

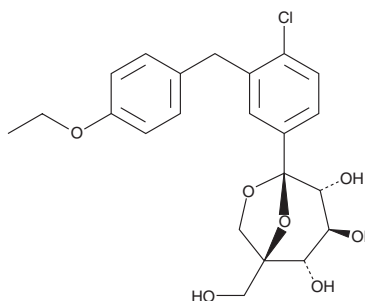
PRODUCT INFORMATION



Ertugliflozin

Item No. 21507

CAS Registry No.: 1210344-57-2
Formal Name: 1,6-anhydro-1-C-[4-chloro-3-[(4-ethoxyphenyl)methyl]phenyl]-5-C-(hydroxymethyl)-β-L-idopyranose
Synonym: PF-04971729
MF: C₂₂H₂₅ClO₇
FW: 436.9
Purity: ≥98%
UV/Vis.: λ_{max}: 224, 278 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

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

Ertugliflozin is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, ertugliflozin should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. Ertugliflozin 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

Ertugliflozin is an inhibitor of sodium-glucose cotransporter 2 (SGLT2; IC₅₀ = 0.927 nM for the human enzyme).¹ It is selective for SGLT2 over SGLT1 (IC₅₀ = 2,050 nM for the human enzyme). Ertugliflozin increases urinary glucose excretion in rats in a dose-dependent manner. Formulations containing ertugliflozin have been used in the treatment of type 2 diabetes.

Reference

1. Mascitti, V., Maurer, T.S., Robinson, R.P., *et al.* Discovery of a clinical candidate from the structurally unique dioxo-bicyclo[3.2.1]octane class of sodium-dependent glucose cotransporter 2 inhibitors. *J. Med. Chem.* **54**(8), 2952-2960 (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, 10/27/2022

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