

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.

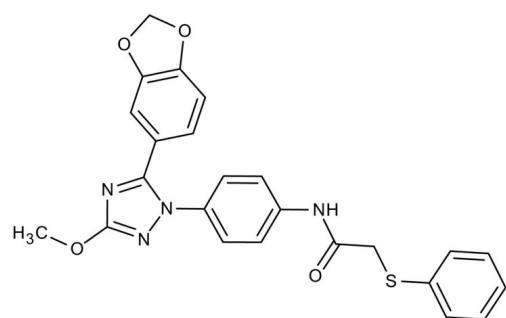
Ordering Information

[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)
Solubility	Soluble in DMSO or ethanol.



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-fr@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