

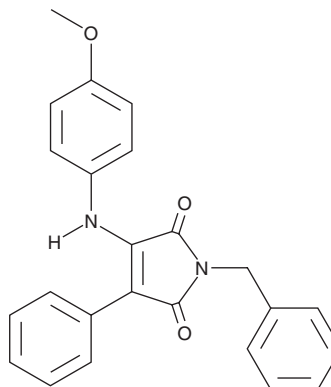
PRODUCT INFORMATION



GSK3987

Item No. 30954

CAS Registry No.: 264206-85-1
Formal Name: 3-[(4-methoxyphenyl)amino]-4-phenyl-1-(phenylmethyl)-1H-pyrrole-2,5-dione
MF: C₂₄H₂₀N₂O₃
FW: 384.4
Purity: ≥98%
UV/Vis.: λ_{max}: 242 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

GSK3987 is supplied as a solid. A stock solution may be made by dissolving the GSK3987 in the solvent of choice, which should be purged with an inert gas. GSK3987 is soluble in organic solvents such as methanol and DMSO. The solubility of GSK3987 in these solvents is approximately 100 and 200 mg/ml, respectively.

Description

GSK3987 is a dual agonist of liver X receptor α (LXRα) and LXRβ.¹ It recruits steroid receptor coactivator 1 (SRC-1) to LXRα and LXRβ in a ligand-sensing assay (LiSA; EC₅₀s = 50 and 40 nM, respectively). GSK3987 is 50-fold selective for LXRα and LXRβ over a panel of nuclear receptors and GSK3β (IC₅₀ = >5,000 nM) but does induce ABCA1 expression in a reporter assay using THP-1 cells (EC₅₀ = 80 nM). It inhibits LPS-induced IL-6 secretion in THP-1 macrophages. GSK3987 increases the expression of the sterol regulatory element binding protein 1c (SREBP-1c) and induces triglyceride accumulation in HepG2 cells in a concentration-dependent manner.

Reference

1. Jaye, M.C., Krawiec, J.A., Campobasso, N., *et al.* Discovery of substituted maleimides as liver X receptor agonists and determination of a ligand-bound crystal structure. *J. Med. Chem.* **48**(17), 5419-5422 (2005).

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, 12/13/2022

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