

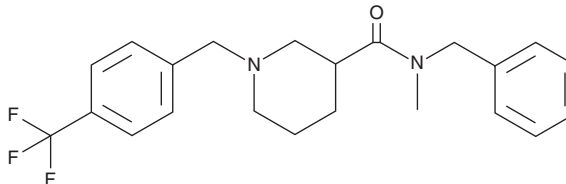
PRODUCT INFORMATION



MDK1088

Item No. 37394

CAS Registry No.: 1350920-22-7
Formal Name: N-methyl-N-(phenylmethyl)-1-[[4-(trifluoromethyl)phenyl]methyl]-3-piperidinecarboxamide
MF: C₂₂H₂₅F₃N₂O
FW: 390.4
Purity: ≥98%
Supplied as: A neat oil
Storage: -20°C
Stability: ≥4 years
Item Origin: Synthetic



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

Laboratory Procedures

MDK1088 is supplied as a neat oil. A stock solution may be made by dissolving the MDK1088 in the solvent of choice, which should be purged with an inert gas. MDK1088 is soluble (≥10 mg/ml) in DMSO and ethanol.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of MDK1088 can be prepared by directly dissolving the neat oil in aqueous buffers. MDK1088 is soluble (≥10 mg/ml) in PBS (pH 7.2). We do not recommend storing the aqueous solution for more than one day.

Description

MDK1088 is an antiprotozoal agent.¹ It inhibits the replication of *T. cruzi* in NIH3T3 cells (IC₅₀ = 0.008 μM).

Reference

1. Germain, A.R., Carmody, L.C., Dockendorff, C., *et al.* Identification of small-molecule inhibitors of *Trypanosoma cruzi* replication. *Bioorg. Med. Chem. Lett.* **21**(23), 7197-7200 (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, 02/17/2025

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