

Mdivi-1

Inhibitor of mitochondrial division

Mdivi-1 is a selective inhibitor of mitochondrial division in yeast and mammalian cells which acts via inhibiting the mitochondrial division dynamin. In cells, Mdivi-1 inhibits apoptosis by inhibiting mitochondrial outer membrane permeabilization. Mdivi-1 causes the rapid (<5 min) reversible and dose-dependent formation of net-like mitochondria in wild-type cells with an $IC_{50} = \sim 10\mu M$. In yeast, time-lapse fluorescence microscopy revealed no detectable mitochondrial division after treatment with Mdivi-1. It inhibits Dnm1 GTPase activity in a dose dependent manner with an estimated IC_{50} of 1-10 μM but is not a general inhibitor of GTPases.

Citations: 21

[View Online »](#)

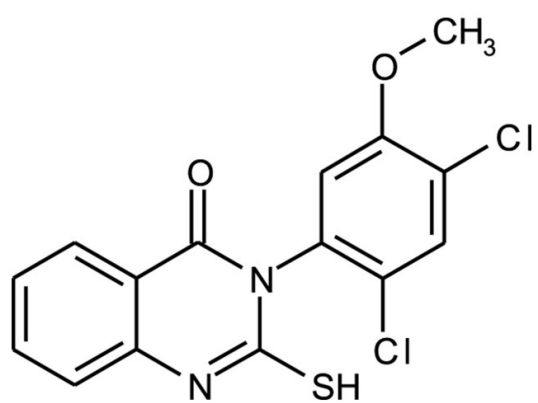
Ordering Information

[Order Online »](#)

BML-CM127-0010	10mg
BML-CM127-0050	50mg

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability As indicated on product label or CoA when stored as recommended.

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Appearance White solid.

CAS 338967-87-6

Couple Target Dynamin, MPTP

Couple Type Inhibitor

Formula $C_{15}H_{10}Cl_2N_2O_2S$

Identity Determined by 1H -NMR.

MW 353.2

Purity $\geq 97\%$ (HPLC)

Solubility Soluble in DMSO (25mg/ml).

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
[info-
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
[info-
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)