

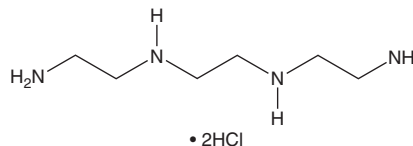
# PRODUCT INFORMATION



## Trientine (hydrochloride)

Item No. 24000

**CAS Registry No.:** 38260-01-4  
**Formal Name:** N<sup>1</sup>,N<sup>2</sup>-bis(2-aminoethyl)-1,2-ethanediamine, dihydrochloride  
**Synonym:** Triethylenetetramine  
**MF:** C<sub>6</sub>H<sub>18</sub>N<sub>4</sub> • 2HCl  
**FW:** 219.2  
**Purity:** ≥98%  
**Supplied as:** A 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

Trientine (hydrochloride) is supplied as a solid. A stock solution may be made by dissolving the trientine (hydrochloride) in water. Trientine (hydrochloride) is slightly soluble in water. We do not recommend storing the aqueous solution for more than one day.

### Description

Trientine is a polyamine chelating agent.<sup>1</sup> It reduces the amount of a copper-albumin complex and increases formation of low molecular weight complexes in a dialysis assay in a concentration-dependent manner. Trientine mobilizes copper *ex vivo* in human sera, increasing the supernatant percentage in an analytical ultracentrifugation experiment. *In vivo*, trientine (40 mg/kg, i.v.) mobilizes copper in the kidneys, but not the liver, to increase urinary excretion in rats. Trientine also decreases nickel accumulation in the liver, kidney, and spleen in mice injected with nickel chloride.<sup>2</sup> Formulations containing trientine have been used to reduce copper accumulation in patients with Wilson's disease.

### References

1. Sarkar, B., Sass-Kortsak, A., Clarke, R., *et al.* A comparative study of *in vitro* and *in vivo* interaction of D-penicillamine and triethylenetetramine with copper. *Proc. R. Soc. Med.* **70(Suppl 3)**, 13-18 (1977).
2. Behari, J.R. Use of liposome entrapped triethylenetetramine (TETA) in the treatment of mice loaded with nickel chloride. *Ind. Health* **25(4)**, 221-224 (1987).

#### 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, 12/15/2022

#### 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