# 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  $\rm O^6$ -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 »

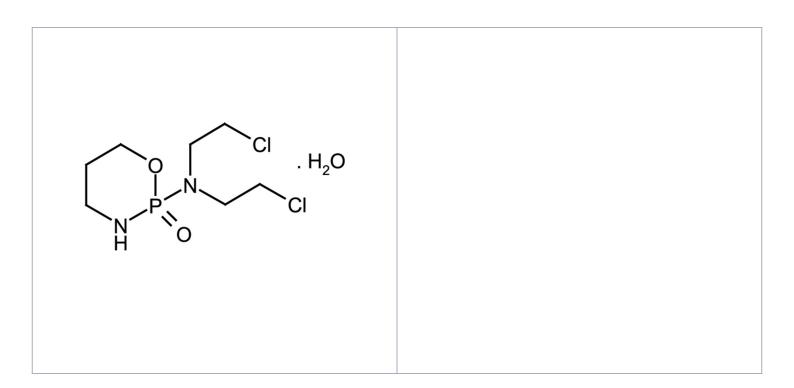
## **Ordering Information**

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, Cytoxan

**Appearance** White to off-white crystalline solid.

**CAS** 6055-19-2

Formula  $C_7H_{15}CI_2N_2O_2P \cdot H_2O$ 

**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

