

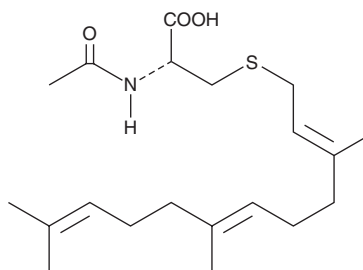
PRODUCT INFORMATION



N-acetyl-S-farnesyl-L-Cysteine

Item No. 63270

CAS Registry No.: 135304-07-3
Formal Name: N-acetyl-S-(3,7,11-trimethyl-2E,6E,10-dodecatrienyl)-L-cysteine
Synonym: AFC
MF: C₂₀H₃₃NO₃S
FW: 367.5
Purity: ≥98%
Supplied as: A solution in methyl acetate
Storage: -20°C
Stability: ≥2 years



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

Laboratory Procedures

N-acetyl-S-farnesyl-L-Cysteine is supplied as a solution in methyl acetate. To change the solvent, simply evaporate the methyl acetate under a gentle stream of nitrogen and immediately add the solvent of choice. Solvents such as ethanol, DMSO, and dimethyl formamide purged with an inert gas can be used. The solubility of N-acetyl-S-farnesyl-L-cysteine in these solvents is approximately 50, 20, and 33 mg/ml, respectively. The solubility of N-acetyl-S-farnesyl-L-cysteine in 0.1 M Na₂CO₃ is approximately 63 mg/ml.

Description

N-acetyl-S-farnesyl-L-Cysteine is a synthetic substrate for the isoprenylated protein methyltransferase (also known as S-adenosylmethionine-dependent methyltransferase).^{1,2} Because it is able to serve as a substrate for the methyltransferase, it effectively functions as an inhibitor of methylation of endogenous isoprenylated proteins.

References

1. Volker, C., Lane, P., Kwee, C., *et al.* A single activity carboxyl methylates both farnesyl and geranylgeranyl cysteine residues. *FEBS Lett.* **295(1-3)**, 189-194 (1991).
2. Pérez-Sala, D., Gilbert, B.A., Tan, E.W., *et al.* Prenylated protein methyltransferases do not distinguish between farnesylated and geranylgeranylated substrates. *Biochem. J.* **284(Pt 3)**, 835-840 (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, 10/31/2023

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