

# Ponasterone A

## Ecdysteroid

A member of the ecdysteroid family. Analog of ecdysone with similar properties to muristerone A. It is a functional, reliable and economical substitute for muristerone A as an inducer for the ecdysone-inducible mammalian expression system. Results show that ponasterone A was able to induce expression of  $\beta$ -galactosidase to levels similar to those obtained with muristerone A induction. Potent compound for ecdysteroid activity; EC<sub>50</sub> = 0.07  $\mu$ M.

Citations: 12

[View Online »](#)

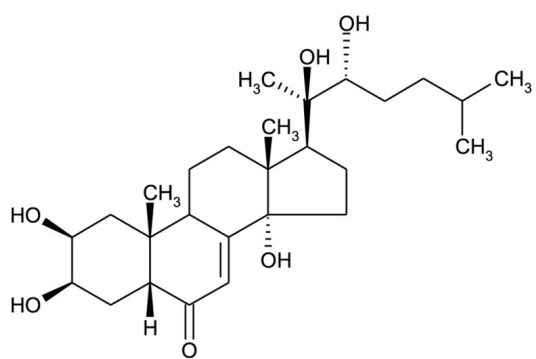
## Ordering Information

[Order Online »](#)

ALX-370-014-M001	1mg
ALX-370-014-M005	5mg

## Manuals, SDS & CofA

[View Online »](#)



## Handling & Storage

<b>Use/Stability</b>	As indicated on product label or CoA when stored as recommended. To prepare a 1mM stock solution (1.1mM) in 100% ethanol: Add 2ml of 100% ethanol to 1mg (2µmol). Gently mix to dissolve the product (5 to 10 min.). Do not heat solution to dissolve the product. Store the 1mM stock solution at +4°C or at -20°C.
<b>Long Term Storage</b>	-20°C
<b>Shipping</b>	Ambient Temperature

## Regulatory Status

RUO - Research Use Only

## Product Details

<b>Alternative Name</b>	25-Deoxy-20-hydroxyecdysone, 25-Deoxyecdysterone, 2β,3β,14α,20R,22R-Pentahydroxy-5β-cholest-7-en-6-one
<b>Appearance</b>	White to off-white powder.
<b>CAS</b>	13408-56-5
<b>Formula</b>	$C_{27}H_{44}O_6$
<b>Identity</b>	Identity determined by NMR.
<b>MI</b>	<b>14:</b> 7588
<b>MW</b>	464.6
<b>Purity</b>	≥95% (HPLC)
<b>Solubility</b>	Soluble in acetic acid, 100% ethanol, methanol or DMSO; insoluble in water.
<b>Source</b>	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)