

## PRODUCT INFORMATION

### ***trans*-hydroxy Glimepiride**

Item No. 34278

**CAS Registry No.:** 600177-94-4

**Formal Name:** 3-ethyl-2,5-dihydro-N-[2-[4-[[[*trans*-4-(hydroxymethyl)cyclohexyl]amino]carbonyl]amino]sulfonyl]phenyl]ethyl]-4-methyl-2-oxo-1H-pyrrole-1-carboxamide

MF:  $C_{24}H_{34}N_4O_6S$

FW: 506.6

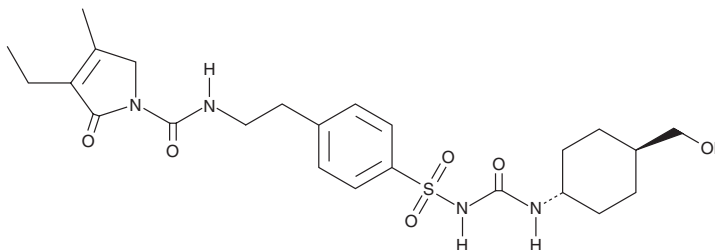
Purity:  $\geq 95\%$ 

**Supplied as:** A solid

**Storage:** -20°C

**Stability:**  $\geq 4$  years

Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.



## Laboratory Procedures

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

### Description

*trans*-hydroxy Glimepiride is an active metabolite of the sulfonylurea glimepiride (Item No. 12090).<sup>1</sup> It is formed from glimepiride primarily in the liver by the cytochrome P450 (CYP) isoform CYP2C9.

## Reference

1. Langtry, H.D. and Balfour, J.A. Glimepiride. A review of its use in the management of type 2 diabetes mellitus. *Drugs* **55**(4), 563-584 (1998).

### 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.

**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**