

PRODUCT INFORMATION



Rosuvastatin lactone

Item No. 33885

CAS Registry No.: 503610-43-3
Formal Name: N-[4-(4-fluorophenyl)-6-(1-methylethyl)-5-
[(1E)-2-[(2S,4R)-tetrahydro-4-hydroxy-6-
oxo-2H-pyran-2-yl]ethenyl]-2-pyrimidinyl]-
N-methyl-methanesulfonamide

Synonyms: (+)-Rosuvastatin lactone,
Rosuvastatin-5S-lactone, RSTL

MF: C₂₂H₂₆FN₃O₅S

FW: 463.5

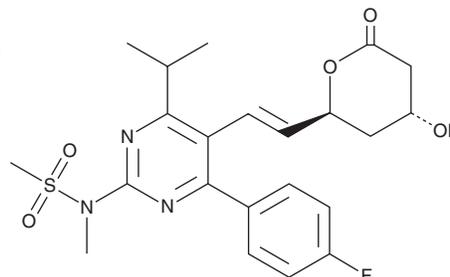
Purity: ≥98%

UV/Vis.: λ_{max}: 245 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

Rosuvastatin lactone is supplied as a solid. A stock solution may be made by dissolving the rosuvastatin lactone in the solvent of choice, which should be purged with an inert gas. Rosuvastatin lactone is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of rosuvastatin lactone in these solvents is approximately 15 and 30 mg/ml, respectively. Rosuvastatin lactone is slightly soluble in ethanol.

Description

Rosuvastatin lactone is the lactone form of the HMG-CoA reductase inhibitor rosuvastatin (Item Nos. 12029 | 18813).¹ It is found *in vivo* following rosuvastatin administration.² Rosuvastatin lactone is an inhibitor of the cytochrome P450 (CYP) isoforms CYP2C9.1 and CYP2C9.3 (IC₅₀s = 8.52 and 4.1 μM, respectively).¹

References

- Shiozawa, A., Yamaori, S., Kamijo, S., *et al.* Effects of acid and lactone forms of statins on S-warfarin 7-hydroxylation catalyzed by human liver microsomes and recombinant CYP2C9 variants (CYP2C9.1 and CYP2C9.3). *Drug Metab. Pharmacokinet.* **36**, 100364 (2021).
- Martin, P.D., Warwick, M.J., Dane, A.L., *et al.* Metabolism, excretion, and pharmacokinetics of rosuvastatin in healthy adult male volunteers. *Clin. Ther.* **25(11)**, 2822-2835 (2003).

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