

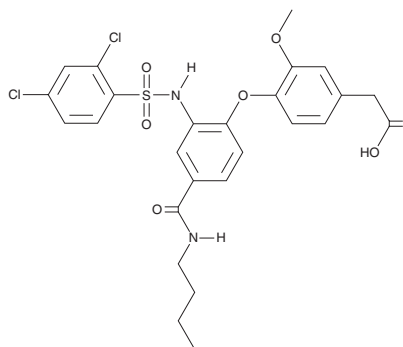
PRODUCT INFORMATION



AMG 009

Item No. 42392

CAS Registry No.: 1027847-67-1
Formal Name: 4-[4-[(butylamino)carbonyl]-2-[[[2,4-dichlorophenyl)sulfonyl]amino]phenoxy]-3-methoxy-benzeneacetic acid
MF: C₂₆H₂₆Cl₂N₂O₇S
FW: 581.5
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

AMG 009 is supplied as a solid. A stock solution may be made by dissolving the AMG 009 in the solvent of choice, which should be purged with an inert gas. AMG 009 is sparingly soluble (1-10 mg/ml) in ethanol and DMSO.

Description

AMG 009 is a dual antagonist of DP₂/CRTH₂ and DP₁ (IC₅₀s = 3 and 12 nM, respectively).¹ It decreases airway resistance in a guinea pig model of airflow obstruction induced by aerosolized prostaglandin D₂ (PGD₂; Item No. 12010) when administered at doses of 3, 10, and 30 mg/kg.

Reference

1. Liu, J., Fu, Z., Wang, Y., *et al.* Discovery and optimization of CRTH2 and DP dual antagonists. *Bioorg. Med. Chem. Lett.* **19(22)**, 6419-6423 (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, 12/11/2024

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