

PRODUCT INFORMATION



rac-trans-4-hydroxy Glyburide

Item No. 33752

CAS Registry No.: 23155-00-2
Formal Name: 5-chloro-N-[2-[4-[[[(*trans*-4-hydroxycyclohexyl)amino]carbonyl]amino]sulfonyl]phenyl]ethyl]-2-methoxy-benzamide

Synonyms: 4-*trans*-hydroxycyclohexyl Glyburide, 4-*trans*-hydroxy Glibenclamide

MF: C₂₃H₂₈ClN₃O₆S

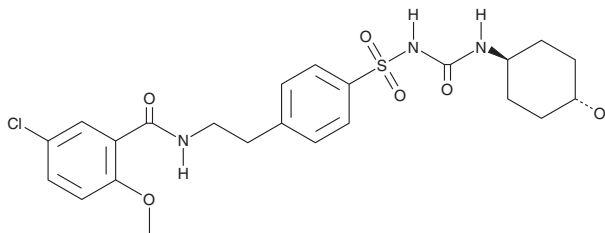
FW: 510.0

Purity: ≥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

rac-trans-4-hydroxy Glyburide is supplied as a solid. A stock solution may be made by dissolving the *rac-trans*-4-hydroxy glyburide in the solvent of choice, which should be purged with an inert gas. *rac-trans*-4-hydroxy Glyburide is slightly soluble in DMSO and methanol.

Description

rac-trans-4-hydroxy Glyburide is an active metabolite of the SUR1/K_{ir}6.2 sulfonylurea inhibitor glyburide (Item No. 15009).^{1,2} It is formed from glyburide by the cytochrome P450 (CYP) isoforms CYP2C8 and CYP2C9.¹ *rac-trans*-4-hydroxy Glyburide inhibits glyburide binding to rat brain synaptosomes at the high and low affinity sites of SUR1/K_{ir}6.2 with IC₅₀ values of 0.95 and 100 nM, respectively.²

References

1. Zharikova, O.L., Fokina, V.M., Nanovskaya, T.N., *et al.* Identification of the major human hepatic and placental enzymes responsible for the biotransformation of glyburide. *Biochem. Pharmacol.* **78(12)**, 1483-1490 (2009).
2. Hill, R.A., Rudra, S., Peng, B., *et al.* Hydroxyl-substituted sulfonylureas as potent inhibitors of specific [³H]glyburide binding to rat brain synaptosomes. *Bioorg. Med. Chem.* **11(9)**, 2099-2113 (2003).

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, 11/07/2022

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