

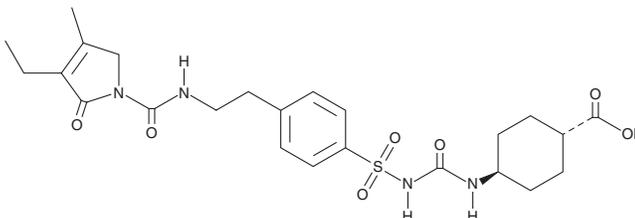
PRODUCT INFORMATION



trans-carboxy Glimepiride

Item No. 34277

CAS Registry No.: 1217739-04-2
Formal Name: *trans*-4-[[[[4-[2-[[[3-ethyl-2,5-dihydro-4-methyl-2-oxo-1H-pyrrol-1-yl]carbonyl]amino]ethyl]phenyl]sulfonyl]amino]carbonyl]amino]-cyclohexanecarboxylic acid
MF: C₂₄H₃₂N₄O₇S
FW: 520.6
Purity: ≥95%
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

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

Description

trans-carboxy Glimepiride is a metabolite of the sulfonyleurea glimepiride (Item No. 12090).¹ It is formed from glimepiride in a two-step process mediated by the cytochrome P450 (CYP) isoform CYP2C9 and cytosolic enzymes.

Reference

1. Noh, K., Kim, E., Jeong, T.C., *et al.* Simultaneous determination of glimepiride and its metabolites in human plasma by liquid chromatography coupled to a tandem mass spectrometry. *Arch. Pharm. Res.* **34**(12), 2073-2078 (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, 11/18/2022

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