

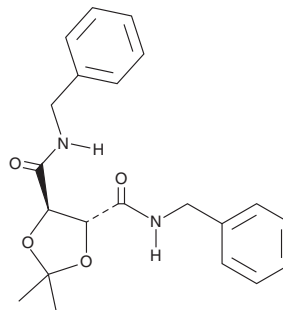
PRODUCT INFORMATION



HOR1-C59

Item No. 45597

CAS Registry No.: 189574-20-7
Formal Name: (4R,5R)-2,2-dimethyl-N⁴,N⁵-bis(phenylmethyl)-1,3-dioxolane-4,5-dicarboxamide
MF: C₂₁H₂₄N₂O₄
FW: 368.4
Purity: ≥98%
Supplied as: A 50 mg/ml solution in ethanol
Storage: -20°C
Stability: ≥4 years



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

Description

HOR1-C59 is an olfactory receptor 5V1 (Or5V1) agonist.¹ It induces G_{α_s}-Gβ_γ dissociation in HEK293 cells overexpressing human Or5V1 (EC₅₀ = 7.12 nM). *In vivo*, HOR1-C59 (50 ng/g) improves glucose tolerance, increases gluconeogenesis, and decreases liver steatosis and epididymal adipocyte size in high-fat diet-fed mice.

Reference

1. Ge, X.Y., Cheng, J., Zhang, L.J., *et al.* Identification of Or5v1/Olfr110 as an oxylipin receptor and anti-obesity target. *Cell*. **189(5)**, 1481-1498 (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, 03/24/2026

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