

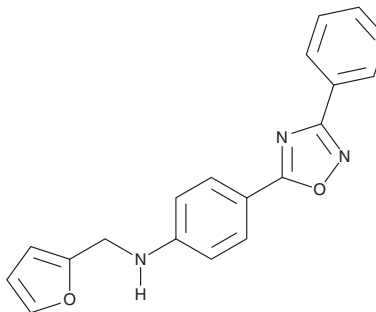
PRODUCT INFORMATION



SKL1223

Item No. 42952

Formal Name: N-(furan-2-ylmethyl)-4-(3-phenyl-1,2,4-oxadiazol-5-yl)aniline
MF: C₁₉H₁₅N₃O₂
FW: 317.3
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

SKL1223 is supplied as a solid. A stock solution may be made by dissolving the SKL1223 in the solvent of choice, which should be purged with an inert gas. SKL1223 is sparingly soluble (1-10 mg/ml) in DMSO and slightly soluble (0.1-1 mg/ml) in ethanol.

Description

SKL1223 is an inhibitor of thioredoxin-interacting protein (TXNIP).¹ It inhibits TXNIP promoter activity in a reporter assay using INS-1 cells expressing human TXNIP (IC₅₀ = 0.64 μM) and decreases the levels of TXNIP in the same cells in a concentration-dependent manner. SKL1223 (5 or 20 μM) induces glucagon secretion from TC-1 mouse lung cancer cells and decreases glucagon-induced glucose production in primary mouse hepatocytes. It decreases fasting blood levels of glucose, serum levels of glucagon and triglycerides, and hepatic levels of triglycerides in mice when administered at a dose of 70 mg/kg per day. SKL1223 decreases blood levels of glucose and serum levels of glucagon in *db/db* mice but not in *Txnip*^{-/-} mice.

Reference

1. Wang, X., Luo, Z., Han, J., *et al.* Orally available small molecule regulates TXNIP expression and glucagon action for the treatment of diabetes. *Research Square* (2024).

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, 05/15/2025

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