

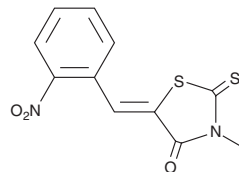
PRODUCT INFORMATION



AKOS B018304

Item No. 35694

CAS Registry No.: 6308-22-1
Formal Name: 5-[(2-nitrophenyl)methylene]-2-thioxo-4-thiazolidinone
Synonym: NSC 43396
MF: C₁₀H₆N₂O₃S₂
FW: 266.3
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

AKOS B018304 is supplied as a solid. A stock solution may be made by dissolving the AKOS B018304 in the solvent of choice, which should be purged with an inert gas. AKOS B018304 is soluble in acetonitrile, methanol, and DMSO.

Description

AKOS B018304 is an arylalkylidene derivative of 1,3-thiazolidin-4-one.¹ It inhibits the binding of ΔFosB to an oligonucleotide containing the activator protein-1 (AP-1) consensus site, Cdk5-AP-1, in a cell-free assay (IC₅₀ = 10.1 μM).² AKOS B018304 (12.5 μM) induces ΔFosB- and ΔFosB/JunD-mediated gene transcription in a reporter assay using Neuro2a cells.

References

1. Jadav, S.S., Sinha, B.N., Hilgenfeld, R., *et al.* Thiazolidone derivatives as inhibitors of chikungunya virus. *Eur. J. Med. Chem.* **89**, 172-178 (2015).
2. Wang, Y., Cesena, T.I., Ohnishi, Y., *et al.* Small molecule screening identifies regulators of the transcription factor ΔFosB. *ACS Chem. Neurosci.* **3**(7), 546-556 (2012).

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, 07/03/2024

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