

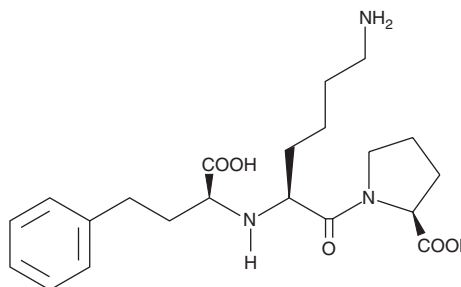
PRODUCT INFORMATION



Lisinopril

Item No. 16833

CAS Registry No.: 76547-98-3
Formal Name: N2-[(1S)-1-carboxy-3-phenylpropyl]-L-lysyl-L-proline
Synonyms: MK-521, MK-522
MF: C₂₁H₃₁N₃O₅
FW: 405.5
Purity: ≥98%
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥4 years



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

Laboratory Procedures

Lisinopril is supplied as a crystalline solid. Lisinopril is sparingly soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. For biological experiments, we suggest that organic solvent-free aqueous solutions of lisinopril be prepared by directly dissolving the lisinopril compound in aqueous buffers. The solubility of lisinopril in PBS (pH 7.2) is approximately 1 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Lisinopril is an angiotensin-converting enzyme (ACE) inhibitor (IC₅₀ = 1.2 nM).¹ It reduces the formation of endothelin-1 (Item No. 24127) and increases nitric oxide (NO) in human vascular endothelial cells when used at a concentration of 0.1 nM.² Lisinopril inhibits the pressor response to angiotensin I in anesthetized rats and dogs (ID₅₀s = 2.3 and 6.5 µg/kg, respectively).¹

References

1. Patchett, A.A., Harris, E., Tristram, E.W., *et al.* A new class of angiotensin-converting enzyme inhibitors. *Nature* **288(5788)**, 280-283 (1980).
2. Desideri, G., Grassi, D., Croce, G., *et al.* Different effects of angiotensin converting enzyme inhibitors on endothelin-1 and nitric oxide balance in human vascular endothelial cells: Evidence of an oxidant-sensitive pathway. *Mediators Inflamm.* 305087 (2008).

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, 12/21/2022

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