

PRODUCT INFORMATION



MK-0941

Item No. 25642

CAS Registry No.: 1137916-97-2
Formal Name: 3-[[6-(ethylsulfonyl)-3-pyridinyl]oxy]-5-[(1S)-2-hydroxy-1-methylethoxy]-N-(1-methyl-1H-pyrazol-3-yl)-benzamide, monomethanesulfonate

MF: C₂₁H₂₄N₄O₆S • CH₃SO₃H

FW: 556.6

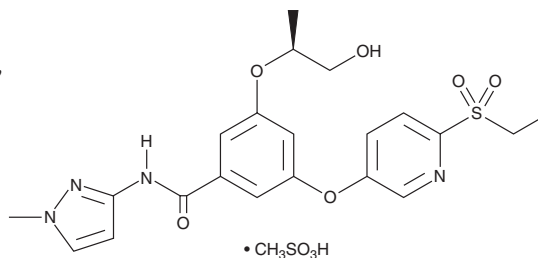
Purity: ≥98%

UV/Vis.: λ_{max}: 214, 276 nm

Supplied as: A crystalline 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

MK-0941 is supplied as a crystalline solid. A stock solution may be made by dissolving the MK-0941 in the solvent of choice, which should be purged with an inert gas. MK-0941 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of MK-0941 in ethanol is approximately 5 mg/ml and approximately 20 mg/ml in DMSO and DMF.

Description

MK-0941 is a glucokinase activator with EC₅₀ values of 0.24 and 0.065 μM at low (2.5 mM) and high (10 mM) concentrations of glucose, respectively.¹ It increases insulin secretion from isolated rat islets of Langerhans and glucose uptake by hepatocytes *in vitro*. MK-0941 reduces blood glucose levels in high-fat diet-fed *db/db* mice and in high-fat diet-fed mice with diabetes induced by streptozotocin (Item No. 13104).

Reference

1. Eiki, J.-i., Nagata, Y., Futamura, M., *et al.* Pharmacokinetic and pharmacodynamic properties of the glucokinase activator MK-0941 in rodent models of type 2 diabetes and healthy dogs. *Mol. Pharmacol.* **80**(6), 1156-1165 (2011).

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, 11/18/2022

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