

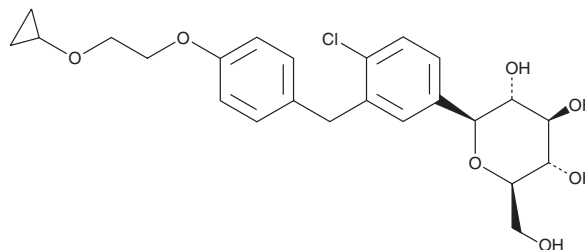
PRODUCT INFORMATION



EGT-1442

Item No. 21359

CAS Registry No.: 1118567-05-7
Formal Name: (1S)-1,5-anhydro-1-C-[4-chloro-3-[[4-[2-(cyclopropyloxy)ethoxy]phenyl]methyl]phenyl]-D-glucitol
Synonym: THR-1442
MF: C₂₄H₂₉ClO₇
FW: 464.9
Purity: ≥95%
UV/Vis.: λ_{max}: 225, 277 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

EGT-1442 is supplied as a crystalline solid. A stock solution may be made by dissolving the EGT-1442 in the solvent of choice, which should be purged with an inert gas. EGT-1442 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of EGT-1442 in these solvents is approximately 30 mg/ml.

EGT-1442 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, EGT-1442 should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. EGT-1442 has a solubility of approximately 0.5 mg/ml in a 1:1 solution of ethanol:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

EGT-1442 is a potent, selective sodium glucose co-transporter 2 (SGLT2) inhibitor with IC₅₀ values of 5.6 μM and 2 nM for human SGLT1 and SGLT2, respectively.¹ It produces a stable urinary excretion of glucose in rats and dogs with ED₅₀ values of 0.38 and 0.09 mg/kg, respectively, and reduces HbA(1c) and blood glucose in *db/db* mice in a concentration dependent manner.

Reference

1. Zhang, W., Welihinda, A., Mechanic, J., *et al.* EGT1442, a potent and selective SGLT2 inhibitor, attenuates blood glucose and HbA1c levels in *db/db* mice and prolongs the survival of stroke-prone rats. *Pharmacol. Res.* **63(4)**, 284-293 (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, 10/19/2022

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