

# Piceatannol

Syk inhibitor. SIRT1 activator.

Originally isolated from *Euphorbia lagascae*. Selective protein tyrosine kinase Syk inhibitor ( $IC_{50} \sim 10 \mu M$ ). In RBL-2H3 cells the selective inhibition of Syk by piceatannol results in inhibition of FcεR1-mediated signaling. Activator of human deacetylase SIRT1.

Citations: 16

[View Online »](#)

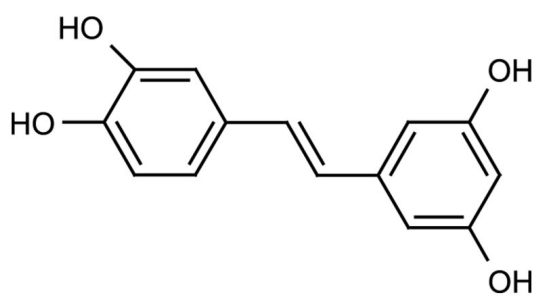
## Ordering Information

[Order Online »](#)

ALX-270-202-M001	1mg
ALX-270-202-M010	10mg

## Manuals, SDS & CofA

[View Online »](#)



## Handling & Storage

<b>Use/Stability</b>	As indicated on product label or CoA when stored as recommended. Unstable in sodium hydroxide.
<b>Handling</b>	Protect from light and oxygen.
<b>Long Term Storage</b>	+4°C
<b>Shipping</b>	Ambient Temperature

## Regulatory Status

RUO - Research Use Only

## Product Details

<b>Alternative Name</b>	3,4,3',5'-Tetrahydroxy-trans-stilbene
<b>Appearance</b>	Off-white to light pink crystalline solid.
<b>CAS</b>	10083-24-6
<b>Couple Target</b>	SIRT, SIRT1, Syk
<b>Couple Type</b>	Activator, Inhibitor
<b>Formula</b>	$C_{14}H_{12}O_4$
<b>MW</b>	244.2
<b>Purity</b>	≥98% (HPLC (UV))
<b>Solubility</b>	Soluble in DMSO (>50mg/ml), methanol (>50mg/ml), dioxan (>50mg/ml), acetonitrile (>50mg/ml); slightly soluble in water (1mg/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-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)