

Compound E

γ -Secretase inhibitor

γ -Secretase is a multimeric and transmembrane aspartyl protease constituted of four subunits: presenilin, nicastrin, Aph-1, and Pen-2. Presenilin is the catalytic subunit of γ -secretase. The carboxyl-terminal fragments (CTFs) of Amyloid precursor protein (APP) and of Notch are two well-known substrates of γ -secretase. Aberrant cleavage of these substrates by γ -secretase is associated with Alzheimer's disease and cancer, respectively. γ -Secretase is, therefore, a promising drug target. Compound E is a cell permeable, potent, selective, and non-competitive inhibitor of γ -secretase. It inhibits the cleavage of both APP and Notch CTFs (IC_{50} =0.3nM for total β -amyloid). At higher concentrations (20-400 μ M), It only weakly affects the activity of presenilin. Compound E was shown to impede ovarian folliculogenesis, promotes neuronal differentiation, and down-regulate thymocyte development.

Citations: 43

[View Online »](#)

Ordering Information

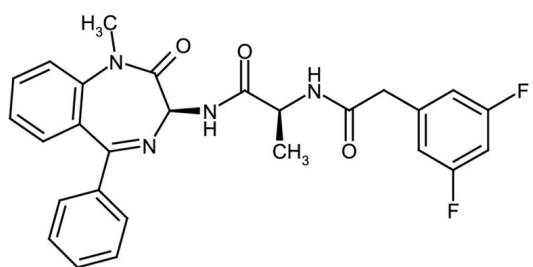
[Order Online »](#)

ALX-270-415-M001	1mg
ALX-270-415-C250	250 μ g

Manuals, SDS & CofA

[View Online »](#)

- Cell-permeable, potent, and non-competitive selective inhibitor of γ -secretase and Notch pathway
- Inducer of neuronal differentiation
- Active in vitro and in vivo



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Stock solutions in DMSO are stable for at least 1 month when stored at -20°C.
Handling	Keep under inert gas.
Short Term Storage	+4°C
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	(2S)-2-[[[(3,5-Difluorophenyl)acetyl]amino]-N-[(3S)-1-methyl-2-oxo-5-phenyl-2,3-dihydro-1H-1,4-benzodiazepin-3-yl]propanamide, γ -Secretase Inhibitor, GSI
Appearance	White to off-white powder.
CAS	209986-17-4
Couple Target	Secretase
Couple Type	Inhibitor
Formula	$C_{27}H_{24}F_2N_4O_3$
Identity	Identity determined by MS and NMR.
MW	490.5
Purity	$\geq 95\%$ (HPLC)
Solubility	Soluble in DMSO (50 mg/ml).



ENZO LIFE SCIENCES,
Last modified: **May 29, 2024**
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-be@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-fr@enzolifesciences.com

France
Phone: +33 472 440 655
info-de@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-uk@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com