

Boc-Gly-Val-Val-CHO

γ -secretase inhibitor

A novel substrate-based (γ -40 site) aldehyde inhibitor of γ -secretase. In whole cells overexpressing amyloid β protein precursor, Boc-GVV-CHO dose-dependently inhibited amyloid β cleavage. γ -secretase is an important therapeutic target for strategies designed to lower amyloid β production.

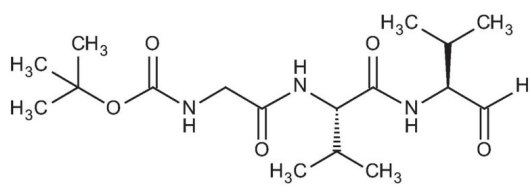
Ordering Information

[Order Online »](#)

BML-PI128-0005	5mg
BML-PI128-0025	25mg

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Store, as supplied, at -20°C for up to 1 year. Store solutions at -20°C for up to 3 months.
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name	Boc-GVV-CHO
Appearance	White to off-white solid.
Formula	$C_{17}H_{31}N_3O_5$
MW	357.5
Purity	≥98% (TLC)
Sequence	Boc-GVV-CHO
Solubility	Soluble in DMSO (>25 mg/ml) or ethanol (10 mg/ml warm).
Source	Synthetic.



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)