

PRODUCT INFORMATION



SQ-609

Item No. 41875

CAS Registry No.: 627052-25-9

Formal Name: 1-[[1-(tricyclo[3.3.1.1^{3,7}]dec-1-ylmethyl)-4-piperidinyl]methyl]-4-piperidinol

MF: C₂₂H₃₈N₂O

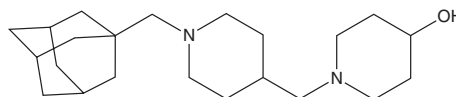
FW: 346.6

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

SQ-609 is supplied as a solid. A stock solution may be made by dissolving the SQ-609 in the solvent of choice, which should be purged with an inert gas. SQ-609 is sparingly soluble (1-10 mg/ml) in ethanol.

Description

SQ-609 is an antibacterial agent.¹ It reduces the proliferation of *M. tuberculosis* in *M. tuberculosis*-infected J774 macrophages (MIC = 4 µg/ml). *In vivo*, SQ-609 (10 mg/kg per day) increases survival time in *M. tuberculosis*-infected female mice.

Reference

1. Bogatcheva, E., Hanrahan, C., Nikonenko, B., *et al.* Identification of SQ609 as a lead compound from a library of dipiperidines. *Bioorg. Med. Chem. Lett.* **21(18)**, 5353-5357 (2011).

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, 10/15/2024

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