

Methotrexate

Dihydrofolate reductase inhibitor

Folic acid antagonist. Potent inhibitor of dihydrofolate reductase. Induces apoptosis. Inhibits DNA synthesis. Anticancer agent.

Citations: 12

[View Online »](#)

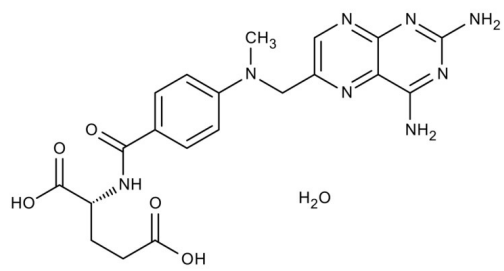
Ordering Information

[Order Online »](#)

ALX-440-045-M050	50mg
ALX-440-045-G001	1g

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended.
Handling	Protect from light.
Long Term Storage	+4°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	(+)-Amethopterin, 4-Amino-10-methylfolic acid, Methylaminopterin, MTX
Appearance	Yellow powder.
CAS	133073-73-1
Couple Target	Dihydrofolate reductase
Couple Type	Inhibitor
Formula	$C_{20}H_{22}N_8O_5 \cdot H_2O$
MI	14: 5985
MW	454.5
Purity	≥98% (HPLC)
Solubility	Soluble in alkaline solution; insoluble in water, 100% ethanol or ether.

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