

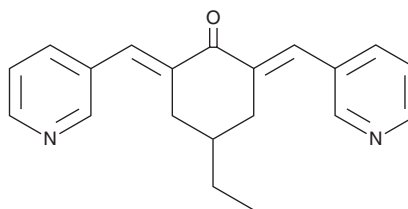
PRODUCT INFORMATION



MCB-613

Item No. 18315

CAS Registry No.: 296792-62-6
Formal Name: 4-ethyl-2,6-bis(3-pyridinylmethylene)-cyclohexanone
MF: C₂₀H₂₀N₂O
FW: 304.4
Purity: ≥95%
UV/Vis.: λ_{max}: 227, 317 nm
Supplied as: A crystalline 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

MCB-613 is supplied as a crystalline solid. A stock solution may be made by dissolving the MCB-613 in the solvent of choice, which should be purged with an inert gas. MCB-613 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of MCB-613 in these solvents is approximately 2 mg/ml.

Description

The p160 steroid receptor coactivator (SRC) family, SRC-1/NCOA1, SRC-2/NCOA2/TIF2/GRIP1, and SRC-3/NCOA3/AIB1/RAC3/ACTR/pCIP, interact with nuclear receptors and other transcription factors to drive target gene expression and integrate growth signaling pathways on which cancer cells rely. MCB-613 is a small molecule stimulator of SRC-1, SRC-2, and SRC-3 that has been used to disrupt cancer cell homeostatic dependence on SRCs.¹ MCB-613 can increase the interaction of SRCs with the coactivators CBP and CARM1 and can induce ER stress coupled to the generation of reactive oxygen species.¹ Overactivation of SRCs by 1-7 μM MCB-613 was shown to selectively and dose dependently induce paraptosis in a variety of cancer cells *in vitro* but not in normal cells.¹ At 20 mg/kg, MCB-613 was also shown to inhibit tumor growth in an MCF-7 breast cancer xenograft model in mice.¹

Reference

1. Wang, L., Yu, Y., Chow, D.-C., *et al.* Characterization of a steroid receptor coactivator small molecule stimulator that overstimulates cancer cells and leads to cell stress and death. *Cancer Cell* **28**, 240-252 (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, 02/15/2024

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