

Cyclophosphamide . monohydrate

DNA alkylating agent

Alkylates DNA, forming DNA-DNA cross-links that result in inhibition of DNA synthesis and cell death. Interestingly the repair enzyme O⁶-alkylguanine-DNA alkyltransferase, long thought to be involved with resistance to methylating and chloroethylating agents, may also contribute to resistance to the cytotoxic and mutagenic effects of cyclophosphamide.

Citations: 6

[View Online »](#)

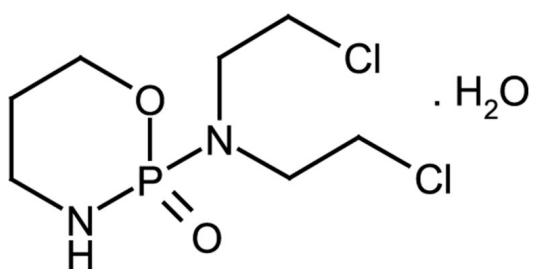
Ordering Information

[Order Online »](#)

ALX-400-051-G001	1g
ALX-400-051-G005	5g

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

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

Long Term Storage -20°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name 2-[Bis(2-chloroethyl)amino]tetrahydro-2H-1,3,2-oxazaphosphorine-2-oxide, Cytosan

Appearance White to off-white crystalline solid.

CAS 6055-19-2

Formula $C_7H_{15}Cl_2N_2O_2P \cdot H_2O$

Identity Determined by IR and UV.

MI 14: 2747

MW 261.1 . 18.0

Purity ≥97% (HPLC)

RTECS RP6157750

Solubility Soluble in water; slightly soluble in 100% ethanol.

Source Synthetic.



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)