

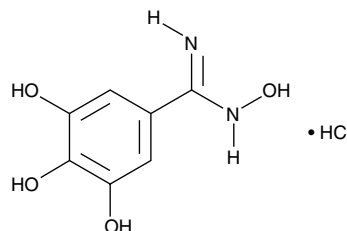
# Product Information



## Trimidox (hydrochloride)

Item No. 10011124

**CAS Registry No.:** 95933-75-8  
**Formal Name:** N,3,4,5-tetrahydroxybenzenecarboximidamide, monohydrochloride  
**Synonym:** VF 233 (hydrochloride)  
**MF:** C<sub>7</sub>H<sub>8</sub>N<sub>2</sub>O<sub>4</sub> • HCl  
**FW:** 220.6  
**Purity:** ≥98%  
**Stability:** ≥2 years at -20°C  
**Supplied as:** A crystalline solid  
**UV/Vis.:** λ<sub>max</sub>: 220, 280 nm



### Laboratory Procedures

For long term storage, we suggest that trimidox be stored as supplied at -20°C. It will be stable for at least two years.

Trimidox is supplied as a crystalline solid. A stock solution may be made by dissolving the trimidox in an organic solvent purged with an inert gas. Trimidox is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of trimidox in these solvents is approximately 20 mg/ml.

Trimidox is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, trimidox should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. Trimidox has a solubility of approximately 0.3 mg/ml in a 1:2 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

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.<sup>1</sup> 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.<sup>1,2</sup> Trimidox inhibits growth of human promyelocytic leukemia HL-60 cells with an IC<sub>50</sub> value of 35 μM.<sup>1</sup>

### 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 & Clinical Pharmacology & Toxicology* **98**, 44-50 (2006).

### Related Products

For a list of related products please visit: [www.caymanchem.com/catalog/10011124](http://www.caymanchem.com/catalog/10011124)

**WARNING: THIS PRODUCT IS FOR LABORATORY RESEARCH ONLY: NOT FOR ADMINISTRATION TO HUMANS. NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.**

#### SAFETY DATA

This material should be considered hazardous until information to the contrary becomes available. Do not ingest, swallow, or inhale. Do not get in eyes, on skin, or on clothing. Wash thoroughly after handling. This information contains some, but not all, of the information required for the safe and proper use of this material. 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

Cayman Chemical Company makes **no warranty or guarantee** of any kind, whether written or oral, expressed or implied, including without limitation, any warranty of fitness for a particular purpose, suitability and merchantability, which extends beyond the description of the chemicals hereof. Cayman **warrants only** to the original customer that the material will meet our specifications at the time of delivery.

Cayman will carry out its delivery obligations with due care and skill. Thus, in no event will Cayman have **any obligation or liability**, whether in tort (including negligence) or in contract, for any direct, indirect, incidental or consequential damages, even if Cayman is informed about their possible existence.

This limitation of liability does not apply in the case of intentional acts or negligence of Cayman, its directors or its employees.

Buyer's **exclusive remedy** and Cayman's sole liability hereunder shall be limited to a refund of the purchase price, or at Cayman's option, the replacement, at no cost to Buyer, of all material that does not meet our specifications.

Said refund or replacement is conditioned on Buyer giving written notice to Cayman within thirty (30) days after arrival of the material at its destination. Failure of Buyer to give said notice within thirty (30) days shall constitute a waiver by Buyer of all claims hereunder with respect to said material.

For further details, please refer to our **Warranty and Limitation of Remedy** located on our website and in our catalog.

Copyright Cayman Chemical Company, 05/06/2013

### Cayman Chemical

#### Mailing address

1180 E. Ellsworth Road  
Ann Arbor, MI  
48108 USA

#### Phone

(800) 364-9897  
(734) 971-3335

#### Fax

(734) 971-3640

#### E-Mail

[custserv@caymanchem.com](mailto:custserv@caymanchem.com)

#### Web

[www.caymanchem.com](http://www.caymanchem.com)