

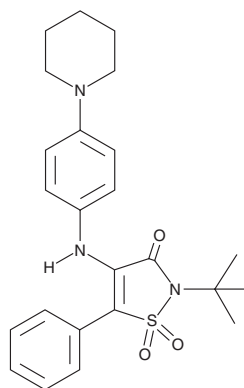
PRODUCT INFORMATION



AZ876

Item No. 34413

CAS Registry No.: 898800-26-5
Formal Name: 2-(1,1-dimethylethyl)-5-phenyl-4-[[4-(1-piperidinyl)phenyl]amino]-3(2H)-isothiazolone, 1,1-dioxide
MF: C₂₄H₂₉N₃O₃S
FW: 439.6
Purity: ≥95%
UV/Vis.: λ_{max}: 254 nm
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

AZ876 is supplied as a solid. A stock solution may be made by dissolving the AZ876 in the solvent of choice, which should be purged with an inert gas. AZ876 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of AZ876 in ethanol is approximately 1 mg/ml and approximately 30 mg/ml in DMSO and DMF.

Description

AZ876 is a liver X receptor (LXR) agonist.^{1,2} It activates LXRA and LXRβ in a transactivation assay in U2OS osteosarcoma cells (EC₅₀s= 6 and 73 nM, respectively).¹ AZ876 reduces phenylephrine-induced increases in cell size and protein synthesis in isolated neonatal rat ventricular myocytes.² It decreases plasma cholesterol and VLDL levels and increases plasma triglyceride and HDL levels, as well as decreases the number and size of aortic root lesions in mice fed a Western diet in an APOE*3-Leiden transgenic mouse model of atherosclerosis when administered at a dose of 20 μmol/kg.¹

References

1. van der Hoorn, J.W.A., Lindén, D., Lindahl, U., *et al.* Low dose of the liver X receptor agonist, AZ876, reduces atherosclerosis in APOE*3Leiden mice without affecting liver or plasma triglyceride levels. *Br. J. Pharmacol.* **162**(7), 1553-1563 (2011).
2. Cannon, M.V., Yu, H., Candido, W.M., *et al.* The liver X receptor agonist AZ876 protects against pathological cardiac hypertrophy and fibrosis without lipogenic side effects. *Eur. J. Heart Fail.* **17**(3), 273-282 (2015).

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, 11/22/2022

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