

# N-Acetyl-S-geranylgeranyl-L-cysteine

## Blocks methyl esterification of geranylgeranylated proteins

AGGC specifically blocks methyl esterification of geranylgeranylated proteins such as Rab proteins. AGGC blocks receptor-mediated signal transduction in human neutrophils ( $IC_{50}=4\text{ }\mu\text{M}$ ). Induces insulin release from HIT-T15 cells by mimicking the action of Rab3A and inhibits  $\beta 2$ -integrin-induced actin polymerization ( $IC_{50}\sim 45\text{ nM}$ ) in neutrophils.

Citations: 4

[View Online »](#)

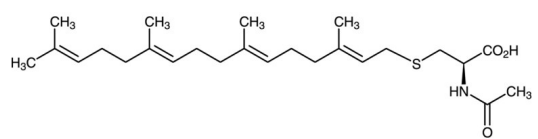
## Ordering Information

[Order Online »](#)

BML-G220-0005	5mg
BML-G220-0025	25mg

## Manuals, SDS & CofA

[View Online »](#)



## Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Store, as supplied, at -20°C for up to 1 year. Store solutions at -20°C for up to 3 months.
Long Term Storage	-20°C
Shipping	Ambient Temperature

## Regulatory Status

RUO - Research Use Only

## Product Details

Appearance	Yellow viscous oil.
CAS	139332-94-8
Couple Target	Methyltransferase
Couple Type	Inhibitor
Formula	$C_{25}H_{41}NO_3S$
MW	435.7
Purity	≥98% (TLC)
Solubility	Soluble in DMSO (>25mg/ml) or 100% ethanol (>25mg/ml).

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-be@enzolifesciences.com](mailto:info-be@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)