

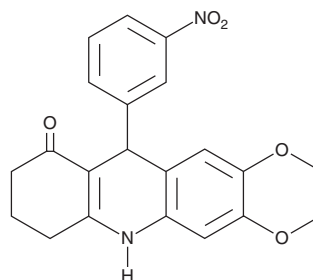
PRODUCT INFORMATION



SSE1806

Item No. 39780

Formal Name: 11-(3-nitrophenyl)-2,3,7,8,9,11-hexahydro-[1,4]dioxino[2,3-b]acridin-10(6H)-one
MF: C₂₁H₁₈N₂O₅
FW: 378.4
Purity: ≥95%
UV/Vis.: λ_{max}: 212 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥2 years



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

Laboratory Procedures

SSE1806 is supplied as a crystalline solid. A stock solution may be made by dissolving the SSE1806 in the solvent of choice, which should be purged with an inert gas. SSE1806 is soluble in DMSO and methanol.

Description

SSE1806 is a microtubule destabilizer.¹ It inhibits tubulin polymerization in a cell-free assay when used at a concentration of 10 μM. SSE1806 inhibits the growth of MCF-7 breast cancer, HCT116 colon cancer, K562 leukemia, and KB-3-1 and paclitaxel-resistant KB-V1 cervical cancer cells (GI₅₀s = 5.02, 2.01, 17.74, 1.29, and 1.74 μM, respectively), as well as induces apoptosis and cell cycle arrest at the G₂/M phase in HCT116 cells.

Reference

1. Firdous, F., Riaz, S., Furqan, M., *et al.* Design, synthesis, and biological evaluation of SSE1806, a microtubule destabilizer that overcomes multidrug resistance. *ACS Med. Chem. Lett.* **14**(10), 1369-1377 (2023).

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, 10/17/2023

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