

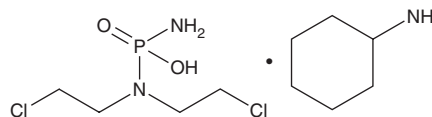
PRODUCT INFORMATION



Phosphoramidate Mustard (cyclohexylammonium salt)

Item No. 34078

CAS Registry No.: 1566-15-0
Formal Name: N,N-bis(2-chloroethyl)-phosphorodiamidic acid, compd. with cyclohexanamine
Synonyms: NSC 69945, PMC
MF: C₄H₁₁Cl₂N₂O₂P • C₆H₁₃N
FW: 320.2
Purity: ≥95%
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years
Item Origin: Synthetic



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

Laboratory Procedures

Phosphoramidate mustard (cyclohexylammonium salt) is supplied as a solid. A stock solution may be made by dissolving the phosphoramidate mustard (cyclohexylammonium salt) in the solvent of choice, which should be purged with an inert gas. Phosphoramidate mustard (cyclohexylammonium salt) is slightly soluble in chloroform and methanol.

Description

Phosphoramidate mustard is an alkylating agent and active metabolite of cyclophosphamide (Item No. 13849).¹ It is formed from cyclophosphamide *via* the ring-opened tautomer of the cytochrome P450 (CYP) isoform-formed intermediate 4-hydroxycyclophosphamide.³ Phosphoramidate mustard induces DNA crosslinking, alkylates guanine in DNA, and increases the production of covalent DNA-protein conjugates in, and is cytotoxic to, HT-1080 human fibrosarcoma cells in a concentration-dependent manner. It is toxic to adult mice and teratogenic to embryos when administered to pregnant dams at a dose of 154 mg/kg on day 11 of gestation.²

References

1. Groehler, A., IV, Villalta, P.W., Campbell, C., *et al.* Covalent DNA-protein cross-linking by phosphoramidate mustard and nornitrogen mustard in human cells. *Chem. Res. Toxicol.* **29(2)**, 190-202 (2016).
2. Gibson, J.E. and Becker, B.A. Teratogenicity of structural truncates of cyclophosphamide in mice. *Teratology* **4(2)**, 141-150 (1971).
3. Huitema, A.D., Mathôt, R.A., Tibben, M.M., *et al.* A mechanism-based pharmacokinetic model for the cytochrome P450 drug-drug interaction between cyclophosphamide and thioTEPA and the autoinduction of cyclophosphamide. *J. Pharmacokinetic. Pharmacodyn.* **28(3)**, 211-230 (2001).

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, 11/07/2022

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