

# PRODUCT INFORMATION

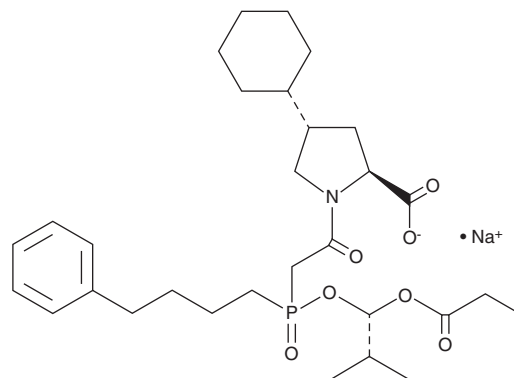


## Fosinopril (sodium salt)

Item No. 21899

**CAS Registry No.:** 88889-14-9  
**Formal Name:** 4S-cyclohexyl-1-[2-[(R)-[(1S)-2-methyl-1-(1-oxopropoxy)propoxy](4-phenylbutyl)phosphinyl]acetyl]-L-proline, monosodium salt

**Synonym:** SQ 28,555  
**MF:** C<sub>30</sub>H<sub>45</sub>NO<sub>7</sub>P • Na  
**FW:** 585.7  
**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

Fosinopril is supplied as a crystalline solid. A stock solution may be made by dissolving the fosinopril (sodium salt) in the solvent of choice. Fosinopril (sodium salt) is soluble in the organic solvent methanol, which should be purged with an inert gas. The solubility of fosinopril (sodium salt) in methanol is approximately 1 mg/ml.

### Description

Fosinopril is a prodrug form of the angiotensin converting enzyme (ACE) inhibitor fosinoprilat.<sup>1</sup> Oral administration of fosinopril inhibits angiotensin I-induced pressor responses in normotensive rats, dogs, and monkeys when administered at doses of 15, 15, and 10 μmol/kg, respectively. Fosinopril (2.5 mg/kg) reduces fractional shortening and decreases left ventricular size in a porcine model of congestive heart failure.<sup>2</sup> Formulations containing fosinopril have been used in the treatment of hypertension and congestive heart failure.

### References

1. DeForrest, J.M., Waldron, T.L., Harvey, C., *et al.* Fosinopril, a phosphinic acid inhibitor of angiotensin I converting enzyme: *In vitro* and preclinical *in vivo* pharmacology. *J. Cardiovasc. Pharmacol.* **14(5)**, 730-736 (1989).
2. McElmurray, J.H., III, Mukherjee, R., New, R.B., *et al.* Angiotensin-converting enzyme and matrix metalloproteinase inhibition with developing heart failure: Comparative effects on left ventricular function and geometry. *J. Pharmacol. Exp. Ther.* **291(2)**, 799-811 (1999).

#### 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, 11/18/2022

#### 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