

Ecdysterone

Natural steroid hormone produced by arthropods; ecdysone agonist.

Ordering Information

[Order Online »](#)

LKT-E0813-M025	25mg
----------------	------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Handling Store under nitrogen.

Long Term Storage -20°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name 20-Hydroxyecdysone, Crustecdysone, Isoinokosterone, Polypodine A

CAS 5289-74-7

Formula $C_{27}H_{44}O_7$

MW 480.64

Purity ≥96%

Solubility Soluble in alcohol.



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com