

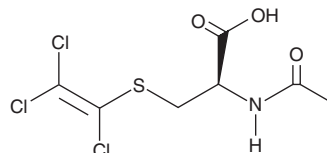
# PRODUCT INFORMATION



## N-acetyl-S-(trichlorovinyl)-L-Cysteine

Item No. 35086

CAS Registry No.: 111348-61-9  
Formal Name: N-acetyl-S-(1,2,2-trichloroethenyl)-L-cysteine  
MF:  $C_7H_8Cl_3NO_3S$   
FW: 292.6  
Purity:  $\geq 90\%$   
Supplied as: A 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

N-acetyl-S-(trichlorovinyl)-L-Cysteine is supplied as a solid. A stock solution may be made by dissolving the N-acetyl-S-(trichlorovinyl)-L-cysteine in the solvent of choice, which should be purged with an inert gas. N-acetyl-S-(trichlorovinyl)-L-Cysteine is soluble in chloroform.

### Description

N-acetyl-S-(trichlorovinyl)-L-Cysteine is a metabolite of the volatile organic compound (VOC) and environmental contaminant tetrachloroethylene (perchloroethylene; PERC).<sup>1,2</sup>

### References

1. Luo, Y.-S., Cichocki, J.A., McDonald, T.J., *et al.* Simultaneous detection of the tetrachloroethylene metabolites S-(1,2,2-trichlorovinyl) glutathione, S-(1,2,2-trichlorovinyl)-L-cysteine, and N-acetyl-S-(1,2,2-trichlorovinyl)-L-cysteine in multiple mouse tissues via ultra-high performance liquid chromatography electrospray ionization tandem mass spectrometry. *J. Toxicol. Environ. Health A* **80**(9), 513-524 (2017).
2. Birner, G., Rutkowska, A., and Dekant, W. N-acetyl-S-(1,2,2-trichlorovinyl)-L-cysteine and 2,2,2-trichloroethanol: Two novel metabolites of tetrachloroethene in humans after occupational exposure. *Drug Metab. Dispos.* **24**(1), 41-48 (1996).

#### 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, 12/09/2022

#### 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