

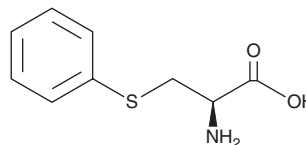
# PRODUCT INFORMATION



## S-Phenylcysteine

Item No. 29734

**CAS Registry No.:** 34317-61-8  
**Formal Name:** S-phenyl-L-cysteine  
**Synonyms:** 3-(phenylthio)-L-Alanine,  
(2R)-2-Amino-3-phenylsulfanylpropanoic acid  
**MF:** C<sub>9</sub>H<sub>11</sub>NO<sub>2</sub>S  
**FW:** 197.3  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 250 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

S-Phenylcysteine is slightly soluble in aqueous solutions. To enhance aqueous solubility, dilute the organic solvent solution into aqueous buffers or isotonic saline. If performing biological experiments, ensure the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. We do not recommend storing the aqueous solution for more than one day.

### Description

S-Phenylcysteine is an adduct that is formed by the binding of benzene oxide to cysteine residues in hemoglobin.<sup>1</sup> It has been found in red blood cells isolated from benzene-exposed rats and mice.

### Reference

1. Bechtold, W.E., Sun, J.S., Birnbaum, L.S., *et al.* S-Phenylcysteine formation in hemoglobin as a biological exposure index to benzene *Arch. Toxicol.* **66**(5), 303-309 (1992).

#### 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, 12/08/2022

#### 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](http://WWW.CAYMANCHEM.COM)