

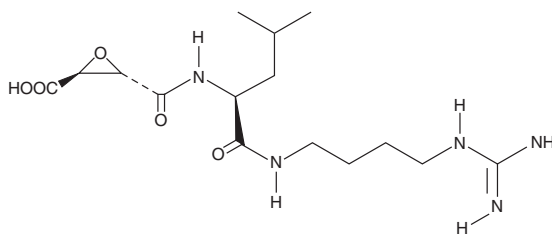
Product Information



E-64

Catalog No. 10007963

CAS Registry No.: 66701-25-5
Formal Name: 3-[[[(1S)-1-[[[4-[(aminoiminomethyl)amino]butyl]amino]carbonyl]-3-methylbutyl]amino]carbonyl]-(2S,3S)-oxiranecarboxylic acid
MF: C₁₅H₂₇N₅O₅
FW: 357.4
Purity: ≥98%
Stability: ≥2 years at -20°C
Supplied as: A crystalline solid



Laboratory Procedures

For long term storage, we suggest that E-64 be stored as supplied at -20°C. It should be stable for at least two years.

E-64 is supplied as a crystalline solid. A stock solution may be made by dissolving the E-64 in an organic solvent purged with an inert gas. E-64 is soluble in organic solvents such as DMSO. The solubility of E-64 in DMSO is approximately 20 mg/ml.

If aqueous stock solutions are required for biological experiments, they can best be prepared by diluting the organic solvent into aqueous buffers or isotonic saline. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. We do not recommend storing the aqueous solution for more than one day.

E-64 is a natural, potent and irreversible inhibitor of cysteine proteases.¹ Its IC₅₀ values for inhibiting cathepsins K, S, and L, *in vitro*, are 1.4, 4.1, and 2.5 nM, respectively.² E-64 also inhibits papain, calpain, and cathepsins B and H, but not serine proteases (*e.g.*, plasmin, trypsin, tissue kallikrein) or aspartic proteases (*e.g.*, pepsin).³ The mode of action involves the formation of a thioether linkage between E-64 and the thiol group of the protease; E-64 does not react with the functional thiol group of non-protease enzymes (*e.g.*, creatine kinase, L-lactate hydrogenase). While E-64 is usually used in biochemical assays, it can also be used in intact cells at higher (μM) concentrations.⁴⁻⁶

References

1. Drahl, C., Cravatt, B.F., and Sorensen, E.J. Protein-reactive natural products. *Natural Products Chemistry* **44**, 5788-5809 (2005).
2. Susa, M., Luong-Nguyen, N.-H., Cappellen, D., *et al.* Human primary osteoclasts: *In vitro* generation and applications as pharmacological and clinical assay. *Journal of Translational Medicine* **2(6)**, (2004).
3. Barrett, A.J., Kembhavi, A.A., Brown, M.A., *et al.* L-trans-Epoxy succinyl-leucylamido(4-guanidino)butane (E-64) and its analogues as inhibitors of cysteine proteinases including cathepsins B, H and L. *Biochem. J.* **201**, 189-198 (1982).
4. Giusti, I., D'Ascenzo, S., Millimaggi, D., *et al.* Cathepsin B mediates the pH-dependent proinvasive activity of tumor-shed microvesicles. *Neoplasia* **10(5)**, 481-488 (2008).
5. Ahlberg, J., Berkenstam, A., Henell, F., *et al.* Degradation of short and long lived proteins in isolated rat liver lysosomes; effects of pH, temperature, and proteolytic inhibitors. *J. Biol. Chem.* **260(9)**, 5847-5854 (1985).
6. Drew, M.E., Banerjee, R., Uffman, E.W., *et al.* Plasmodium food vacuole plasmepsins are activated by falcipains. *J. Biol. Chem.* **283(19)**, 12870-12876 (2008).

Related Products

Lactacystin - Cat. No. 70980 • Clasto-Lactacystin β-lactone - Cat. No. 70988 • Salinosporamide A - Cat. No. 10007311 • Salinosporamide B - Cat. No. 10007563

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

MATERIAL 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 Material 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, 01/04/2010

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

Web

www.caymanchem.com