

PRODUCT INFORMATION

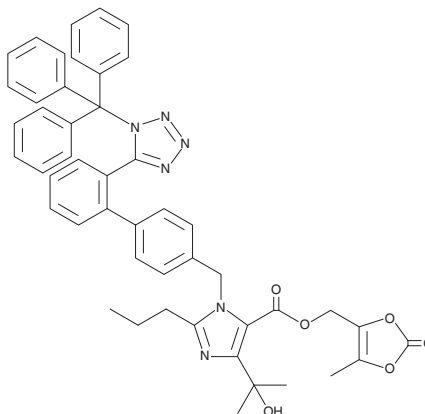


Tritylomesartan Medoxomil

Item No. 34854

CAS Registry No.: 144690-92-6
Formal Name: 4-(1-hydroxy-1-methylethyl)-2-propyl-1-[[2'-[1-(triphenylmethyl)-1H-tetrazol-5-yl][1,1'-biphenyl]-4-yl]methyl]-1H-imidazole-5-carboxylic acid, (5-methyl-2-oxo-1,3-dioxol-4-yl)methyl ester

MF: $C_{48}H_{44}N_6O_6$
FW: 800.9
Purity: $\geq 98\%$
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

Tritylomesartan medoxomil is supplied as a solid. A stock solution may be made by dissolving the tritylomesartan medoxomil in the solvent of choice, which should be purged with an inert gas. Tritylomesartan medoxomil is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of tritylomesartan medoxomil in these solvents is approximately 1 mg/ml.

Description

Tritylomesartan medoxomil is an intermediate in the synthesis of olmesartan medoxomil (Item No. 11614), a prodrug form of the angiotensin II receptor 1 (AT_1) antagonist olmesartan (Item No. 23412).¹

Reference

1. Hanumantha Rao, B., Subramanyeswara Rao, I.V., Ravi Kanth, V., *et al.* A competent and commercially viable process for the synthesis of the anti-hypertensive drug olmesartan medoxomil. *Sci. Pharm.* **83**(3), 465-478 (2015).

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