

PRODUCT INFORMATION

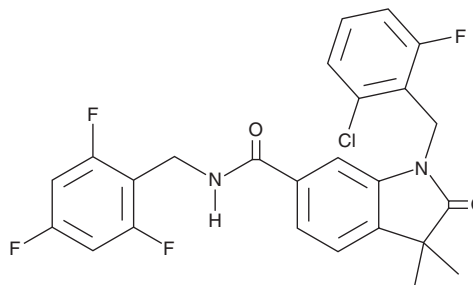


C53

Item No. 37354

CAS Registry No.: 2259624-71-8
Formal Name: 1-[(2-chloro-6-fluorophenyl)methyl]-2,3-dihydro-3,3-dimethyl-2-oxo-N-[(2,4,6-trifluorophenyl)methyl]-1H-indole-6-carboxamide

Synonym: STING Agonist-12
MF: C₂₅H₁₉ClF₄N₂O₂
FW: 490.9
Purity: ≥98%
UV/Vis.: λ_{max}: 219 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

C53 is supplied as a crystalline solid. A stock solution may be made by dissolving the C53 in the solvent of choice, which should be purged with an inert gas. C53 is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of C53 in these solvents is approximately 2 and 1 mg/ml, respectively.

Description

C53 is a STING agonist.¹ It activates human wild-type STING (EC₅₀ = 185 nM), as well as STING containing the polymorphisms R232, H232, or HAQ in HEK293 cells in an interferon-stimulated gene 54 (ISG54) luciferase reporter assay when used at a concentration of 10 μM. C53 (10 μM) induces the expression of *IFN-α*, *IFN-β*, *TNF-α*, chemokine (C-X-C motif) ligand 10 (*CXCL10*), *IL-12p40*, and *IL-6* in isolated human monocyte-derived dendritic cells (hMO-DCs). It also increases the expression of the dendritic cell markers *CD83* and *CD86* in hMO-DCs when used at a concentration of 10 μM.

Reference

1. Pryde, D.C., Middya, S., Banerjee, M., *et al.* The discovery of potent small molecule activators of human STING. *Eur. J. Med. Chem.* **209**, 112869 (2021).

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, 08/05/2022

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