

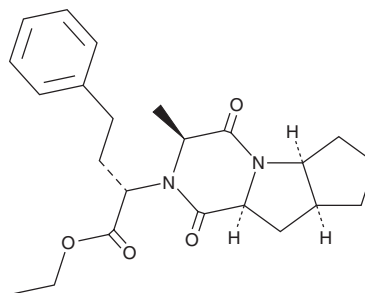
PRODUCT INFORMATION



Ramipril Diketopiperazine

Item No. 34917

CAS Registry No.: 108731-95-9
Formal Name: (αS,3S,5aS,8aS,9aS)-decahydro-3-methyl-1,4-dioxo-α-(2-phenylethyl)-2H-cyclopenta[4,5]pyrrolo[1,2-a]pyrazine-2-acetic acid, ethyl ester
MF: C₂₃H₃₀N₂O₄
FW: 398.5
Purity: ≥95%
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

Ramipril diketopiperazine is supplied as a solid. A stock solution may be made by dissolving the ramipril diketopiperazine in the solvent of choice, which should be purged with an inert gas. Ramipril diketopiperazine is soluble in organic solvents such as ethanol and DMSO.

Description

Ramipril diketopiperazine is a potential impurity found in commercial preparations of ramipril.¹

Reference

1. Orlandini, S., Gotti, R., Giannini, I., *et al.* Development of a capillary electrophoresis method for the assay of ramipril and its impurities: An issue of cis-trans isomerization. *J. Chromatogr. A* **1218(18)**, 2611-2617 (2011).

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, 12/08/2022

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