

PRODUCT INFORMATION

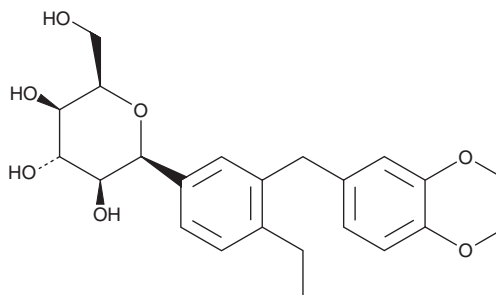


Licogliflozin (mixed isomers)

Item No. 35772

CAS Registry No.: 2730989-98-5
Formal Name: (2S,3S,4S,5R,6R)-2-(3-((2,3-dihydrobenzo[b][1,4]dioxin-6-yl)methyl)-4-ethylphenyl)-6-(hydroxymethyl)tetrahydro-2H-pyran-3,4,5-triol

Synonym: LIK-066
MF: C₂₃H₂₈O₇
FW: 416.5
Purity: ≥90% (mixture of isomers)
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

Licogliflozin (mixed isomers) is supplied as a solid. A stock solution may be made by dissolving the licogliflozin (mixed isomers) in the solvent of choice, which should be purged with an inert gas. Licogliflozin (mixed isomers) is soluble in DMSO.

Description

Licogliflozin (mixed isomers) is a mixture of isomers of the sodium-glucose cotransporter 1 (SGLT1) and SGLT2 inhibitor licogliflozin.

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, 12/13/2022

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