

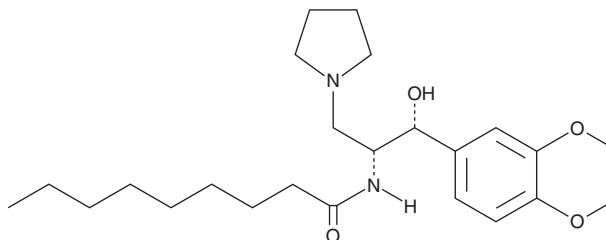
PRODUCT INFORMATION



Genz-123346

Item No. 28500

CAS Registry No.: 491833-30-8
Formal Name: N-[(1R,2R)-2-(2,3-dihydro-1,4-benzodioxin-6-yl)-2-hydroxy-1-(1-pyrrolidinylmethyl)ethyl]-nonanamide
MF: C₂₄H₃₈N₂O₄
FW: 418.6
Purity: ≥98%
Supplied as: A crystalline 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

Genz-123346 is supplied as a crystalline solid. A stock solution may be made by dissolving the genz-123346 in the solvent of choice, which should be purged with an inert gas. Genz-123346 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of genz-123346 in these solvents is approximately 30 mg/ml.

Description

Genz-123346 is a glucosylceramide synthase inhibitor (IC₅₀ = 14 nM).¹ It reduces hepatic levels of glucosylceramide and triglycerides as well as blood levels of alanine aminotransferase (ALT), hemoglobin A1c, and non-fasting glucose in *ob/ob* mice when administered at a dose of 120 mg/kg per day.² Genz-123346 (125 mg/kg per day) also decreases hepatic fat mass and reduces the size and number of hepatic lipid droplets in diet-induced obese mice.

References

1. Zhao, H., Przybylska, M., Wu, I.-H., *et al.* Inhibiting glycosphingolipid synthesis improves glycemic control and insulin sensitivity in animal models of type 2 diabetes. *Diabetes* **56(5)**, 1210-1218 (2007).
2. Zhao, H., Przybylska, M., Wu, I.-H., *et al.* Inhibiting glycosphingolipid synthesis ameliorates hepatic steatosis in obese mice. *Hepatology* **50(1)**, 85-93 (2009).

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, 12/08/2022

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