

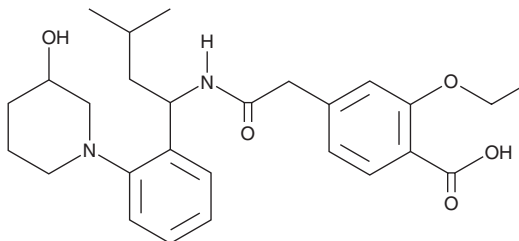
PRODUCT INFORMATION



3-hydroxy Repaglinide

Item No. 45941

CAS Registry No.: 874908-14-2
Formal Name: 2-ethoxy-4-[2-[[1-[2-(3-hydroxy-1-piperidiny)phenyl]-3-methylbutyl]amino]-2-oxoethyl]-benzoic acid
MF: C₂₇H₃₆N₂O₅
FW: 468.6
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

3-hydroxy Repaglinide is supplied as a solid. A stock solution may be made by dissolving the 3-hydroxy repaglinide in the solvent of choice, which should be purged with an inert gas. 3-hydroxy Repaglinide is slightly soluble (0.1-1 mg/ml) in acetonitrile.

Description

3-hydroxy Repaglinide is a previously proposed metabolite of repaglinide (Item No. 19387).¹ It was previously proposed to be a metabolite formed by the cytochrome P450 (CYP) isoform CYP2C8, but it was later shown that CYP2C8 metabolizes repaglinide to 4-hydroxy repaglinide.

Reference

1. Sharma, R., Balesano, A., Doran, A.C., *et al.* Human cytochrome P4502C8 metabolizes repaglinide to 4'-hydroxyrepaglinide, not 3'-hydroxyrepaglinide. *Drug Metab. Dispos.* **53(11)**, 100177 (2025).

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, 06/26/2026

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