

**α-**

# Hydroxyfarnesylphosphonic acid

## Farnesyltransferase inhibitor

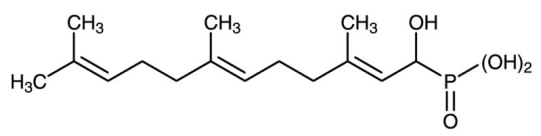
Potent inhibitor of farnesyltransferase in isolated enzyme assays ( $IC_{50}$ =30 nM) and in whole cells. Inhibition of farnesyltransferase is selective over geranylgeranyltransferase I and II ( $IC_{50}$ =35.8  $\mu$ M and 67  $\mu$ M respectively).  $\alpha$ -Hydroxyfarnesylphosphonic acid inhibited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

Citations: 3

[View Online »](#)

Manuals, SDS & CofA

[View Online »](#)



## Handling & Storage

<b>Use/Stability</b>	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.
<b>Handling</b>	Do not evaporate to dryness. Compound decomposes out of solution.
<b>Long Term Storage</b>	-20°C
<b>Shipping</b>	Blue Ice

## Regulatory Status

RUO - Research Use Only

## Product Details

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