# C8 ceramine, D-erythro

Ceramide-activated protein kinase activator

 ${
m C_8}$  ceramine is an analog of ceramide in which the amide is substituted with an amine. This substitution renders it inert to ceramidases. It is highly effective at induction of apoptosis in U937 cells. Ceramine induces peak DNA fragmentation in 6 hrs compared to 18 hrs for  ${
m C_8}$  ceramide.

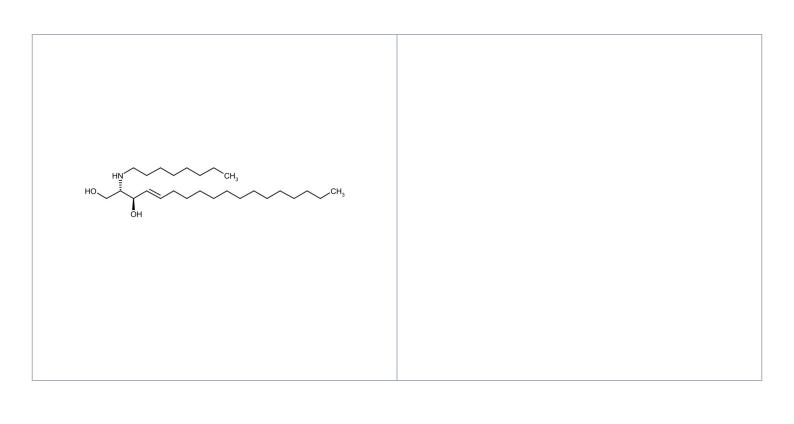
# **Ordering Information**

Order Online »

**BML-SL145-0005** 5mg

Manuals, SDS & CofA

View Online »



## **Handling & Storage**

Use/Stability As indicated on product label or CoA when stored as recommended. Store, as supplied,

at -20°C for up to 1 year. Store solutions at -20°C for up to 3 months.

Long Term Storage -20°C

**Shipping** Ambient Temperature

### Regulatory Status RUO - Research Use Only

### **Product Details**

Appearance White solid.

CAS 170926-06-4

Couple Target Ceramide-activated protein kinase

Couple Type Activator

Formula  $C_{26}H_{53}NO_2$ 

**MW** 411.7

Purity ≥98% (TLC)

**Solubility** Soluble in ethanol (25 mg/ml) or DMSO (25 mg/ml).

Last modified: May 29, 2024

