

# Farnesylthioacetic acid

## Methyltransferase inhibitor

FTA potently and specifically inhibits methyl esterification of farnesylated proteins such as the  $\gamma$  subunit of transducin (70% at 10 $\mu$ M). It does not act as a substrate for methyl transferase.

Citations: 3

[View Online »](#)

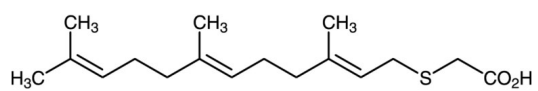
## Ordering Information

[Order Online »](#)

BML-G222-0005	5mg
BML-G222-0025	25mg

## Manuals, SDS & CofA

[View Online »](#)



## Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Solutions are stable for up to 6 months when stored at -20°C.
Handling	Protect from light.
Long Term Storage	-20°C
Shipping	Ambient Temperature

## Regulatory Status

RUO - Research Use Only

## Product Details

Alternative Name	FTA
Appearance	Yellow viscous oil.
CAS	135784-48-4
Couple Target	Methyltransferase
Couple Type	Inhibitor
Formula	$C_{17}H_{28}O_2S$
MW	296.5
Purity	≥98% (TLC)
Solubility	Soluble in DMSO (>25mg/ml) or 100% ethanol (>25mg/ml).
Source	Synthetic.



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)