

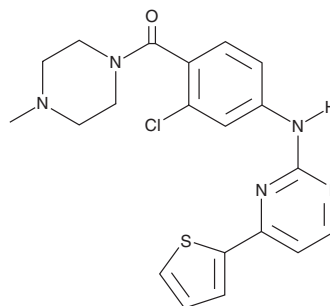
PRODUCT INFORMATION



GSK3 β Inhibitor XVIII

Item No. 21198

CAS Registry No.: 1139875-74-3
Formal Name: [2-chloro-4-[[4-(2-thienyl)-2-pyrimidinyl]amino]phenyl](4-methyl-1-piperazinyl)-methanone
MF: C₂₀H₂₀ClN₅OS
FW: 413.9
Purity: $\geq 98\%$
UV/Vis.: λ_{max} : 288 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

GSK3 β inhibitor XVIII is supplied as a crystalline solid. A stock solution may be made by dissolving the GSK3 β inhibitor XVIII in the solvent of choice, which should be purged with an inert gas. GSK3 β inhibitor XVIII is soluble in the organic solvent DMSO.

Description

GSK3 β inhibitor XVIII is an inhibitor of glycogen synthase kinase 3 β (GSK3 β ; IC₅₀ = 64 nM).¹ It induces expression of a luciferase reporter in HEK293 cells expressing wild-type Wnt (EC₅₀ = ~1.25 μ M) but not in cells expressing a β -catenin/TCF binding site-mutant Wnt. GSK3 β inhibitor XVIII increases β -catenin accumulation in the nucleus of HEK293 cells. It induces proliferation of R7T1 β -cells (EC₅₀ = 1,143 nM), an effect that can be blocked by the COX and Wnt signaling inhibitor sulindac (Item No. 10004386), and rat primary β -cells.

Reference

1. Wang, W., Walker, J.R., Wang, X., *et al.* Identification of small-molecule inducers of pancreatic β -cell expansion. *Proc. Nat. Acad. Sci. USA* **106**(5), 1427-1432 (2009).

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, 11/16/2022

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