

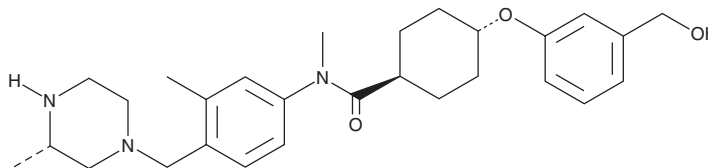
PRODUCT INFORMATION



DS-3801b

Item No. 42548

CAS Registry No.: 1369412-66-7
Formal Name: *trans*-4-[3-(hydroxymethyl)phenoxy]-N-methyl-N-[3-methyl-4-[[[(3S)-3-methyl-1-piperazinyl]methyl]phenyl]-cyclohexanecarboxamide
MF: C₂₈H₃₉N₃O₃
FW: 465.6
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

DS-3801b is supplied as a solid. A stock solution may be made by dissolving the DS-3801b in the solvent of choice, which should be purged with an inert gas. DS-3801b is soluble (≥10 mg/ml) in ethanol and DMSO.

Description

DS-3801b is an agonist of the motilin receptor.¹ It induces the production of inositol monophosphate (IP₁) in COS-7 cells expressing the human or rabbit motilin receptor (EC₅₀s = 5.3 and 1.3 nM, respectively). DS-3801b induces contractions in isolated rabbit duodenal muscle strips (EC₅₀ = 7.3 nM).

Reference

1. Toda, N., Shida, T., Takano, R., *et al.* Discovery of DS-3801b, a non-macrolide GPR38 agonist with N-methylanilide structure. *Bioorg. Med. Chem. Lett.* **59**, 128554 (2022).

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, 02/26/2025

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