

# SH-5

## Akt inhibitor

Inhibitor of Akt (PKB) activation without affecting activation of the upstream kinase PDK-1, or other kinases downstream of Ras such as MAPK. Induces apoptosis and kills a variety of cancer cell lines containing high levels of active Akt and depending upon Akt for survival.

Citations: 7

[View Online »](#)

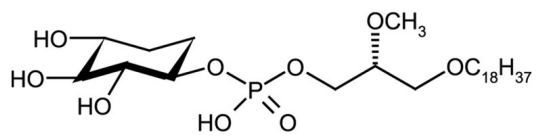
## Ordering Information

[Order Online »](#)

ALX-270-349-M001	1mg
------------------	-----

## Manuals, SDS & CofA

[View Online »](#)



## Handling & Storage

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

**Long Term Storage** +4°C

**Shipping** Ambient Temperature

**Regulatory Status** RUO - Research Use Only

## Product Details

**Alternative Name** D-3-Deoxy-2-O-methyl-myo-inositol 1-[(R)-2-methoxy-3-(octadecyloxy)propyl hydrogen phosphate]

**Appearance** White to off-white solid.

**Couple Target** Akt

**Couple Type** Inhibitor

**Formula**  $C_{29}H_{59}O_{10}P$

**Identity** Determined by 1H-NMR, 31P-NMR, 13C-NMR and MS.

**MW** 598.7

**Purity** ≥98% (TLC)

**Solubility** Soluble in DMSO (10mg/ml) or methanol; slightly soluble in water.

**Last modified: May 29, 2024**



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)