

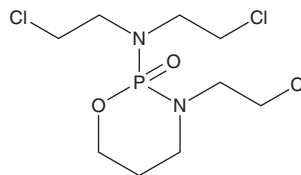
# PRODUCT INFORMATION



## Trofosfamide

Item No. 21874

**CAS Registry No.:** 22089-22-1  
**Formal Name:** N,N,3-tris(2-chloroethyl)tetrahydro-2H-1,3,2-oxazaphosphorin-2-amine, 2-oxide  
**Synonym:** NSC 109723  
**MF:** C<sub>9</sub>H<sub>18</sub>Cl<sub>3</sub>N<sub>2</sub>O<sub>2</sub>P  
**FW:** 323.6  
**Purity:** ≥95%  
**Supplied as:** A crystalline solid  
**Storage:** -20°C  
**Stability:** ≥4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

### Laboratory Procedures

Trofosfamide is supplied as a crystalline solid. A stock solution may be made by dissolving the trofosfamide in the solvent of choice, which should be purged with an inert gas. Trofosfamide is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of trofosfamide in ethanol and DMF is approximately 50 mg/ml and approximately 30 mg/ml in DMSO.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of trofosfamide can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of trofosfamide in PBS, pH 7.2, is approximately 10 mg/ml. We do not recommend storing the aqueous solution for more than one day.

### Description

Trofosfamide is a nitrogen mustard alkylating agent and a derivative of cyclophosphamide (Item No. 13849) that has antiproliferative activities.<sup>1,2</sup> It increases the number of micronucleated polychromatic red blood cells isolated from mouse bone marrow, indicating cytogenetic activity, when administered at doses of 40, 80, and 160 mg/kg.<sup>3</sup> Trofosfamide (100 mg/kg) inhibits proliferation of isolated splenocytes in the allogenic mixed lymphocyte reaction.<sup>1</sup> It reduces sheep red blood cell-induced increases in the number of antibody-producing spleen cells isolated from rats when administered at a dose of 40 mg/kg.<sup>2</sup>

### References

1. Devlin, R.G., Schwartz, N.L., Baronowsky, P.E., *et al.* Inhibition of cellular immune reactions by cyclophosphamide analogues ifosfamide and trofosfamide. I. Mixed lymphocyte reaction. *Proc. Soc. Exp. Biol. Med.* **145(2)**, 389-391 (1974).
2. Harrison, E.F. and Fuquay, M.E. Immunosuppressive properties of cyclophosphamide analogues. *Proc. Soc. Exp. Biol. Med.* **139(3)**, 957-963 (1972).
3. Wild, D. Cytogenetic effects in the mouse of 17 chemical mutagens and carcinogens evaluated by the micronucleus test. *Mutat. Res.* **56(3)**, 319-327 (1978).

#### WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

#### SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the complete Safety Data Sheet, which has been sent via email to your institution.

#### WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 10/10/2022

#### CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD  
ANN ARBOR, MI 48108 · USA

**PHONE:** [800] 364-9897  
[734] 971-3335

**FAX:** [734] 971-3640

CUSTSERV@CAYMANCHEM.COM  
WWW.CAYMANCHEM.COM