

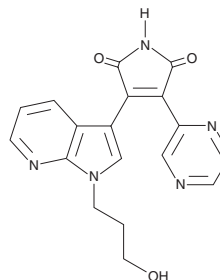
PRODUCT INFORMATION



GSK3 β Inhibitor XI

Item No. 23940

CAS Registry No.: 626604-39-5
Formal Name: 3-[1-(3-hydroxypropyl)-1H-pyrrolo[2,3-b]pyridin-3-yl]-4-(2-pyrazinyl)-1H-pyrrole-2,5-dione
Synonym: Glycogen Synthase Kinase 3 β Inhibitor XI
MF: C₁₈H₁₅N₅O₃
FW: 349.3
Purity: $\geq$ 98%
Supplied as: A solid
Storage: -20°C
Stability: $\geq$ 4 years



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

Laboratory Procedures

GSK3 β inhibitor XI is supplied as a solid. A stock solution may be made by dissolving the GSK3 β inhibitor XI in the solvent of choice, which should be purged with an inert gas. GSK3 β inhibitor XI is soluble in the organic solvent DMSO at a concentration of approximately 5 mg/ml.

Description

GSK3 β inhibitor XI is a potent inhibitor of glycogen synthase kinase 3 β (GSK3 β ; K_i = 25 nM).¹ It is selective for GSK3 β over a panel of 79 kinases at a concentration of 10 μ M. GSK3 β inhibitor XI inhibits GSK3 β in HEK293 cells (EC₅₀ = 32 nM).

Reference

1. O'Neill, D.J., Shen, L., Prouty, C., *et al.* Design, synthesis, and biological evaluation of novel 7-azaindolyheteroaryl-maleimides as potent and selective glycogen synthase kinase-3 β (GSK-3 β) inhibitors. *Bioorg. Med. Chem.* **12**(12), 3167-3185 (2004).

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, 10/29/2024

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