

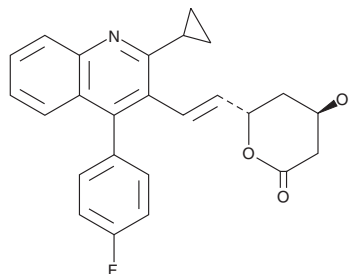
PRODUCT INFORMATION



Pitavastatin lactone

Item No. 21785

CAS Registry No.: 141750-63-2
Formal Name: 6S-[(1E)-2-[2-cyclopropyl-4-(4-fluorophenyl)-3-quinoliny]ethenyl] tetrahydro-4R-hydroxy-2H-pyran-2-one
MF: C₂₅H₂₂FNO₃
FW: 403.5
Purity: ≥98%
Supplied as: A solid
Storage: 4°C
Stability: ≥4 years



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

Laboratory Procedures

Pitavastatin lactone is supplied as a solid. A stock solution may be made by dissolving the pitavastatin lactone in the solvent of choice, which should be purged with an inert gas. Pitavastatin lactone is soluble in the organic solvent chloroform.

Description

Pitavastatin lactone is a major phase 2 metabolite of the HMG-CoA reductase inhibitor pitavastatin (Item No. 15414).^{1,2} Pitavastatin lactone is formed when pitavastatin undergoes glucuronidation by the UDP-glucuronosyltransferase (UGT) isoforms UGT1A1, UGT1A3, or UGT2B7 to form pitavastatin glucuronide, which then undergoes non-enzymatic conversion to pitavastatin lactone. It can be retroconverted to pitavastatin via hydrolysis.

References

1. Fujino, H., Yamada, I., Shimada, S., *et al.* Metabolic fate of pitavastatin, a new inhibitor of HMG-CoA reductase: Human UDP-glucuronosyltransferase enzymes involved in lactonization. *Xenobiotica* **33**(1), 27-41 (2003).
2. Aoki, T., Nishimura, H., Nakagawa, S., *et al.* Pharmacological profile of a novel synthetic inhibitor of 3-hydroxy-3-methylglutaryl-coenzyme A reductase. *Arzneimittelforschung* **47**(8), 904-909 (1997).

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, 12/14/2022

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