Boc-Gly-Val-Val-CHO

y-secretase inhibitor

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

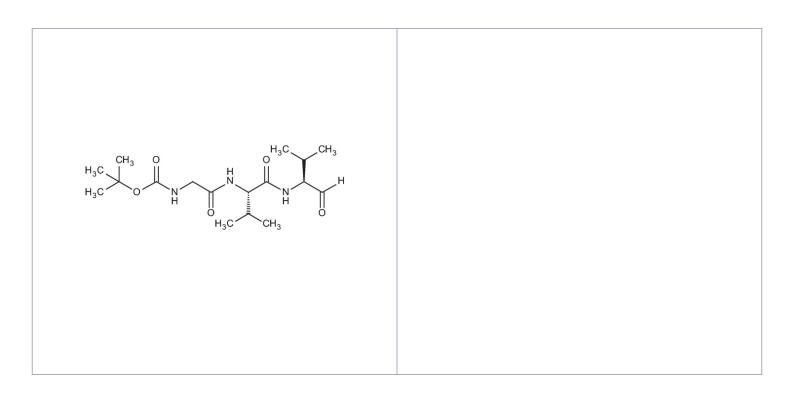
Ordering Information

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.

Last modified: May 29, 2024