

FCCP

MPTP inhibitor

A potent reversible inhibitor of mitochondrial oxidative phosphorylation. FCCP is a useful tool for depolarizing mitochondrial membrane potential. Treatment of cells with FCCP at varying concentrations leads to partial (100 nM) or complete (10 μ M) depolarization and apoptosis.

Citations: 19

[View Online »](#)

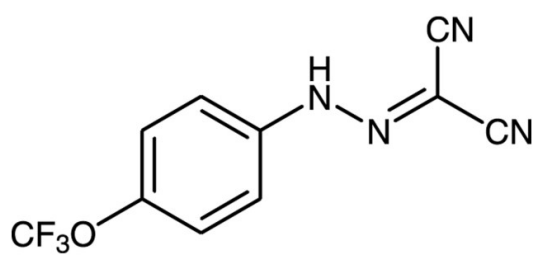
Ordering Information

[Order Online »](#)

BML-CM120-0010	10mg
BML-CM120-0050	50mg

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

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

Long Term Storage Ambient

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Carbonylcyanide-4-(trifluoromethoxy)-phenylhydrazine

Appearance Yellow solid.

CAS 370-86-5

Couple Target MPTP

Couple Type Inhibitor

Formula $C_{10}H_5F_3N_4O$

Identity Determined by ¹H-NMR.

MW 254.2

Purity ≥98% (HPLC)

Solubility Soluble in DMSO

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)