

S-Farnesyl-L-cysteine methyl ester

Multidrug resistance transporter activator

Stimulates the multidrug resistance transporter ATPase activity (4-5-fold at 10-20 μ M) and competes for drug binding.

Citations: 1

[View Online »](#)

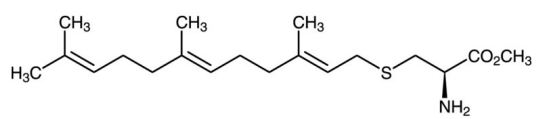
Ordering Information

[Order Online »](#)

BML-G201-0005	5mg
BML-G201-0025	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 under an inert atmosphere in solvent that has been purged with nitrogen or argon at -20°C for up to one month.
Long Term Storage	-20°C
Shipping	Dry Ice

Regulatory Status

RUO - Research Use Only

Product Details

Appearance	Colorless oil.
CAS	125741-64-2
Couple Target	Multidrug resistance transporter
Couple Type	Activator
Formula	$C_{19}H_{33}NO_2S$
MW	339.5
Purity	≥98% (TLC)
Solubility	Soluble in DMSO (25 mg/ml) and ethanol (25 mg/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-
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)