

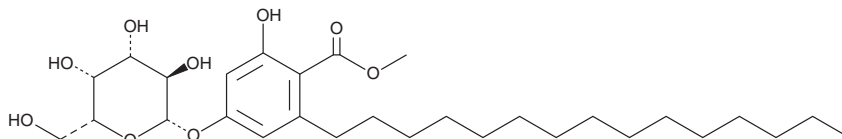
PRODUCT INFORMATION



Aquastatin C

Item No. 44783

CAS Registry No.: 1344044-32-1
Formal Name: 4-(β-D-galactopyranosyloxy)-2-hydroxy-6-pentadecyl-benzoic acid, methyl ester
MF: C₂₉H₄₈O₉
FW: 540.7
Purity: ≥95%
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years
Item Origin: Semi-synthetic



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

Laboratory Procedures

Aquastatin C is supplied as a solid. A stock solution may be made by dissolving the aquastatin C in the solvent of choice, which should be purged with an inert gas. Aquastatin C is soluble in methanol and DMSO.

Description

Aquastatin C is a fungal metabolite that has been found in *Sporothrix*.¹ Unlike aquastatin A (Item No. 33253), aquastatin C does not inhibit bacterial enoyl-ACP reductases, including *S. aureus* enoyl-acyl carrier protein reductase (FabI; IC₅₀ = >100 μM), and is not active against *S. aureus*, methicillin-resistant *S. aureus* (MRSA), or *S. pneumoniae* (MICs = >128 for all).

Reference

1. Kwon, Y.-J., Sohn, M.-J., and Kim, W.-G. Aquastatin C, a new glycoaromatic derivative from *Sporothrix* sp. FN611. *J. Antibiot. (Tokyo)* **64**(2), 213-216 (2011).

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, 11/25/2025

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