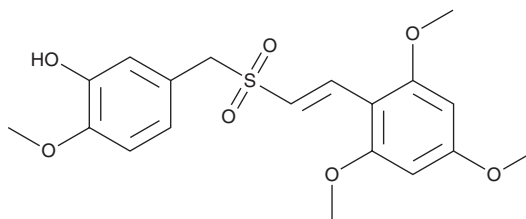


PRODUCT INFORMATION



ON-013100
Item No. 43947

CAS Registry No.: 865783-95-5
Formal Name: 2-methoxy-5-[[[(1E)-2-(2,4,6-trimethoxyphenyl)ethenyl]sulfonyl]methyl]-phenol
MF: C₁₉H₂₂O₇S
FW: 394.4
Purity: ≥95%
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

ON-013100 is supplied as a solid. A stock solution may be made by dissolving the ON-013100 in the solvent of choice, which should be purged with an inert gas. ON-013100 is soluble (≥10 mg/ml) in DMSO.

Description

ON-013100 is a cell cycle inhibitor.¹ It induces cell cycle arrest at the G₂/M phase in Granta 519 mantle cell lymphoma cells when used at a concentration of 1 μM. ON-13100 decreases levels of murine double minute 2 (MDM2), cyclin D1, and cyclin dependent kinase 4 (Cdk4) and increases levels of cyclin B1 in Granta 519 cells. It induces apoptosis in Granta 519 cells when used at a concentration of 0.01 μM. ON-013100 (1 or 10 μM), in combination with doxorubicin (Item No. 15007) or vincristine (Item No. 11764), induces cytotoxicity in Granta 519 and Z-138C mantle cell lymphoma cells.

Reference

1. Park, I.-W., Reddy, M.V.R., Reddy, E.P., *et al.* Evaluation of novel cell cycle inhibitors in mantle cell lymphoma. *Oncogene* 26(38), 5635-5642 (2007).

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, 08/13/2025

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