

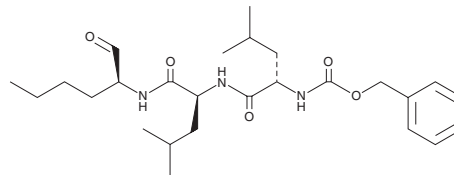
PRODUCT INFORMATION



γ -Secretase Inhibitor I

Item No. 45205

CAS Registry No.: 133407-83-7
Formal Name: N-[(phenylmethoxy)carbonyl]-L-leucyl-N-[(1S)-1-formylpentyl]-L-leucinamide
Synonyms: GSI-I, Z-Leu-Leu-Nle-CHO, Z-LLNle-CHO
MF: C₂₆H₄₁N₃O₅
FW: 475.6
Purity: \geq 95%
Supplied as: A solid
Storage: -20°C
Stability: \geq 4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

γ -Secretase inhibitor I is supplied as a solid. A stock solution may be made by dissolving the γ -Secretase inhibitor I in the solvent of choice, which should be purged with an inert gas. γ -Secretase inhibitor I is sparingly soluble (1-10 mg/ml) in ethanol and soluble (\geq 10 mg/ml) in DMSO.

Description

γ -Secretase inhibitor I is a γ -secretase and proteasome inhibitor.¹ It reduces growth of BT474, T47D, SK-BR-3, and MDA-MB-468 breast cancer cells (ED₅₀s = 2.5, 2.4, 1.6, and 1.4 μ M, respectively) and decreases levels of the γ -secretase cleavage product Notch1 intracellular domain (N1ICD) in the same cells. γ -Secretase inhibitor I reduces proteasome activity in MCF-7 and MDA-MB-231 breast cancer cells when used at concentrations ranging from 1 to 4 μ M.

Reference

1. Han, H., Ma, I., Hendzel, M.J., *et al.* The cytotoxicity of γ -secretase inhibitor I to breast cancer cells is mediated by proteasome inhibition, not by γ -secretase inhibition. *Breast Cancer Res.* **11(4)**, R57 (2005).

WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the [complete](#) Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 04/01/2026

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM