

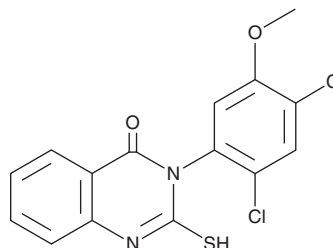
PRODUCT INFORMATION



Mdivi 1

Item No. 15559

CAS Registry No.: 338967-87-6
Formal Name: 3-(2,4-dichloro-5-methoxyphenyl)-2,3-dihydro-2-thioxo-4(1H)-quinazolinone
Synonym: Mitochondrial Division Inhibitor 1
MF: C₁₅H₁₀Cl₂N₂O₂S
FW: 353.2
Purity: ≥98%
UV/Vis.: λ_{max}: 223, 295 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

Mdivi 1 is supplied as a crystalline solid. A stock solution may be made by dissolving the mdivi 1 in the solvent of choice, which should be purged with an inert gas. Mdivi 1 is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of mdivi 1 in these solvents is approximately 20 mg/ml.

Mdivi 1 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, mdivi 1 should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. Mdivi 1 has a solubility of approximately 0.1 mg/ml in a 1:6 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Mdivi 1 is a mitochondrial division inhibitor.^{1,2} It inhibits the GTPase activity of yeast, but not human, dynamin-1 (Dnm1; IC₅₀ = ~1-10 μM for the recombinant yeast enzyme) and human dynamin-related protein 1 (DRP1) in A549 lung cancer cells when used at a concentration of 50 μM. Mdivi 1 (50 μM) inhibits mitochondrial fragmentation and DRP1 self-assembly induced by staurosporine (Item No. 81590) in COS cells overexpressing DRP1.¹ It induces apoptosis in, and inhibits colony formation of, A549 cells when used at concentrations of 25 and 50 μM, respectively.² Mdivi 1 (1 μM) decreases doxorubicin-induced increases in infarct-to-risk ratios in a Langendorff isolated perfused rat heart model of ischemia-reperfusion injury.³ It reduces the severity of lung injury in a mouse model of LPS-induced acute lung injury when administered at a dose of 20 mg/kg.⁴

References

1. Cassidy-Stone, A., Chipuk, J.E., Ingeman, E., *et al.* Chemical inhibition of the mitochondrial division dynamin reveals its role in Bax/Bak-dependent mitochondrial outer membrane permeabilization. *Dev. Cell* **14**(2), 193-204 (2008).
2. Wu, D., Dasgupta, A., Chen, K.-H., *et al.* Identification of novel dynamin-related protein 1 (Drp1) GTPase inhibitors: Therapeutic potential of Drpitor1 and Drpitor1a in cancer and cardiac ischemia-reperfusion injury. *FASEB J.* **34**(1), 1447-1464 (2020).
3. Gharanei, M., Hussain, A., Janneh, O., *et al.* Attenuation of doxorubicin-induced cardiotoxicity by mdivi-1: A mitochondrial division/mitophagy inhibitor. *PLoS One* **8**(10), e77713 (2013).
4. Deng, S., Zhang, L., Mo, Y., *et al.* Mdivi-1 attenuates lipopolysaccharide-induced acute lung injury by inhibiting MAPKs, oxidative stress and apoptosis. *Pulm. Pharmacol. Ther.* **62**, 101918 (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, 02/01/2024

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