

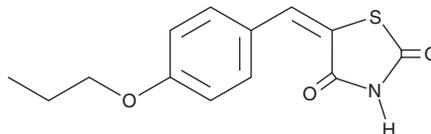
PRODUCT INFORMATION



SMI-16a

Item No. 29456

CAS Registry No.: 587852-28-6
Formal Name: 5-[(4-propoxyphenyl)methylene]-
2,4-thiazolidinedione
MF: C₁₃H₁₃NO₃S
FW: 263.3
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

SMI-16a is supplied as a solid. A stock solution may be made by dissolving the SMI-16a in the solvent of choice, which should be purged with an inert gas. SMI-16a is soluble in DMSO.

Description

SMI-16a is a Pim-1 kinase inhibitor (IC₅₀ = 63 nM).¹ It is selective for Pim-1 over a panel of 60 kinases. SMI-16a (5 μM) inhibits phosphorylation of the Pim-1 target protein Bad in DU145-Pim cells and inhibits the growth of PC3, DU145, LNCaP, K562, and MV4-11 cancer cells. It induces apoptosis and cell cycle arrest at the G₁ phase in DU145 cells.

Reference

1. Beharry, Z., Zemskova, M., Mahajan, S., *et al.* Novel benzylidene-thiazolidine-2,4-diones inhibit Pim protein kinase activity and induce cell cycle arrest in leukemia and prostate cancer cells. *Mol. Cancer Ther.* **8**(6), 1473-1483 (2009).

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, 10/25/2022

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