

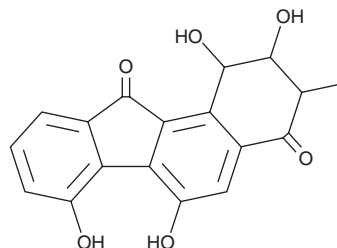
PRODUCT INFORMATION



Fluostatin B

Item No. 37771

CAS Registry No.: 158906-40-2
Formal Name: 2,3-dihydro-1,2,6,7-tetrahydroxy-3-methyl-1H-benzo[a]fluorene-4,11-dione
MF: C₁₈H₁₄O₆
FW: 326.3
Purity: ≥90%
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years
Item Origin: Bacterium/*Streptomyces* sp. TA-3391



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

Laboratory Procedures

Fluostatin B is supplied as a solid. A stock solution may be made by dissolving the fluostatin B in the solvent of choice, which should be purged with an inert gas. Fluostatin B is soluble in DMSO and slightly soluble in acetone and ethyl acetate.

Description

Fluostatin B is a fluorenone originally found in *Streptomyces*.¹ It inhibits dipeptidyl peptidase 3 (DPP-3; IC₅₀ = 24 µg/ml).

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, 02/15/2023

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