

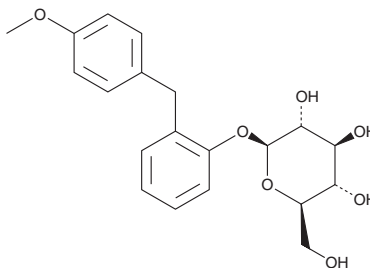
PRODUCT INFORMATION



Sergliflozin A

Item No. 14342

CAS Registry No.: 360775-96-8
Formal Name: 2-[(4-methoxyphenyl)methyl]
phenyl, β -D-glucopyranoside
MF: $C_{20}H_{24}O_7$
FW: 376.4
Purity: $\geq 98\%$
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

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

Description

Sergliflozin A is an inhibitor of sodium-glucose cotransporter 2 (SGLT2; $K_i = 2.39$ nM for the human transporter) and an active metabolite of sergliflozin etabonate.¹ It is formed from sergliflozin etabonate by esterases and selectively inhibits SGLT2 over SGLT1 ($K_i = 708$ nM for the human transporter). Co-infusion of sergliflozin A and glucose reduces maximal renal glucose reabsorption rates in rats.

Reference

1. Katsuno, K., Fujimori, Y., Takemura, Y., *et al.* Sergliflozin, a novel selective inhibitor of low-affinity sodium glucose cotransporter (SGLT2), validates the critical role of SGLT2 in renal glucose reabsorption and modulates plasma glucose level. *J. Pharmacol. Exp. Ther.* **320**(1), 323-330 (2007).

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, 09/23/2024

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