

PRODUCT INFORMATION

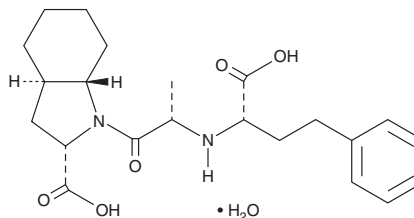


Trandolaprilat (hydrate)

Item No. 29577

CAS Registry No.: 951393-55-8
Formal Name: (2S,3aR,7aS)-1-[(2S)-2-[[[(1S)-1-carboxy-3-phenylpropyl]amino]-1-oxopropyl]octahydro-1H-indole-2-carboxylic acid, monohydrate

Synonym: RU 44403
MF: C₂₂H₃₀N₂O₅ • H₂O
FW: 420.5
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

Trandolaprilat (hydrate) is supplied as a solid. A stock solution may be made by dissolving the trandolaprilat (hydrate) in the solvent of choice, which should be purged with an inert gas. Trandolaprilat (hydrate) is slightly soluble in DMSO and methanol (warmed and sonicated).

Description

Trandolaprilat is an angiotensin-converting enzyme (ACE) inhibitor.¹ It inhibits angiotensin I-induced contraction in isolated rat thoracic aorta (IC₅₀ = 6.9 nM) and potentiates bradykinin-induced contraction in isolated guinea pig ileum with a 50% augmentation concentration (AC₅₀) value of 2.7 nM. Trandolaprilat induces relaxation of norepinephrine- or potassium chloride-precontracted isolated rat thoracic aorta with intact endothelium in a concentration-dependent manner. Trandolaprilat inhibits angiotensin I-induced pressor responses in rats and dogs (ID₅₀s = 9.9 and 7.2 µg/kg, respectively) and potentiates the bradykinin-induced depressor response in rats (ED₅₀ = 4.1 µg/kg).²

References

1. Sekiguchi, N., Ishii, Y., Fujikura, H., *et al.* Pharmacological action of (-)-(2S,3aR,7aS)-1-[(S)-N-[(S)-1-carboxyl-3-phenylpropyl]alanyl]hexahydro-2-indolinecarboxylic acid (trandolaprilat) in isolated smooth muscle preparations. *Gen. Pharmacol.* **24**(3), 585-590 (1993).
2. Brown, N.L., Badel, M.-Y., Benzoni, F., *et al.* Angiotensin-converting enzyme inhibition, anti-hypertensive activity and hemodynamic profile of trandolapril (RU 44570). *Eur. J. Pharmacol.* **148**(1), 79-91 (1988).

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