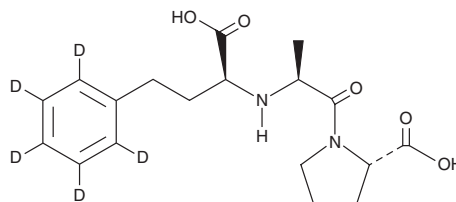


PRODUCT INFORMATION



Enalaprilat-d₅ Item No. 45856

CAS Registry No.: 349554-00-3
Formal Name: N-[(1S)-1-carboxy-3-(phenyl-d₅)propyl]-L-alanyl-L-proline
MF: C₁₈H₁₉D₅N₂O₅
FW: 353.4
Chemical Purity: ≥98% (Enalaprilat)
Deuterium Incorporation: ≥99% deuterated forms (d₁-d₅); ≤1% d₀
Supplied as: A 1 mg/ml solution in methanol
Storage: -20°C
Stability: ≥4 years



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

Laboratory Procedures

Enalaprilat-d₅ is intended for use as an internal standard for the quantification of enalaprilat (Item No. 17603) 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).

Description

Enalaprilat is an angiotensin-converting enzyme (ACE) inhibitor (IC₅₀ = 5.8 nM) and an active metabolite of the prodrug enalapril (Item No. 16041).^{1,2} It reduces the pressor response induced by angiotensin I in hypertensive rats when administered at a dose of 10 mg/kg.²

References

1. Redelinghuys, P., Nchinda, A.T., and Sturrock, E.D. Development of domain-selective angiotensin I-converting enzyme inhibitors. *Ann. N. Y. Acad. Sci.* **1056**, 160-175 (2005).
2. Mencil, J.J., Regan, J.R., Barton, J., et al. Angiotensin converting enzyme inhibitors. 10. Aryl sulfonamide substituted N-[1-carboxy-3-phenylpropyl]-L-alanyl-L-proline derivatives as novel antihypertensives. *J. Med. Chem.* **33**(6), 1606-1615 (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, 06/22/2026

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