

# DAPT

## $\gamma$ -Secretase inhibitor

Cell permeable inhibitor of  $\gamma$ -secretase ( $IC_{50}$ =115nM for total  $\beta$ -amyloid,  $IC_{50}$ =200nM for  $\beta$ -amyloid 1-42). It reduces  $A\beta$  levels *in vivo* in plasma and cerebrospinal fluid in young and aged Tg2576 mice. It blocks the proteolytic processing of neurotrophin receptor like death domain protein (NRADD) Does not inhibit presenilinase. Antagonizes Notch signaling through inhibition of Notch processing. DAPT treatment can influence hematopoietic cell fate decisions and enhances neuronal differentiation in embryonic stem cell-derived embryoid bodies independent of sonic hedgehog (shh) signaling.

Citations: 20

[View Online »](#)

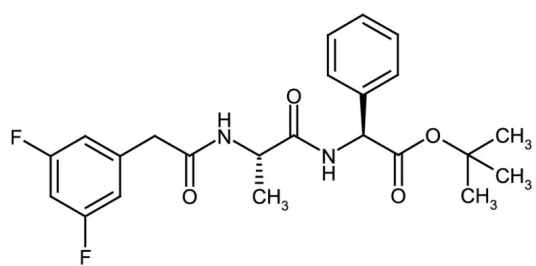
## Ordering Information

[Order Online »](#)

ALX-270-416-M005	5mg
ALX-270-416-M025	25mg

## Manuals, SDS & CofA

[View Online »](#)



## Handling & Storage

<b>Use/Stability</b>	As indicated on product label or CoA when stored as recommended. Stock solutions are stable for up to 3 months when stored at -20°C.
<b>Handling</b>	Protect from light. Packaged under inert gas. After reconstitution, prepare aliquots and store at -20°C.
<b>Long Term Storage</b>	-20°C
<b>Shipping</b>	Ambient Temperature

## Regulatory Status

RUO - Research Use Only

## Product Details

<b>Alternative Name</b>	N-[N-(3,5-Difluorophenacetyl-L-alanyl)]-S-phenylglycine t-butyl ester, $\gamma$ -Secretase inhibitor IX
<b>Appearance</b>	White to off-white solid.
<b>CAS</b>	208255-80-5
<b>Couple Target</b>	Secretase
<b>Couple Type</b>	Inhibitor
<b>Formula</b>	$C_{23}H_{26}F_2N_2O_4$
<b>Identity</b>	Determined by NMR.
<b>MW</b>	432.5
<b>Purity</b>	≥98% (HPLC)
<b>Solubility</b>	Soluble in 100% ethanol, DMSO (20mg/ml) or dichloromethane.



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-  
fr@enzolifesciences.com](mailto:info-fr@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)