

α-

Hydroxyfarnesylphosphonic acid

Farnesyltransferase inhibitor

Potent inhibitor of farnesyltransferase in isolated enzyme assays (IC_{50} =30 nM) and in whole cells. Inhibition of farnesyltransferase is selective over geranylgeranyltransferase I and II (IC_{50} =35.8 μ M and 67 μ M respectively). α -Hydroxyfarnesylphosphonic acid inhibited (1 μ M) Ras processing in Ha-ras-transformed NIH 3T3 fibroblasts.

Citations: 3

[View Online »](#)

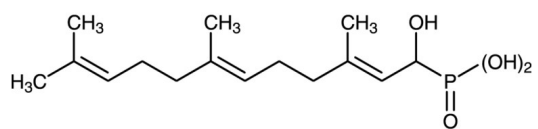
Ordering Information

[Order Online »](#)

BML-G228-0001	1mg
---------------	-----

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Stable for at least 1 year after receipt when stored, as supplied, at -20°C.
Handling	Do not evaporate to dryness. Compound decomposes out of solution.
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Appearance	Liquid.
CAS	140633-12-1
Couple Target	Farnesyltransferase
Couple Type	Inhibitor
Formula	$C_{15}H_{27}O_4P$
Formulation	Solution in ethanol.
MW	302.4
Purity	≥98% (TLC)
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)