

# N-Acetyl-S-farnesyl-L-cysteine

## Methyltransferase inhibitor

Members of the Ras superfamily of G proteins,  $\gamma$ -subunits of heterotrimeric G proteins and the  $\alpha$ -subunit of cGMP phosphodiesterase from retinal rods are all carboxyl methylated at carboxy-terminal S-farnesylcysteine residues. AFC specifically inhibits the S-farnesylcysteine methyltransferase ( $K_m = 20 \mu\text{M}$ ) in cell free extracts and in whole cells. It inhibits carboxyl methylation of p21ras platelet Rap15 and the  $\gamma$ -subunit of transducin. It blocks receptor-mediated signal transduction in platelets and neutrophils not by inhibiting the methyltransferase, but possibly by inhibiting capacitative  $\text{Ca}^{2+}$  entry.

Citations: 6

[View Online »](#)

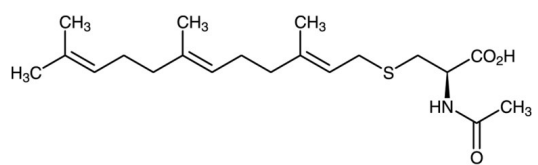
## Ordering Information

[Order Online »](#)

BML-G200-0005	5mg
BML-G200-0025	25mg

## Manuals, SDS & CofA

[View Online »](#)



## Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended.
Handling	Keep under inert gas. Protect from light.
Long Term Storage	-20°C
Shipping	Ambient Temperature

## Regulatory Status

RUO - Research Use Only

## Product Details

Alternative Name	AFC
Appearance	Pale yellow oil.
CAS	135304-07-3
Couple Target	Methyltransferase
Couple Type	Inhibitor
Formula	$C_{20}H_{33}NO_3S$
MW	367.5
Purity	≥98% (HPLC)
Solubility	Soluble in DMSO (25mg/ml), methanol or 100% ethanol (25mg/ml).
Source	Synthetic.

Last modified: May 29, 2024



ENZO LIFE SCIENCES,  
INC.  
Phone: 800.942.0430  
[info-  
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office  
ENZO LIFE SCIENCES  
(ELS) AG  
Phone: +41 61 926 8989  
[info-  
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands  
& Luxembourg  
Phone: +32 3 466 0420  
[info-  
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

France  
Phone: +33 472 440 655  
[info-  
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany  
Phone: +49 7621 5500 526  
[info-  
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland  
Phone (UK customers):  
0845 601 1488  
Phone: +44 1392 825900  
[info-  
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)