

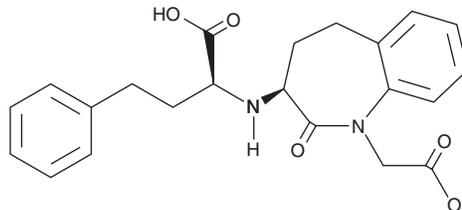
PRODUCT INFORMATION



Benazeprilat

Item No. 21796

CAS Registry No.: 86541-78-8
Formal Name: (3S)-3-[[[(1S)-1-carboxy-3-phenylpropyl]amino]-2,3,4,5-tetrahydro-2-oxo-1H-1-benzazepine-1-acetic acid
Synonym: CGS 14831
MF: C₂₂H₂₄N₂O₅
FW: 396.4
Purity: ≥98%
UV/Vis.: λ_{max}: 239 nm
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

Benazeprilat is supplied as a crystalline solid. A stock solution may be made by dissolving the benazeprilat in the solvent of choice, which should be purged with an inert gas. Benazeprilat is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of benazeprilat in these solvents is approximately 0.5 and 0.16 mg/ml, respectively.

Description

Benazeprilat is an inhibitor of angiotensin-converting enzyme (ACE; IC₅₀ = 0.28 nM in plasma from dog) and the active metabolite of the prodrug benazepril (Item No. 23804).¹ Benazeprilat is produced in the liver following cleavage of the ester group from benazepril. Formulations containing the prodrug benazepril have been used to treat hypertension in human and veterinary medicine.

Reference

1. Toutain, P.L. and Lefèbvre, H.P. Pharmacokinetics and pharmacokinetic/pharmacodynamic relationships for angiotensin-converting enzyme inhibitors. *J. Vet. Pharmacol. Ther.* **27(6)**, 515-525 (2004).

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, 01/03/2023

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