

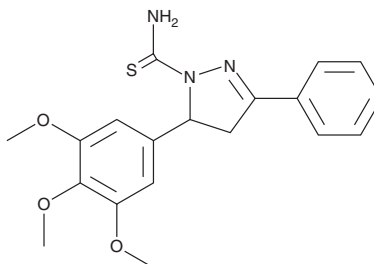
PRODUCT INFORMATION



SSE15206

Item No. 42961

CAS Registry No.: 1370046-40-4
Formal Name: 4,5-dihydro-3-phenyl-5-(3,4,5-trimethoxyphenyl)-1H-pyrazole-1-carbothioamide
MF: C₁₉H₂₁N₃O₃S
FW: 371.5
Purity: ≥98%
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years



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

Laboratory Procedures

SSE15206 is supplied as a solid. A stock solution may be made by dissolving the SSE15206 in the solvent of choice, which should be purged with an inert gas. SSE15206 is sparingly soluble (1-10 mg/ml) in DMSO.

Description

SSE15206 is an inhibitor of tubulin polymerization.¹ It inhibits microtubule polymerization in A549 cells when used at concentrations of 1.25 and 2.5 μ M. SSE15206 (0.5 μ M) induces cell cycle arrest at the G₂/M phase in HCT116 cells. It inhibits the growth of a panel of nine breast, lung, colorectal, cervical, and ovarian cancer cells (GI₅₀s = 0.145-1.269 μ M) and MOLM-13, MV4-11, and K562 leukemia cells (GI₅₀s = 0.264, 0.632, and 0.776 μ M, respectively).

Reference

1. Manzoor, S., Bilal, A., Khan, S., *et al.* Identification and characterization of SSE15206, a microtubule depolymerizing agent that overcomes multidrug resistance. *Sci. Rep.* **8**(1), 3305 (2018).

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, 03/14/2025

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