

PRODUCT INFORMATION

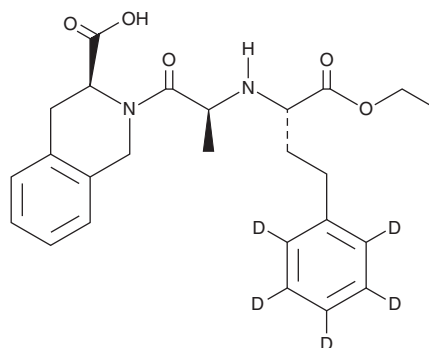


Quinapril-d₅ Item No. 30735

CAS Registry No.: 1279029-79-6
Formal Name: 2-[(2S)-2-[[[(1S)-1-(ethoxycarbonyl)-3-phenyl-d₅-propyl]amino]-1-oxopropyl]-1,2,3,4-tetrahydro-3-isoquinolinecarboxylic acid

Synonym: CI-906-d₅
MF: C₂₅H₂₅D₅N₂O₅
FW: 443.6
Chemical Purity: ≥98% (Quinapril)

Deuterium Incorporation: ≥99% deuterated forms (d₁-d₅); ≤1% d₀
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

Quinapril-d₅ is intended for use as an internal standard for the quantification of quinapril (Item No. 21439) 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).

Quinapril-d₅ is supplied as a solid. A stock solution may be made by dissolving the quinapril-d₅ in the solvent of choice, which should be purged with an inert gas. Quinapril-d₅ is soluble in the organic solvent methanol.

Description

Quinapril is a prodrug form of the angiotensin converting enzyme (ACE) inhibitor quinaprilat.^{1,2} *In vivo*, quinapril (3 mg/kg) reduces mean arterial pressure in renal hypertensive and spontaneously hypertensive rats.² It inhibits angiotensin I-induced pressor responses in normotensive rats and dogs. Quinapril (10-200 mg/kg) prevents left ventricular heart failure in CHF 14.6 cardiomyopathic hamsters. Formulations containing quinapril have been used in the treatment of hypertension, heart failure, and diabetic nephropathy.

References

1. Vago, T., Bevilacqua, M., Conci, F., *et al.* Angiotensin converting enzyme binding sites in human heart and lung: Comparison with rat tissues. *Br. J. Pharmacol.* **107**(3), 821-825 (1992).
2. Kaplan, H.R., Taylor, D.G., and Olson, S.C. Quinapril: Overview of preclinical data. *Clin. Cardiol.* **13**(6 Suppl. 7), VII6-VII12 (1990).

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