

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 β -galactosidase to levels similar to those obtained with muristerone A induction. Potent compound for ecdysteroid activity; EC₅₀ = 0.07 μ M.

Citations: 12

[View Online »](#)

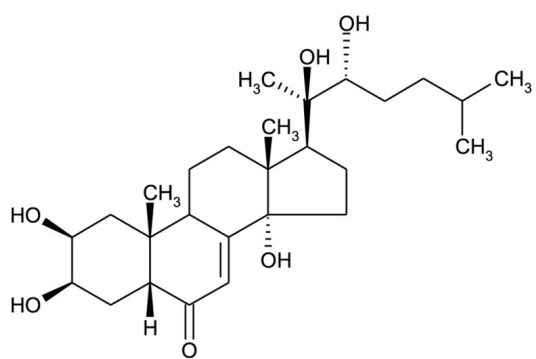
Ordering Information

[Order Online »](#)

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

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	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.
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

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