

PRODUCT INFORMATION

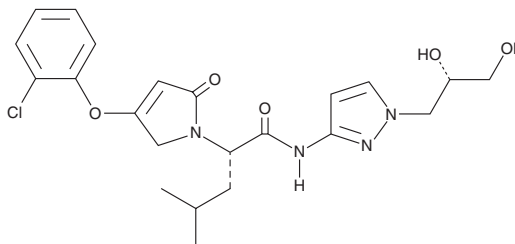


Sinogliatin

Item No. 43205

CAS Registry No.: 1191996-10-7
Formal Name: 4-(2-chlorophenoxy)-N-[1-[(2S)-2,3-dihydroxypropyl]-1H-pyrazol-3-yl]-2,5-dihydro- α S-(2-methylpropyl)-2-oxo-1H-pyrrole-1-acetamide

Synonym: HMS5552
MF: C₂₂H₂₇ClN₄O₅
FW: 462.9
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

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

Description

Sinogliatin is an activator of glucokinase.¹ It increases the hepatic levels of glucokinase and pancreatic levels of insulin in a rat model of diabetes induced by streptozotocin (Item No. 13104) and a high-fat diet (HFD) when administered at doses of 10 or 30 mg/kg per day. Sinogliatin (10 or 30 mg/kg) decreases blood levels of glucose in an oral glucose tolerance test (OGTT) in the same model. Formulations containing sinogliatin have been used in the treatment of type 2 diabetes.

Reference

1. Wang, P., Liu, H., Chen, L., *et al.* Effects of a novel glucokinase activator, HMS5552, on glucose metabolism in a rat model of type 2 diabetes mellitus. *J. Diabetes Res.* 5812607 (2017).

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, 04/09/2025

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