

# 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 »](#)

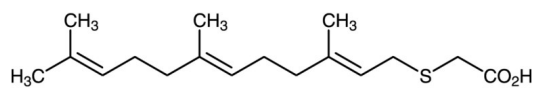
## Ordering Information

[Order Online »](#)

BML-G222-0005	5mg
---------------	-----

## Manuals, SDS & CofA

[View Online »](#)



## Handling & Storage

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

## Regulatory Status

RUO - Research Use Only

## Product Details

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