

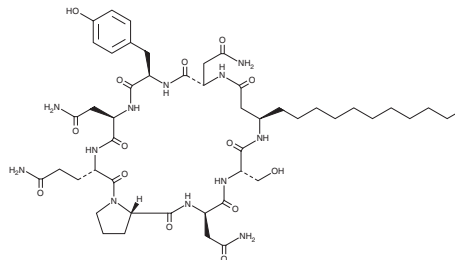
PRODUCT INFORMATION



Iturin A-2

Item No. 38911

CAS Registry No.: 83785-07-3
Formal Name: cyclo[(3R)-3-aminotetradecanoyl-L-asparaginyl-D-tyrosyl-D-asparaginyl-L-glutaminyll-L-prolyl-D-asparaginyl-L-seryl]
Synonym: C14 Iturin
MF: C₄₈H₇₄N₁₂O₁₄
FW: 1,043.2
Purity: ≥90%
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years
Item Origin: Bacterium/*Bacillus* sp.



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

Laboratory Procedures

Iturin A-2 is supplied as a solid. A stock solution may be made by dissolving the Iturin A-2 in the solvent of choice, which should be purged with an inert gas. Iturin A-2 is soluble in methanol, DMSO, and dimethyl formamide.

Description

Iturin A-2 is a cyclic lipopeptide that has been found in *B. subtilis* and has diverse biological activities.¹⁻⁴ It is active against the plant pathogenic fungus *P. expansum* (MIC = 8 µg/disk).¹ Iturin A-2 induces cytotoxicity in MCF-7 and BT474 breast cancer and HeLa cervical cancer cells (IC₅₀s = 66.81, 95.04, and 77.5 µg/ml, respectively).² It reduces southern corn leaf blight lesion area induced by *B. maydis* in greenhouse studies when applied at concentrations of 100 and 300 mg/kg AI and reduces the incidence of southern corn leaf blight in field studies at 300 and 500 mg/kg AI.³ Iturin A-2 (12.5 µg/ml) inhibits cell division, but not nuclear division, in isolated fertilized starfish eggs.⁴

References

1. Bland, J.M. The first synthesis of a member of the iturin family, the antifungal cyclic lipopeptide, iturin-A2. *J. Org. Chem.* **61**(16), (1996).
2. Jiang, J., Zhang, H., Zhang, C., *et al.* Production, purification and characterization of 'iturin A-2' a lipopeptide with antitumor activity from Chinese sauerkraut bacterium *Bacillus velezensis* T701. *Int. J. Pept. Res. Ther.* **27**, 2135-2147 (2007).
3. Ye, Y.F., Li, Q.-Q., Fu, G., *et al.* Identification of antifungal substance (iturin A2) produced by *Bacillus subtilis* B47 and its effect on southern corn leaf blight. *J. Integr. Agric.* **11**(1), 90-99 (2012).
4. Tsuchimori, N., Ikegami, S., Miyashiro, S., *et al.* Isolation of iturin A-2 as an inhibitor of cytokinesis of starfish embryos. *Comp. Biochem. Physiol. Part C* **84**(2), 381-384 (1986).

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, 10/11/2023

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