

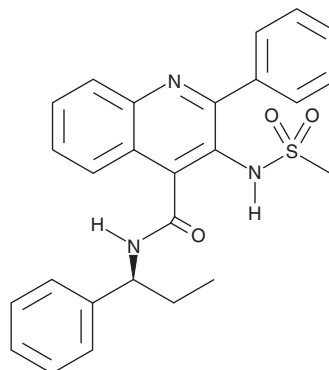
PRODUCT INFORMATION



MLE-4901

Item No. 23587

CAS Registry No.: 941690-55-7
Formal Name: 3-[(methylsulfonyl)amino]-2-phenyl-N-[(1S)-1-phenylpropyl]-4-quinolinecarboxamide
Synonyms: AZD 2624, AZD 4901
MF: C₂₆H₂₅N₃O₃S
FW: 459.6
Purity: ≥98%
UV/Vis.: λ_{max}: 236 nm
Supplied as: A crystalline 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

MLE-4901 is supplied as a crystalline solid. A stock solution may be made by dissolving the MLE-4901 in the solvent of choice, which should be purged with an inert gas. MLE-4901 is soluble in organic solvents such as DMSO and dimethyl formamide (DMF). The solubility of MLE-4901 in these solvents is approximately 20 and 30 mg/ml, respectively.

MLE-4901 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, MLE-4901 should first be dissolved in DMF and then diluted with the aqueous buffer of choice. MLE-4901 has a solubility of approximately 0.25 mg/ml in a 1:3 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

MLE-4901 is an antagonist of the neurokinin-3 (NK₃) receptor.¹ It inhibits human liver microsomal CYP3A4/5 with apparent IC₅₀ values of 7.1 and 19.8 μM in midazolam 1'-hydroxylation and testosterone 6β-hydroxylation assays, respectively.

Reference

1. Li, Y., Zhou, D., Ferguson, S.S., *et al.* *In vitro* assessment of metabolic drug-drug interaction potential of AZD2624, neurokinin-3 receptor antagonist, through cytochrome P450 enzyme identification, inhibition, and induction studies. *Xenobiotica* **40(11)**, 721-729 (2010).

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, 11/08/2022

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