

PRODUCT INFORMATION



1H-1-ethyl Candesartan Cilexetil

Item No. 17912

CAS Registry No.: 914613-35-7
Formal Name: 2-ethoxy-1-[[[2'-(1-ethyl-1H-tetrazol-5-yl)][1,1'-biphenyl]-4-yl]methyl]-1H-benzimidazole-7-carboxylic acid, 1-[[[(cyclohexyloxy)carbonyl]oxy]ethyl ester

MF: C₃₅H₃₈N₆O₆

FW: 638.7

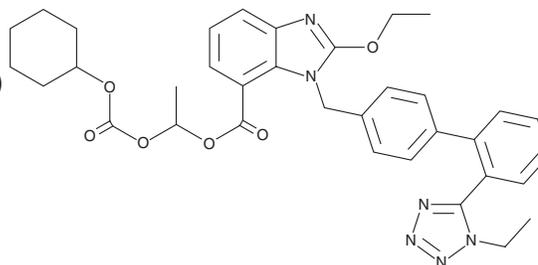
Purity: ≥98%

UV/Vis.: λ_{max}: 254 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

1H-1-ethyl Candesartan cilexetil is supplied as a crystalline solid. A stock solution may be made by dissolving the 1H-1-ethyl candesartan cilexetil in the solvent of choice, which should be purged with an inert gas. 1H-1-ethyl Candesartan cilexetil is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of 1H-1-ethyl candesartan cilexetil in ethanol is approximately 3 mg/ml and approximately 30 mg/ml in DMSO and DMF.

1H-1-ethyl Candesartan cilexetil is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, 1H-1-ethyl candesartan cilexetil should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. 1H-1-ethyl Candesartan cilexetil has a solubility of approximately 0.5 mg/ml in a 1:1 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

1H-1-ethyl Candesartan cilexetil is a potential impurity found in bulk preparations of candesartan cilexetil.^{1,2} It is a degradation product formed by thermal, photolytic, oxidative, or hydrolytic stress during storage.

References

1. Raman, B., Sharma, B.A., Mahale, G., *et al.* Investigation and structural elucidation of a process related impurity in candesartan cilexetil by LC/ESI-ITMS and NMR. *J. Pharm. Biomed. Anal.* **56(2)**, 256-263 (2011).
2. Mehta, S., Shah, R.P., Priyadarshi, R., *et al.* LC and LC-MS/TOF studies on stress degradation behaviour of candesartan cilexetil. *J. Pharm. Biomed. Anal.* **52(3)**, 345-354 (2010).

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, 10/05/2022

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