGR5 contributes to hepatic cystogenesis in rodents with polycystic liver diseases through cyclic adenosine monophosphate/Gas signaling. *Hepatology* **66**(4), 1197-1218 (2017).
2. Lei, Y., Li, G., Li, J., *et al.* Investigation of the potential role of TGR5 in pancreatic cancer by a comprehensive molecular experiments and the liquid chromatography mass spectrometry (LC-MS) based metabolomics. *Discov. Oncol.* **13**(46), (2022).
3. Zhao, Q., Ren, H., Yang, N., *et al.* *Bifidobacterium pseudocatenuatum*-mediated bile acid metabolism to prevent rheumatoid arthritis via the gut-Joint axis. *Nutrients* **15**(2), 255 (2023).

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, 12/06/2023

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