

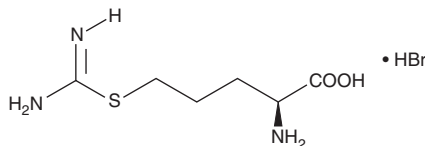
# PRODUCT INFORMATION



## Thioarginine (hydrobromide)

Item No. 80240

**CAS Registry No.:** 2108196-23-0  
**Formal Name:** 5-[(aminoiminomethyl)thio]-L-norvaline, monohydrobromide  
**MF:**  $C_6H_{13}N_3O_2S \cdot HBr$   
**FW:** 272.2  
**Purity:**  $\geq 95\%$   
**Supplied as:** A crystalline solid  
**Storage:**  $-20^\circ C$   
**Stability:**  $\geq 4$  years



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

### Laboratory Procedures

Thioarginine (hydrobromide) is supplied as a crystalline solid. Aqueous solutions of thioarginine (hydrobromide) can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of thioarginine (hydrobromide) in PBS (pH 7.2) is approximately 3 mg/ml. We do not recommend storing the aqueous solution for more than one day.

### Description

Thioarginine is a colorimetric substrate for arginase.<sup>1</sup> Arginase binds to and hydrolyzes thioarginine to release 2-amino-5-mercaptovaleric acid, which reacts with DTNB (Ellman's reagent) to form 2-nitro-5-thiobenzoate and can be quantified by colorimetric detection at 412 nm as a measure of arginase activity.

### Reference

1. Han, S.H. and Viola, R.E. A spectrophotometric assay of arginase. *Anal. Biochem.* **295(1)**, 117-119 (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, 04/08/2024

#### 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