

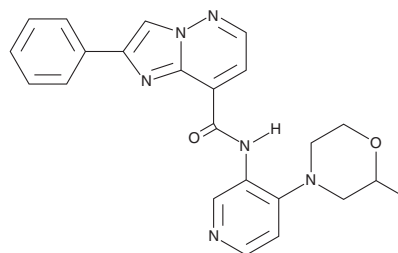
PRODUCT INFORMATION



GSK3β Inhibitor XIII

Item No. 45310

CAS Registry No.: 2227316-74-5
Formal Name: N-[4-(2-methyl-4-morpholinyl)-3-pyridinyl]-2-phenyl-imidazo[1,2-b]pyridazine-8-carboxamide
Synonyms: Glycogen Synthase Kinase β Inhibitor XIII, Glycogen Synthase Kinase β Inhibitor 13, GSK3β Inhibitor 13
MF: C₂₃H₂₂N₆O₂
FW: 414.5
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

GSK3β inhibitor XIII is supplied as a solid. A stock solution may be made by dissolving the GSK3β inhibitor XIII in the solvent of choice, which should be purged with an inert gas. GSK3β inhibitor XIII is slightly soluble (0.1-1 mg/ml) in acetonitrile.

Description

GSK3β inhibitor XIII is an inhibitor of glycogen synthase kinase β (GSK3β; IC₅₀ = 0.73 nM).¹ It is selective for GSK3β over cyclin-dependent kinase 2 (Cdk2), Cdk5, AP2-associated protein kinase 1 (AAK1), Pim-1, and PKCθ (IC₅₀s = 1.4, 6.4, 0.2, 0.38, and 0.36 μM, respectively) and a panel of 30 additional kinases (IC₅₀s = >2 μM for all). *In vivo*, GSK3β inhibitor XIII (30 mg/kg) reduces brain tau phosphorylation in a triple-transgenic mouse model of Alzheimer's disease.

Reference

1. Hartz, R.A., Ahuja, V.T., Sivaprakasam, P., *et al.* Design, structure-activity relationships, and *in vivo* evaluation of potent and brain-penetrant imidazo[1,2-b]pyridazines as glycogen synthase kinase-3β (gsk-3β) inhibitors. *J. Med. Chem.* **66**(6), 4321-4252 (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, 04/13/2026

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