

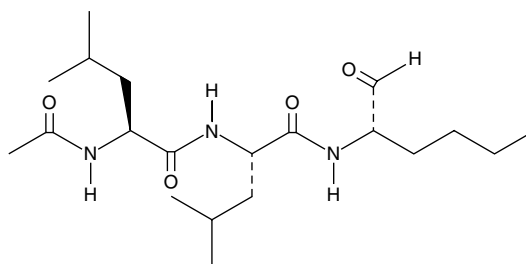
# Product Information



## Calpain Inhibitor I

Item No. 14921

**CAS Registry No.:** 110044-82-1  
**Formal Name:** N-acetyl-L-leucyl-N-[(1S)-1-formylpentyl]-L-leucinamide  
**Synonyms:** Ac-Leu-Leu-Nle-Aldehyde, ALLN, MG 101  
**MF:** C<sub>20</sub>H<sub>37</sub>N<sub>3</sub>O<sub>4</sub>  
**FW:** 383.5  
**Purity:** ≥95%  
**Stability:** ≥2 years at -20°C  
**Supplied as:** A crystalline solid



### Laboratory Procedures

For long term storage, we suggest that calpain inhibitor I be stored as supplied at -20°C. It should be stable for at least two years.

Calpain inhibitor I is supplied as a crystalline solid. A stock solution may be made by dissolving the calpain inhibitor I in the solvent of choice. Calpain inhibitor I is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide, which should be purged with an inert gas. The solubility of calpain inhibitor I in these solvents is approximately 20 mg/ml.

Calpain inhibitor I is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, calpain inhibitor I should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. Calpain inhibitor I has a solubility of approximately 0.5 mg/ml in a 1:1 solution of ethanol:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Calpain inhibitor I is a synthetic tripeptide aldehyde that acts as a potent inhibitor of cysteine proteases including calpain I (K<sub>i</sub> = 190 nM), calpain II (K<sub>i</sub> = 220 nM), cathepsin B (K<sub>i</sub> = 150 nM), and cathepsin L (K<sub>i</sub> = 500 pM).<sup>1</sup> At 10-100 μM calpain inhibitor I dose-dependently prevents the degradation of IκBα and IκBβ by the ubiquitin-proteasome complex, which blocks activation of NFκB and the production of TNF and IL-1β, suggesting a potential therapeutic effect for inflammatory diseases.<sup>2</sup> At 10 μM, calpain inhibitor I can inhibit nitric oxide production by activated macrophages by interfering with transcription of the inducible nitric oxide synthase gene.<sup>3</sup>

### References

1. Sasaki, T., Kishi, M., Saito, M., *et al.* Inhibitory effect of di- and tripeptidyl aldehydes on calpains and cathepsins. *J. Enzyme Inhib.* **3(3)**, 195-201 (1990).
2. Haas, M., Page, S., Page, M., *et al.* Effect of proteasome inhibitors on monocytic IκB-α and -β depletion, NF-κB activation, and cytokine production. *J. Leukoc. Biol.* **63(3)**, 395-404 (1998).
3. Griscavage, J.M., Wilk, S., and Ignarro, L.J. Serine and cysteine proteinase inhibitors prevent nitric oxide production by activated macrophages by interfering with transcription of the inducible NO synthase gene. *Biochem. Biophys. Res. Commun.* **215(2)**, 721-729 (1995).

### Related Products

For a list of related products please visit: [www.caymanchem.com/catalog/14921](http://www.caymanchem.com/catalog/14921)

**WARNING: THIS PRODUCT IS FOR LABORATORY RESEARCH ONLY: NOT FOR ADMINISTRATION TO HUMANS. NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.**

#### SAFETY DATA

This material should be considered hazardous until information to the contrary becomes available. Do not ingest, swallow, or inhale. Do not get in eyes, on skin, or on clothing. Wash thoroughly after handling. This information contains some, but not all, of the information required for the safe and proper use of this material. 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

Cayman Chemical Company makes **no warranty or guarantee** of any kind, whether written or oral, expressed or implied, including without limitation, any warranty of fitness for a particular purpose, suitability and merchantability, which extends beyond the description of the chemicals hereof. Cayman **warrants only** to the original customer that the material will **meet our specifications at the time of delivery**.

Cayman will carry out its delivery obligations with due care and skill. Thus, in no event will Cayman have **any obligation or liability**, whether in tort (including negligence) or in contract, for any direct, indirect, incidental or consequential damages, even if Cayman is informed about their possible existence.

This limitation of liability does not apply in the case of intentional acts or negligence of Cayman, its directors or its employees.

Buyer's **exclusive remedy** and Cayman's sole liability hereunder shall be limited to a **refund** of the purchase price, or at Cayman's option, the **replacement**, at no cost to Buyer, of all material that does not meet our specifications.

Said refund or replacement is conditioned on Buyer giving written notice to Cayman within thirty (30) days after arrival of the material at its destination. Failure of Buyer to give said notice within thirty (30) days shall constitute a waiver by Buyer of all claims hereunder with respect to said material.

For further details, please refer to our **Warranty and Limitation of Remedy located on our website and in our catalog**.

Copyright Cayman Chemical Company, 09/05/2013

### Cayman Chemical

#### Mailing address

1180 E. Ellsworth Road  
Ann Arbor, MI  
48108 USA

#### Phone

(800) 364-9897  
(734) 971-3335

#### Fax

(734) 971-3640

#### E-Mail

[custserv@caymanchem.com](mailto:custserv@caymanchem.com)

#### Web

[www.caymanchem.com](http://www.caymanchem.com)