# SecinH3

Cell-permeable cytohesin family small GEF antagonist.

A Sec7 domain-binding selective antagonist against cytohesin family small GEFs. SecinH3 inhibits large GEFs only at much higher concentrations and is shown to effectively inhibit insulin signaling.

## **Ordering Information**

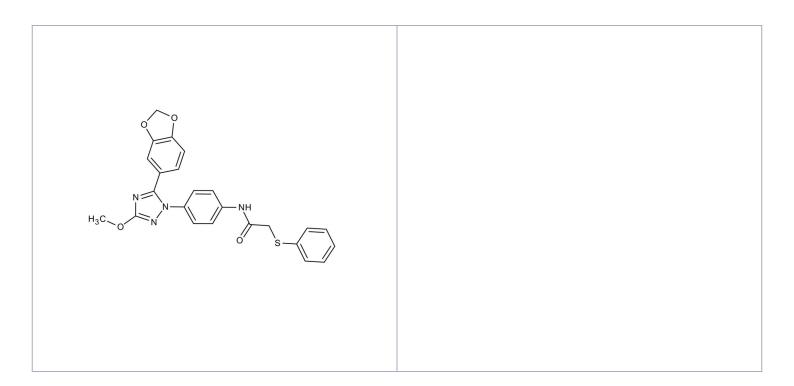
Order Online »

ENZ-CHM114-0005

5mg

Manuals, SDS & CofA

**View Online »** 



### **Handling & Storage**

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

Short Term Storage -20°C

**Long Term Storage** -80°C

**Shipping** Ambient Temperature

#### Regulatory Status RUO - Research Use Only

#### **Product Details**

Alternative Name N-(4-(5-(1,3-Benzodioxol-5-yl)-3-methoxy-1H-1,2,4-triazol-

1-yl)phenyl)-2-(phenylthio)acetamide, Cytohesin inhibitor I

**Appearance** White powder solid.

**CAS** 853625-60-2

Couple Target Cytohesin, Insulin

Couple Type Inhibitor

Formula  $C_{24}H_{20}N_4O_4S$ 

**MW** 460.5

Purity ≥99% (HPLC)

**Soluble** in DMSO or ethanol.

Last modified: May 29, 2024

