

Akt inhibitor VIII

Potent inhibitor of Akt1, Akt2 and Akt3 activity.

A cell-permeable compound that potently, selectively and reversibly inhibits Akt1/Akt2/Akt3 activity.

Citations: 4

[View Online »](#)

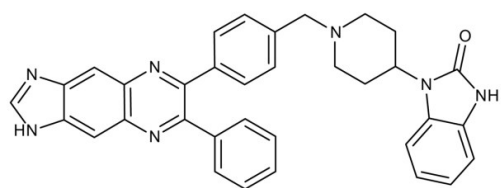
Ordering Information

[Order Online »](#)

ENZ-CHM125-0001	1mg
ENZ-CHM125-0005	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	Akti-1/2, Isozyme-selective Akt inhibitor VIII, 1,3-Dihydro-1-(1-((4-(6-phenyl-1H-imidazo[4,5-g]quinoxalin-7-yl)phenyl)methyl)-4-piperidiny)-2H-benzimidazol-2-one, Imidazoquinoxaline 16h
Appearance	Off-white to yellow powder.
CAS	612847-09-3
Couple Target	Akt
Couple Type	Inhibitor
Formula	$C_{34}H_{29}N_7O$
MW	551.6
Purity	≥98% (HPLC)
Solubility	Soluble in DMSO.

Last modified: May 29, 2024



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