

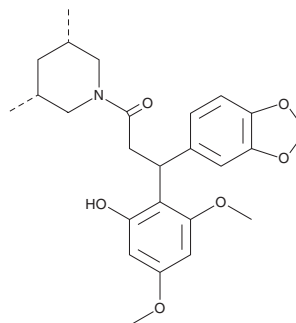
PRODUCT INFORMATION



ML-209

Item No. 30164

CAS Registry No.: 1334526-14-5
Formal Name: *rel*-3-(1,3-benzodioxol-5-yl)-1-[(3*R*,5*S*)-3,5-dimethyl-1-piperidiny]-3-(2-hydroxy-4,6-dimethoxyphenyl)-1-propanone
MF: C₂₅H₃₁NO₆
FW: 441.5
Purity: ≥98%
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

ML-209 is supplied as a solid. A stock solution may be made by dissolving the ML-209 in the solvent of choice, which should be purged with an inert gas. ML-209 is soluble in the organic solvent DMSO at a concentration of approximately 100 mM.

Description

ML-209 is an antagonist of retinoic acid receptor-related orphan receptor γ t (ROR γ t; IC₅₀ = 1.1 μ M in a reporter assay).¹ It inhibits ROR γ t-induced transcription in HEK293T cells expressing the human receptor (IC₅₀ = 300 nM). ML-209 inhibits ROR γ t-dependent differentiation of isolated naïve cord blood human CD4⁺ T cells into Th17 T helper cells.

Reference

1. Huh, J.R., Englund, E.E., Wang, H., *et al.* Identification of potent and selective diphenylpropanamide ROR γ inhibitors. *ACS Med. Chem. Lett.* **4**(1), 79-84 (2013).

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, 12/12/2022

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