

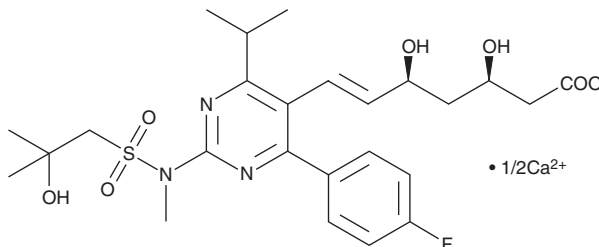
PRODUCT INFORMATION



Rosuvastatin EP Impurity A (calcium salt)

Item No. 35686

CAS Registry No.: 1714147-47-3
Formal Name: 7-[4-(4-fluorophenyl)-2-[[[2-hydroxy-2-methylpropyl)sulfonyl]methylamino]-6-(1-methylethyl)-5-pyrimidinyl]-3,5-dihydroxy-6-heptenoic acid, hemicalcium salt
MF: C₂₅H₃₃FN₃O₇S • 1/2Ca
FW: 558.7
Purity: ≥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

Rosuvastatin EP impurity A (calcium salt) is supplied as a solid. A stock solution may be made by dissolving the rosuvastatin EP impurity A (calcium salt) in the solvent of choice, which should be purged with an inert gas. Rosuvastatin EP impurity A (calcium salt) is soluble in methanol and DMSO.

Description

Rosuvastatin EP impurity A is a potential impurity in commercial preparations of the HMG-CoA reductase inhibitor rosuvastatin.¹

Reference

1. Lee, Y.H., Viji, M., Lee, E., *et al.* Synthesis and characterization of Rosuvastatin calcium impurity A; a HMG-CoA reductase inhibitor. *Tetrahedron Lett.* **58(26)**, 2614-2617 (2017).

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, 01/02/2024

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