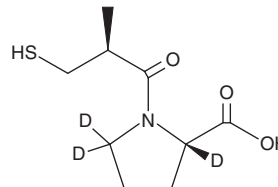


# PRODUCT INFORMATION



## Captopril-d<sub>3</sub> Item No. 25235

**CAS Registry No.:** 1356383-38-4  
**Formal Name:** (S)-1-((S)-3-mercapto-2-methylpropanoyl)pyrrolidine-2-carboxylic-2,5,5-d<sub>3</sub> acid  
**MF:** C<sub>9</sub>H<sub>12</sub>D<sub>3</sub>NO<sub>3</sub>S  
**FW:** 220.3  
**Chemical Purity:** ≥98% (Captopril)  
**Deuterium Incorporation:** ≥99% deuterated forms (d<sub>1</sub>-d<sub>3</sub>); ≤1% d<sub>0</sub>  
**Supplied as:** A 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

Captopril-d<sub>3</sub> is intended for use as an internal standard for the quantification of captopril (Item No. 15313) by GC- or LC-MS. The accuracy of the sample weight in this vial is between 5% over and 2% under the amount shown on the vial. If better precision is required, the deuterated standard should be quantitated against a more precisely weighed unlabeled standard by constructing a standard curve of peak intensity ratios (deuterated versus unlabeled).

Captopril-d<sub>3</sub> is supplied as a solid. A stock solution may be made by dissolving the captopril-d<sub>3</sub> in the solvent of choice, which should be purged with an inert gas. Captopril-d<sub>3</sub> is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of captopril-d<sub>3</sub> in these solvents is approximately 30 mg/ml.

### Description

Captopril is a first generation nonpeptidic ACE inhibitor (IC<sub>50</sub> = 6.3 nM) that was designed based on bradykinin-potentiating peptides isolated from the venom of *B. jararaca*, a pit viper native to Brazil.<sup>1-3</sup> It does not exhibit a domain preference for binding either the C- or N-terminal active sites of the somatic form of ACE.<sup>1</sup> Acute and chronic administration of captopril reduces blood pressure in spontaneously hypertensive Wistar-Kyoto rats and does not induce hypotension in salt-repleted normotensive Wistar-Kyoto rats.<sup>4</sup> Captopril is also a competitive and reversible inhibitor of leukotriene (LTA<sub>4</sub>) hydrolase, which results in the disruption of LTB<sub>4</sub> synthesis with an IC<sub>50</sub> value of 14 μM.<sup>5</sup>

### References

1. Dalkas, G.A., Marchand, D., Galleyrand, J.C., *et al.* Study of a lipophilic captopril analogue binding to angiotensin I converting enzyme. *J. Pept. Sci.* **16(2)**, 91-97 (2010).
2. Redelinghuys, P., Nchinda, A.T., and Sturrock, E.D. Development of domain-selective angiotensin I-converting enzyme inhibitors. *Ann. N.Y. Acad. Sci.* **1056(1)**, 160-175 (2005).
3. Cushman, D.W. and Ondetti, M.A. Design of angiotensin converting enzyme inhibitors. *Nat. Med.* **5(10)**, 1110-1113 (1999).
4. Rubin, B., Antonaccio, M.J., and Horovitz, Z.P. Captopril (SQ 14,225) (D-3-mercapto-2-methylpropanoyl-L-proline): A novel orally active inhibitor of angiotensin-converting enzyme and antihypertensive agent. *Prog. Cardiovasc. Dis.* **21(3)**, 183-194 (1978).
5. Orning, L., Krivi, G., Bild, G., *et al.* Inhibition of leukotriene A<sub>4</sub> hydrolase/aminopeptidase by captopril. *J. Biol. Chem.* **266(25)**, 16507-16511 (1991).

#### 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, 10/31/2022

#### 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