

PRODUCT INFORMATION



Zaragozic Acid A

Item No. 17452

CAS Registry No.: 142561-96-4

Formal Name: (7S)-11-acetate 5-[(2E,4S,6S)-4,6-dimethyl-2-octenoate]-2,7-anhydro-3,4-di-C-carboxy-8,9,10,12,13-pentadeoxy-10-methylene-12-(phenylmethyl)-L-erythro-L-glycero-D-altro-7-trideculo-7,4-furanosonic acid

Synonyms: L-694,599, Squalostatins S₁

MF: C₃₅H₄₆O₁₄

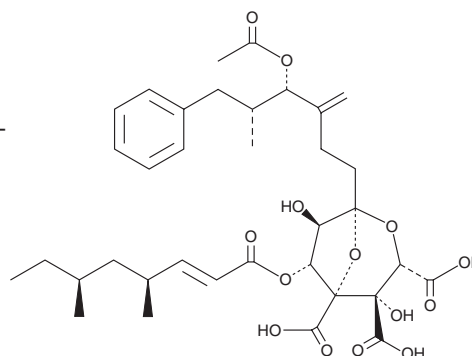
FW: 690.7

Purity: ≥95%

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

Zaragozic acid A is supplied as a solid. A stock solution may be made by dissolving the zaragozic acid A in the solvent of choice, which should be purged with an inert gas. Zaragozic acid A is soluble in organic solvents such as ethanol, methanol, DMSO, and dimethyl formamide.

Description

Squalene synthase catalyzes the first committed step in cholesterol synthesis, mediating the reductive dimerization of farnesyl pyrophosphate (Item No. 63250) to produce squalene.¹ Zaragozic acid A is a fungal metabolite that acts as a reversible competitive inhibitor of squalene synthase ($K_i = 78 \text{ pM}$ *in vitro*).² It dose-dependently reduces cholesterol synthesis in HepG2 cells ($IC_{50} = 6 \text{ }\mu\text{M}$) and inhibits hepatic cholesterol synthesis in mice ($ED_{50} = 0.2 \text{ mg/kg}$).² Zaragozic acid A also inhibits farnesyl transferase (FTase) and geranylgeranyl transferase type I (GGTase I) (IC_{50} s = 216 and 50 nM, respectively).³

References

1. Charlton-Menys, V. and Durrington, P.N. Squalene synthase inhibitors: Clinical pharmacology and cholesterol-lowering potential. *Drugs* **67**(1), 11-16 (2007).
2. Bergstrom, J.D., Kurtz, M.M., Rew, D.J., et al. Zaragozic acids: A family of fungal metabolites that are picomolar competitive inhibitors of squalene synthase. *Proc. Natl. Acad. Sci. USA* **90**(1), 80-84 (1993).
3. Gibbs, J.B., Pompliano, D.L., Mosser, S.D., et al. Selective inhibition of farnesyl-protein transferase blocks ras processing *in vivo*. *J. Biol. Chem.* **268**, 7617-7620 (1993).

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, 11/08/2022

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