

CGP-37157

MPTP inhibitor

Potent, selective inhibitor of mitochondrial $\text{Na}^+\text{Ca}^{2+}$ exchange. IC_{50} = 0.4 μM . Blocks mitochondrial permeability transition.

Citations: 11

[View Online »](#)

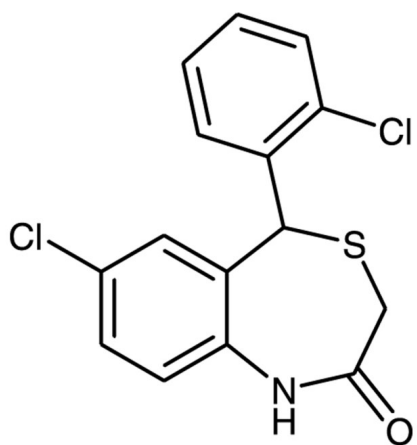
Ordering Information

[Order Online »](#)

BML-CM119-0005	5mg
----------------	-----

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

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

Long Term Storage Ambient

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Appearance	White solid.
CAS	75450-34-9
Couple Target	MPTP
Couple Type	Inhibitor
Formula	$C_{15}H_{11}Cl_2NOS$
MW	324.2
Purity	≥98% (TLC)
Solubility	Soluble in DMSO.

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

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

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

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