

PRODUCT INFORMATION

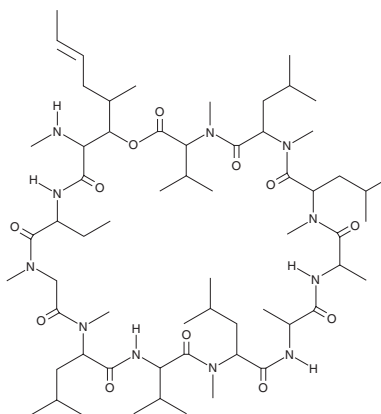


Isocyclosporin A

Item No. 19462

CAS Registry No.: 59865-16-6
Formal Name: (11→1)-lactone-(2S,3R,4R,6E)-3-hydroxy-4-methyl-2-(methylamino)-6-octenoyl-(2S)-2-aminobutanoyl-N-methylglycyl-N-methyl-L-leucyl-L-valyl-N-methyl-L-leucyl-L-alanyl-D-alanyl-N-methyl-L-leucyl-N-methyl-L-leucyl-N-methyl-L-valine

MF: C₆₂H₁₁₁N₁₁O₁₂
FW: 1,202.6
Purity: ≥98%
Supplied as: A powder
Storage: -20°C
Stability: ≥4 years



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

Laboratory Procedures

Isocyclosporin A is supplied as a powder. A stock solution may be made by dissolving the isocyclosporin A in the solvent of choice, which should be purged with an inert gas. Isocyclosporin A is soluble in ethanol, methanol, DMSO, and dimethyl formamide.

Description

Isocyclosporin A is an isomer of cyclosporin A (Item No. 12088) that forms upon acid hydrolysis or upon ionization during mass spectrometry (MS).^{1,2} A method to differentiate the two isomers during MS has been described.¹

References

1. Cirigliano, A. M., and Cabrera, G. M. Differentiation of cyclosporin A from isocyclosporin A by liquid chromatography/electrospray ionization mass spectrometry with post-column addition of divalent metal salt. *Rapid Commun. Mass Spectrom.* **28(5)**, 465-470 (2014).
2. Oliyai, R., and Stella, V. J. Kinetics and mechanism of isomerization of cyclosporin A. *Pharm. Res.* **9(5)**, 617-622 (1992).

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, 09/30/2022

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