

# 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 O<sup>6</sup>-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: 7

[View Online »](#)

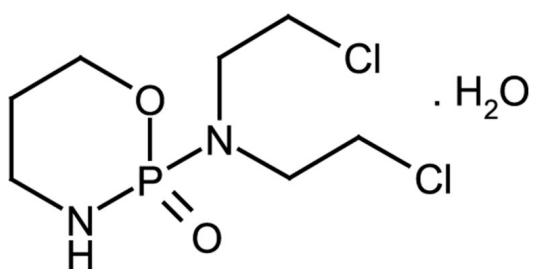
## Ordering Information

[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.

**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-  
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

Germany  
Phone: +49 7621 5500 526  
[info-  
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)

UK & Ireland  
Phone (UK customers):  
0845 601 1488  
Phone: +44 1392 825900  
[info-  
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)