

# Herboxidiene

## Spliceosome inhibitor

A novel polyketide fermentation product produced by *Streptomyces chromofuscus*, originally discovered by screening for herbicidal activity. Potent and selective inhibitor of spliceosome subunit SF3b. Specifically targets SAP155, a subunit of SF3b responsible for pre-mRNA splicing. Displays anti-angiogenic activity via down-regulation of VEGFR-2 and HIF-1 $\alpha$ . Spliceosome inhibitors have attracted enormous attention due to their potential use in cancer treatment.

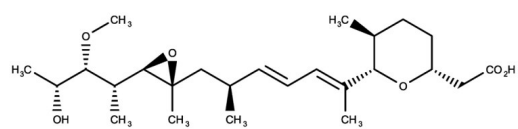
### Ordering Information

[Order Online »](#)

ENZ-CHM487-0005	5mg
-----------------	-----

### Manuals, SDS & CofA

[View Online »](#)



## Handling & Storage

**Use/Stability** As indicated on product label or CoA when stored as recommended.

**Long Term Storage** -20°C

**Shipping** Ambient Temperature

**Regulatory Status** RUO - Research Use Only

## Product Details

**Alternative Name** GEX1A

**Appearance** White solid.

**CAS** 142861-00-5

**Couple Target** Splicing factor 3B

**Couple Type** Inhibitor

**Formula**  $C_{25}H_{42}O_6$

**Identity** Determined by NMR.

**MW** 438.60

**Purity** ≥97% (HPLC)

**Solubility** Soluble in DMSO or ethanol.



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)