

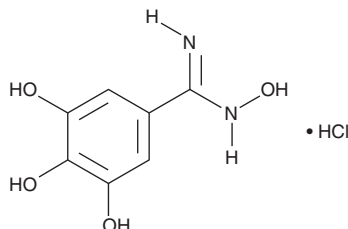
PRODUCT INFORMATION



Trimidox (hydrochloride)

Item No. 10011124

CAS Registry No.: 95933-75-8
Formal Name: N,3,4,5-tetrahydroxybenzenecarboximidamide, monohydrochloride
Synonym: CF 233
MF: C₇H₈N₂O₄ • HCl
FW: 220.6
Purity: ≥98%
UV/Vis.: λ_{max}: 220, 280 nm
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

Trimidox (hydrochloride) is supplied as a crystalline solid. A stock solution may be made by dissolving the trimidox (hydrochloride) in the solvent of choice, which should be purged with an inert gas. Trimidox (hydrochloride) is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of trimidox (hydrochloride) in these solvents is approximately 0.25, 2, and 1 mg/ml, respectively.

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 trimidox (hydrochloride) can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of trimidox (hydrochloride) in PBS (pH 7.2) is approximately 5 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Ribonucleotide reductase, the rate-limiting enzyme for *de novo* DNA synthesis, is a common target for chemotherapy. Its increased activity in cancer cells is associated with malignant transformation and proliferation.¹ Trimidox is a specific ribonucleotide reductase inhibitor that reduces levels of dGTP and dCTP in HL-60 cells, inducing apoptosis *via* activation of caspases without altering the cell cycle distribution.^{1,2} Trimidox inhibits growth of human promyelocytic leukemia HL-60 cells with an IC₅₀ value of 35 μM.¹

References

1. Szekeres, T., Fritzer, M., Strobl, H., *et al.* Synergistic growth inhibitory and differentiating effects of trimidox and tiazofurin in human promyelocytic leukemia HL-60 cells. *Blood* **84**(12), 4316-4321 (1994).
2. Kanno, S.I., Uwai, K., Tomizawa, A., *et al.* Trimidox induces apoptosis via cytochrome c release in NALM-6 human B cell leukaemia cells. *Basic Clin. Pharmacol. Toxicol.* **98**(1), 44-50 (2006).

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, 01/31/2025

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