

# Methotrexate

Dihydrofolate reductase inhibitor

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

Citations: 12

[View Online »](#)

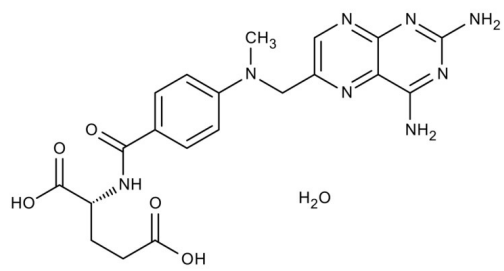
## Ordering Information

[Order Online »](#)

ALX-440-045-M050	50mg
ALX-440-045-M100	100mg
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](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)