

PRODUCT INFORMATION



(3R,5R)-Rosuvastatin (calcium salt)

Item No. 35381

CAS Registry No.: 1422515-55-6
Formal Name: (3R,5R,6E)-7-[4-(4-fluorophenyl)-6-(1-methylethyl)-2-[methyl(methylsulfonyl)amino]-5-pyrimidinyl]-3,5-dihydroxy-6-heptenoic acid, hemicalcium salt

MF: C₂₂H₂₈FN₃O₆S • 1/2Ca

FW: 500.6

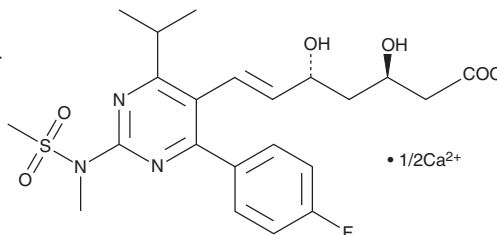
Purity: ≥95%

UV/Vis.: λ_{max}: 242, 310 nm

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

(3R,5R)-Rosuvastatin (calcium salt) is supplied as a solid. A stock solution may be made by dissolving the (3R,5R)-rosuvastatin (calcium salt) in the solvent of choice, which should be purged with an inert gas. (3R,5R)-Rosuvastatin (calcium salt) is soluble in the organic solvent ethanol.

Description

(3R,5R)-Rosuvastatin is a potential impurity found in bulk preparations of the HMG-CoA reductase inhibitor rosuvastatin.¹ It is a degradation product formed by acid or base hydrolysis, thermal, photolytic, hydrolytic, or oxidative stress, or during storage.

Reference

1. Koppala, S., Rayam, P., Reddy, V.R., *et al.* Identification, isolation, characterization and quantification of a new impurity in rosuvastatin calcium tablet dosage form. *Anal. Chem. Ind. J.* **16(10)**, 417-432 (2016).

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, 12/12/2022

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