

# PRODUCT INFORMATION



## DL-Cystathionine (hydrochloride)

Item No. 17819

**Formal Name:** S-(2-amino-2-carboxyethyl)-homocysteine, dihydrochloride

**Synonym:** DL-Allocystathionine

**MF:** C<sub>7</sub>H<sub>14</sub>N<sub>2</sub>O<sub>4</sub>S • 2HCl

**FW:** 295.2

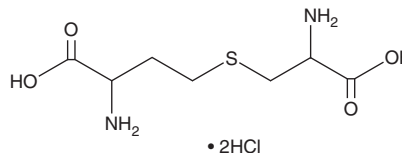
**Purity:** ≥95%

**Supplied as:** A solid

**Storage:** -20°C

**Stability:** ≥4 years

**Special Conditions:** Hygroscopic



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

### Laboratory Procedures

DL-Cystathionine (hydrochloride) is supplied as a solid. Aqueous solutions of DL-cystathionine (hydrochloride) can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of DL-cystathionine (hydrochloride) in PBS (pH 7.2) is approximately 0.25 mg/ml. We do not recommend storing the aqueous solution for more than one day.

### Description

DL-Cystathionine is an intermediate in the synthesis of cysteine that is produced from homocysteine and serine by the action of cystathionine-β-synthase (CBS). CBS can also mediate the condensation of cysteine with homocysteine to produce cystathionine and hydrogen sulfide. An excess of cystathionine in the urine, resulting from a congenital deficiency of cystathionase or an acquired deficiency of vitamin B<sub>6</sub>, is a hallmark of cystathioninuria.

#### 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, 09/06/2023

#### 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