

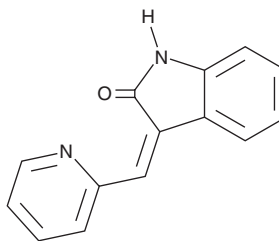
PRODUCT INFORMATION



GSK3 β Inhibitor 1

Item No. 42107

CAS Registry No.: 187325-53-7
Formal Name: (3Z)-1,3-dihydro-3-(2-pyridinylmethylene)-2H-indol-2-one
Synonym: Glycogen Synthase Kinase 3 β Inhibitor I
MF: C₁₄H₁₀N₂O
FW: 222.2
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 1 is supplied as a solid. A stock solution may be made by dissolving the GSK3 β inhibitor 1 in the solvent of choice, which should be purged with an inert gas. GSK3 β inhibitor 1 is sparingly soluble (1-10 mg/ml) in DMSO and slightly soluble (0.1-1 mg/ml) in ethanol.

Description

GSK3 β inhibitor 1 is an inhibitor of glycogen synthase kinase 3 β (GSK3 β ; IC₅₀ = 4.19 nM).¹ *In vivo*, GSK3 β inhibitor 1 (50 mg/kg) decreases blood glucose levels in an oral glucose tolerance test (OGTT) in a rat model of diabetes induced by a high-fat diet and streptozotocin (STZ; Item No. 13104).

Reference

1. Lozinskaya, N.A., Babkov, D.A., Zaryanova, E.V., *et al.* Synthesis and biological evaluation of 3-substituted 2-oxindole derivatives as new glycogen synthase kinase 3 β inhibitors. *Bioorg. Med. Chem.* **27(9)**, 1804-1817 (2019).

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