

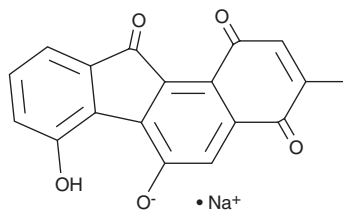
PRODUCT INFORMATION



Fluostatin A (sodium salt)

Item No. 38913

CAS Registry No.: 209673-93-8
Formal Name: 6,7-dihydroxy-3-methyl-1H-benzo[a]fluorene-1,4,11-trione, monosodium salt
MF: C₁₈H₉O₅ • Na
FW: 328.3
Purity: ≥98%
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years
Item Origin: Synthetic



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

Laboratory Procedures

Fluostatin A (sodium salt) is supplied as a solid. A stock solution may be made by dissolving the fluostatin A (sodium salt) in the solvent of choice, which should be purged with an inert gas. Fluostatin A (sodium salt) is soluble in DMSO. Fluostatin A (sodium salt) is also soluble in water. We do not recommend storing the aqueous solution for more than one day.

Description

Fluostatin A is a fluorenone originally isolated from *Streptomyces*.¹ It selectively inhibits dipeptidyl peptidase 3 (DPP-3) over DPP-1, DPP-2, and DPP-4 (IC₅₀s = 0.44, >100, >100, and >100 µg/ml, respectively).

Reference

1. Akiyama, T., Harada, S., Kojima, F., *et al.* Fluostatins A and B, new inhibitors of dipeptidyl peptidase III, produced by *Streptomyces* sp. TA-3391. I. Taxonomy of producing strain, production, isolation, physico-chemical properties and biological properties. *J. Antibiot. (Tokyo)* **51(6)**, 553-559 (1998).

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, 10/10/2023

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