

Metyrapone

Cytochrome P450 inhibitor

Inhibits cytochrome P-450-mediated prostaglandin ω/ω -1 hydroxylase activity.

Citations: 1

[View Online »](#)

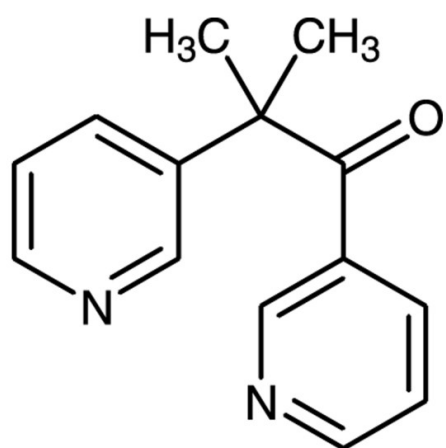
Ordering Information

[Order Online »](#)

BML-EI256-0200	200mg
----------------	-------

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 54-36-4

Couple Target Cytochrome P450

Couple Type Inhibitor

Formula $C_{14}H_{14}N_2O$

MW 226.34

Purity $\geq 99\%$

Solubility Soluble in water (25 mg/ml), DMSO (25 mg/ml) or ethanol (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)