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} = ~10µM. 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 1-10µM but is not a general inhibitor of GTPases.

Citations: 21

View Online »

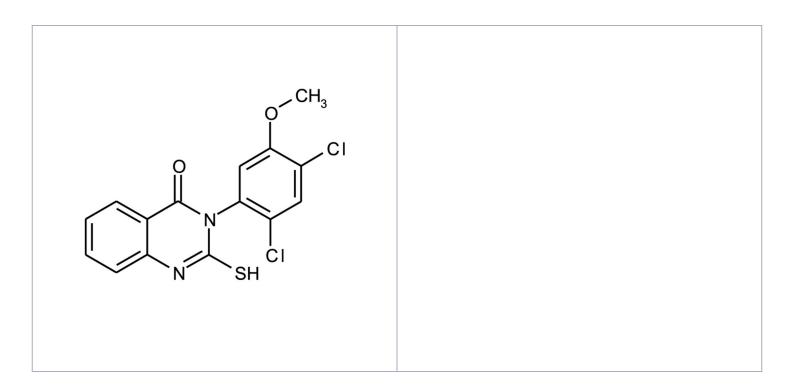
Ordering Information

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 $C_{15}H_{10}CI_2N_2O_2S$

Identity Determined by 1H-NMR.

MW 353.2

Purity ≥97% (HPLC)

Soluble in DMSO (25mg/ml).

Last modified: May 29, 2024