

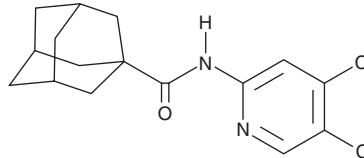
# PRODUCT INFORMATION



## Insulin Sensitizer C59

Item No. 42129

CAS Registry No.: 2761446-81-3  
Formal Name: N-(4,5-dichloro-2-pyridinyl)-  
tricyclo[3.3.1.1<sup>3,7</sup>]decane-1-carboxamide  
MF: C<sub>16</sub>H<sub>18</sub>Cl<sub>2</sub>N<sub>2</sub>O  
FW: 325.2  
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

Insulin sensitizer C59 is supplied as a solid. A stock solution may be made by dissolving the insulin sensitizer C59 in the solvent of choice, which should be purged with an inert gas. Insulin sensitizer C59 is soluble (≥10 mg/ml) in DMSO and sparingly soluble (1-10 mg/ml) in ethanol.

### Description

Insulin sensitizer C59 is a potentiator of insulin-induced plasma membrane translocation of glucose transporter 4 (Glut4).<sup>1</sup> It enhances insulin-stimulated increases in plasma membrane levels of Glut4 in a reporter assay using C2C12 cells when used at a concentration of 10 μM. Insulin sensitizer C59 (10 mg/kg per day) decreases fasting blood levels of glucose, fasting plasma levels of insulin, and body weight in mice fed a high-fat diet. It potentiates insulin-induced increases in the plasma membrane translocation of Glut4 in the forelimbs, hind limbs, brown adipose tissues, and heart but not brain or liver in mice in a model of diet-induced obesity.

### Reference

1. Yin, T.C., Van Vranken, J.G., Srivastava, D., *et al.* Insulin sensitization by small molecules enhancing GLUT4 translocation. *Cell Chem. Biol.* **30(8)**, 933-942 (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, 11/01/2024

#### 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