

Finasteride

5 α -reductase inhibitor

Inhibitor of 5 α -reductase, the enzyme which converts testosterone to the more potent androgen, 5 α -dihydrotestosterone. Used as a treatment in benign prostatic hyperplasia and prostate cancer. Also used to treat male pattern hair loss.

Citations: 8

[View Online »](#)

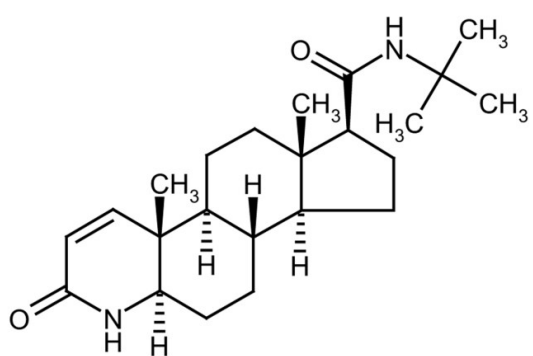
Ordering Information

[Order Online »](#)

ALX-270-491-M100	100mg
ALX-270-491-M500	500mg

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

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

Long Term Storage +4°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name MK-906

Appearance White to off-white crystalline powder.

CAS 98319-26-7

Couple Target Reductase

Couple Type Inhibitor

Formula $C_{23}H_{36}N_2O_2$

Identity Determined by IR.

MI 14: 4082

MW 372.5

Purity ≥98% (Assay (UV))

RTECS CL5245000

Solubility Soluble in DMSO, 100% ethanol, methanol or propanol.

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)