

PRODUCT INFORMATION

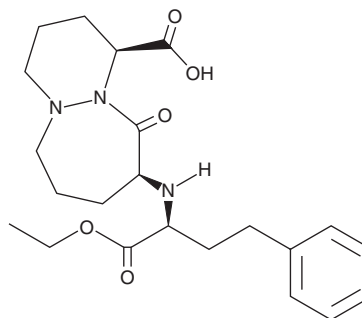


Cilazapril

Item No. 29777

CAS Registry No.: 88768-40-5
Formal Name: (1S,9S)-9-[[[(1S)-1-(ethoxycarbonyl)-3-phenylpropyl]amino]octahydro-10-oxo-6H-pyridazino[1,2-a][1,2]diazepine-1-carboxylic acid

Synonym: Ro 31-2848
MF: C₂₂H₃₁N₃O₅
FW: 417.5
Purity: ≥98%
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

Cilazapril is supplied as a solid. A stock solution may be made by dissolving the cilazapril in the solvent of choice, which should be purged with an inert gas. Cilazapril is slightly soluble in chloroform, methanol, and ethanol.

Description

Cilazapril is a prodrug form of the angiotensin converting enzyme (ACE) inhibitor cilazaprilat.¹ *In vivo*, cilazapril (0.1 mg/kg) inhibits plasma ACE activity and inhibits the angiotensin I-induced pressor response in anesthetized rats and cats. It decreases systolic blood pressure in spontaneously hypertensive rats when administered at a dose of 30 mg/kg. Cilazapril (10 mg/kg, p.o.) decreases blood pressure in volume-depleted renal hypertensive dogs. Formulations containing cilazapril have been used in the treatment of hypertension.

Reference

1. Waterfall, J.F. A review of the preclinical cardiovascular pharmacology of cilazapril, a new angiotensin converting enzyme inhibitor. *Br. J. Clin. Pharmacol.* **27(Suppl 2)**, 139S-150S (1989).

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, 10/28/2022

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