

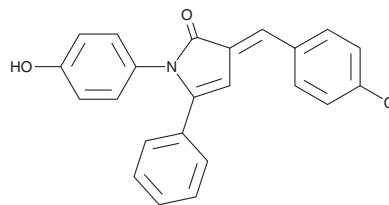
PRODUCT INFORMATION



inS3-54-A18

Item No. 39678

CAS Registry No.: 328998-53-4
Formal Name: 3-[(4-chlorophenyl)methylene]-1,3-dihydro-1-(4-hydroxyphenyl)-5-phenyl-2H-pyrrol-2-one
MF: C₂₃H₁₆ClNO₂
FW: 373.8
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

inS3-54-A18 is supplied as a solid. A stock solution may be made by dissolving the inS3-54-A18 in the solvent of choice, which should be purged with an inert gas. inS3-54-A18 is soluble in acetonitrile, DMSO, and methanol.

Description

inS3-54-A18 is an inhibitor of STAT3.¹ It selectively inhibits the binding of STAT3 to DNA over STAT1 at 25 μM. inS3-54-A18 (20 μM) inhibits STAT3 activity in a reporter assay using MDA-MB-231 breast cancer cells. It induces apoptosis in, and reduces migration and invasion of, A549 non-small cell lung cancer (NSCLC) and MDA-MB-231 cells when used at concentrations of 5 or 10 μM. inS3-54-A18 (200 mg/kg) reduces tumor volume in an A549 mouse xenograft model.

Reference

- Huang, W., Dong, Z., Chen, Y., *et al.* Small-molecule inhibitors targeting the DNA-binding domain of STAT3 suppress tumor growth, metastasis and STAT3 target gene expression *in vivo*. *Oncogene* **35**(6), 783-792 (2016).

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, 11/16/2023

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