

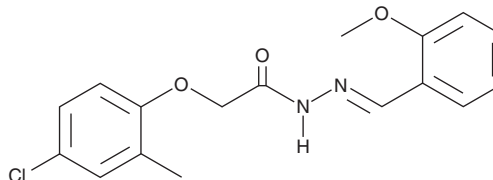
PRODUCT INFORMATION



Ani 9

Item No. 29733

CAS Registry No.: 356102-14-2
Formal Name: 2-(4-chloro-2-methylphenoxy)-acetic acid, 2-[(2-methoxyphenyl)methylene]hydrazide
MF: C₁₇H₁₇ClN₂O₃
FW: 332.8
Purity: ≥98%
Supplied as: A solid
Storage: -20°C
Stability: ≥2 years



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

Laboratory Procedures

Ani 9 is supplied as a solid. A stock solution may be made by dissolving the ani 9 in the solvent of choice, which should be purged with an inert gas. Ani 9 is soluble in the organic solvent DMSO at a concentration of approximately 25 mg/ml.

Description

Ani 9 is an inhibitor of the calcium-activated chloride channel (CaCC) anoctamin 1 (ANO1) that blocks ATP-induced apical membrane chloride currents in ANO1-expressing Fischer rat thyroid (FRT) cells (IC₅₀ = 77 nM).¹ It is selective for ANO1 over ANO2 at 10 μM and the cystic fibrosis transmembrane conductance regulator (CFTR) and epithelial sodium channel (ENaC) at 30 μM. Ani 9 inhibits ATP-induced quenching of a fluorescent intracellular halide sensor in PC3 prostate cancer cells and Capan-1 pancreatic carcinoma cells in a concentration-dependent manner.

Reference

1. Seo, Y., Lee, H.K., Park, J., *et al.* Ani9, A novel potent small-molecule ANO1 inhibitor with negligible effect on ANO2. *PLoS One* **11(5)**, e0155771 (2016).

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, 01/30/2020

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