

Chaetomelic acid A

Farnesyltransferase inhibitor

Potent inhibitor of farnesyltransferase in isolated enzyme assays (IC_{50} =55nM) but is inactive in whole cells. Inhibition of farnesyltransferase is selective over geranylgeranyltransferase I and II (IC_{50} =92 μ M and 34 μ M respectively).

Citations: 2

[View Online »](#)

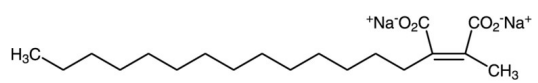
Ordering Information

[Order Online »](#)

BML-G229-0005	5mg
BML-G229-0025	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	White to off-white solid.
CAS	148796-51-4
Couple Target	Farnesyltransferase
Couple Type	Inhibitor
Formula	$C_{19}H_{32}O_4 \cdot 2Na$
MW	370.4
MeltingPoint	>315°C
Purity	≥98% (TLC)
Solubility	Soluble in warm water (2mg/ml, slightly cloudy).
Source	Synthetic.



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

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

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

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

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