

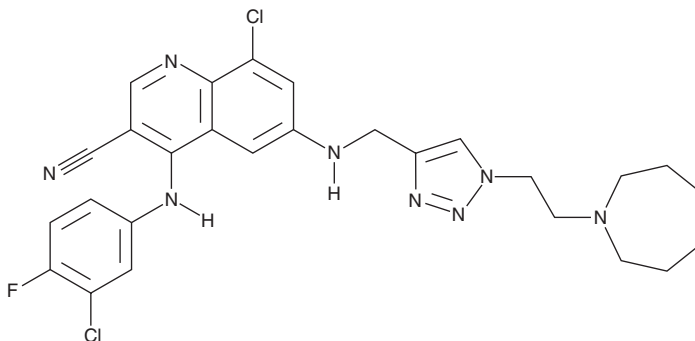
PRODUCT INFORMATION



Cot Inhibitor 1

Item No. 43368

CAS Registry No.: 915365-57-0
Formal Name: 8-chloro-4-[(3-chloro-4-fluorophenyl)amino]-6-[[[1-[2-(hexahydro-1H-azepin-1-yl)ethyl]-1H-1,2,3-triazol-4-yl)methyl]amino]-3-quinolinecarbonitrile
MF: C₂₇H₂₇Cl₂FN₈
FW: 553.5
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

Cot inhibitor 1 is supplied as a solid. A stock solution may be made by dissolving the cot inhibitor 1 in the solvent of choice, which should be purged with an inert gas. Cot inhibitor 1 is soluble (≥10 mg/ml) in DMSO.

Description

Cot inhibitor 1 is an inhibitor of MAP3K8 (IC₅₀ = 28 nM).¹ It inhibits LPS-induced increases in TNF-α levels in isolated human whole blood (IC₅₀ = 5.4 μM).

Reference

1. Wu, J., Green, N., Hotchandani, R., *et al.* Selective inhibitors of tumor progression loci-2 (Tpl2) kinase with potent inhibition of TNF-α production in human whole blood. *Bioorg. Med. Chem. Lett.* **19(13)**, 3485-3488 (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, 04/29/2025

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