

Ketoconazole

Cytochrome P450 inhibitor

Inhibits cytochrome P-450-dependent steps in the biosynthesis of steroid hormones *in vivo*. Antimetastatic and antineoplastic activity. Antipsoriatic activity. Orally active 5-lipoxygenase and thromboxane synthase inhibitor.

Citations: 2

[View Online »](#)

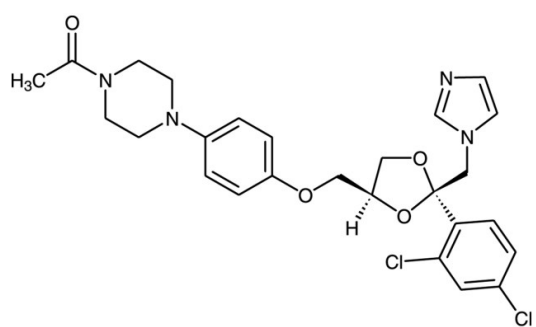
Ordering Information

[Order Online »](#)

BML-EI107-0050	50mg
BML-EI107-0500	500mg

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

CAS 65277-42-1

Couple Target Cytochrome P450, Lipoxygenase, Thromboxane synthase

Couple Type Inhibitor

MW 531.4

Purity ≥99% (HPLC)

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-
be@enzolifesciences.com](mailto:info-be@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)