

PRODUCT INFORMATION

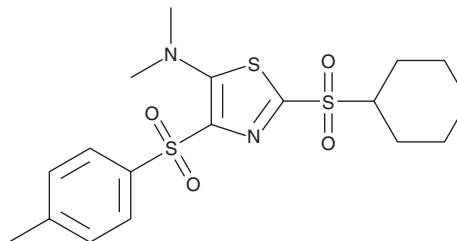


M04

Item No. 35454

CAS Registry No.: 875158-73-9
Formal Name: 2-(cyclohexylsulfonyl)-N,N-dimethyl-4-[(4-methylphenyl)sulfonyl]-5-thiazolamine

MF: C₁₈H₂₄N₂O₄S₃
FW: 428.6
Purity: ≥98%
UV/Vis.: λ_{max}: 230, 328 nm
Supplied as: A crystalline 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

M04 is supplied as a crystalline solid. A stock solution may be made by dissolving the M04 in the solvent of choice, which should be purged with an inert gas. M04 is soluble in chloroform.

Description

M04 is an agonist of stimulator of interferon genes (STING).¹ It induces IFN reporter gene expression in HEK293T cells expressing wild-type human STING but not in HEK293T cells expressing the R71H-G230A-R293Q (HAQ) human STING variant nor in mouse RAW 264.7 cells, indicating allelic- and species-dependent activities, at 75 μM. M04 induces the production of TNF-α, IL-10, IL-1β, and IL-12p70 in human peripheral blood mononuclear cells (PBMCs). It increases the percentage of isolated human monocyte-derived dendritic cells expressing the MHC class II cell surface receptor HLA-DR and the co-stimulatory molecules CD40, CD80, and CD86, as well as increases T cell cross-priming of the same cells in an *ex vivo* assay, when used at a concentration of 50 μM.

Reference

1. Abraham, J., Botto, S., Mizuno, N., *et al.* Characterization of a novel compound that stimulates STING-mediated innate immune activity in an allele-specific manner. *Front. Immunol.* **11**, 1430 (2020).

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, 12/09/2021

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