

PRODUCT INFORMATION



Atorvastatin (calcium salt)

Item No. 10493

CAS Registry No.: 134523-03-8
Formal Name: 2-(4-fluorophenyl)-βR,δR-dihydroxy-5-(1-methylethyl)-3-phenyl-4-[(phenylamino)carbonyl]-1H-pyrrole-1-heptanoic acid, hemicalcium salt

MF: C₃₃H₃₅FN₂O₅ • 1/2Ca

FW: 578.7

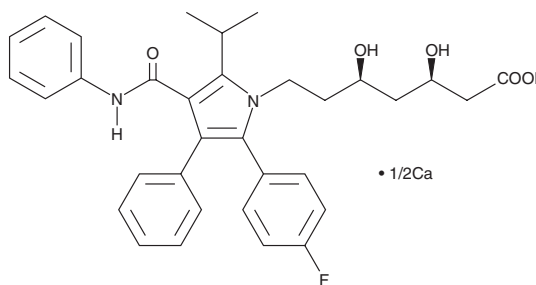
Purity: ≥98%

UV/Vis.: λ_{max}: 247 nm

Supplied as: A crystalline solid

Storage: -20°C

Stability: ≥2 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

Atorvastatin (calcium salt) is supplied as a crystalline solid. A stock solution may be made by dissolving the atorvastatin (calcium salt) in the solvent of choice. Atorvastatin (calcium salt) is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF), which should be purged with an inert gas. The solubility of atorvastatin (calcium salt) in these solvents is approximately 0.5, 15, and 25 mg/ml, respectively.

Atorvastatin (calcium salt) is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, atorvastatin (calcium salt) should first be dissolved in DMF and then diluted with the aqueous buffer of choice. Atorvastatin (calcium salt) has a solubility of approximately 0.1 mg/ml in a 1:9 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Atorvastatin is an inhibitor of HMG-CoA reductase, the rate-limiting enzyme in the mevalonate pathway of cholesterol synthesis, that has IC₅₀ values of 73, 102, and 0.6 nM for HepG2 cells, human fibroblasts, and rat hepatocytes, respectively.¹ Formulations containing atorvastatin have been used in the treatment of hypercholesterolemia and certain dyslipidemias.

Reference

1. Shaw, M.K., Newton, R.S., Sliskovic, D.R., *et al.* Hep-G2 cells and primary rat hepatocytes differ in their response to inhibitors of HMG-CoA reductase. *Biochem. Biophys. Res. Commun.* **170**(2), 726-734 (1990).

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, 05/08/2018

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