

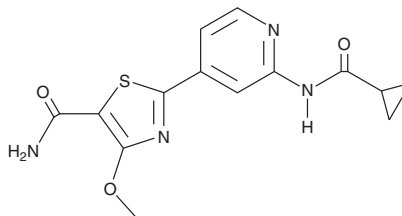
PRODUCT INFORMATION



GSK3 β Inhibitor 2

Item No. 43327

CAS Registry No.: 1702428-31-6
Formal Name: 2-[2-[(cyclopropylcarbonyl)amino]-4-pyridinyl]-4-methoxy-5-thiazolecarboxamide
Synonyms: Glycogen Synthase Kinase 3 β Inhibitor 2, HY 130795
MF: C₁₄H₁₄N₄O₃S
FW: 318.4
Purity: $\geq 95\%$
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

GSK3 β inhibitor 2 is supplied as a solid. A stock solution may be made by dissolving the GSK3 β inhibitor 2 in the solvent of choice, which should be purged with an inert gas. GSK3 β Inhibitor 2 is sparingly soluble (1-10 mg/ml) in DMSO.

Description

GSK3 β inhibitor 2 is glycogen synthase kinase 3 β (GSK3 β) inhibitor (IC₅₀ = 1.1 nM).¹ It decreases brain levels of phosphorylated tau in the 3xTg mouse model of Alzheimer's disease when administered at a dose of 30 mg/kg as a nanosuspension

Reference

1. Sivaprakasam, P., Han, X., Civiello, R.L., *et al.* Discovery of new acylaminopyridines as GSK-3 inhibitors by a structure guided in-depth exploration of chemical space around a pyrrolopyridinone core. *Bioorg. Med. Chem. Lett.* **25**(9), 1856-1863 (2015).

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, 04/28/2025

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