# 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**

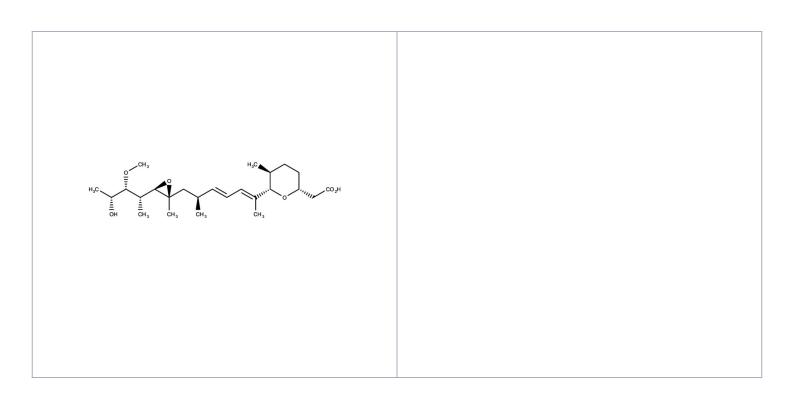
Order Online »

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.

Last modified: February 27, 2025