

PRODUCT INFORMATION



AM4668

Item No. 43419

CAS Registry No.: 1011531-27-3
Formal Name: β S-[4-[[4-methyl-2-[4-(trifluoromethyl)phenyl]-5-thiazolyl]methoxy]phenyl]-3-isoxazolepropanoic acid

MF: C₂₄H₁₉F₃N₂O₄S

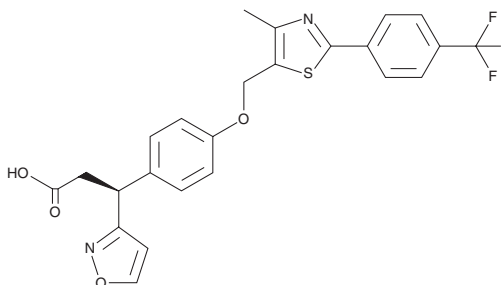
FW: 488.5

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

AM4668 is supplied as a solid. A stock solution may be made by dissolving the AM4668 in the solvent of choice, which should be purged with an inert gas. AM4668 is slightly soluble (0.1-1 mg/ml) in acetonitrile and DMSO.

Description

AM4668 is an agonist of free fatty acid receptor 1 (FFAR1/GPR40).¹ It induces inositol-1,4,5-triphosphate (IP₃) production in A9 cells expressing human FFAR1/GPR40 (EC₅₀ = 3.6 nM). AM4668 induces insulin secretion from pancreatic islets isolated from FFAR1 knock-in mice. *In vivo*, AM4668 (10 mg/kg) reduces blood glucose levels in FFAR1 knock-in mice.

Reference

1. Liu, J.J., Wang, Y., Ma, Z., *et al.* Optimization of GPR40 agonists for type 2 diabetes. *ACS Med. Chem. Lett.* 5(5), 517-521 (2014).

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, 06/17/2025

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