

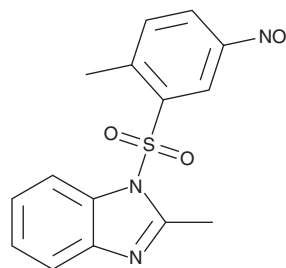
PRODUCT INFORMATION



BI-6015

Item No. 46432

CAS Registry No.: 93987-29-2
Formal Name: 2-methyl-1-[(2-methyl-5-nitrophenyl)sulfonyl]-
1H-benzimidazole
MF: $C_{15}H_{13}N_3O_4S$
FW: 331.3
Purity: $\geq 98\%$
Supplied as: A solid
Storage: $-20^{\circ}C$
Stability: ≥ 4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

BI-6015 is supplied as a solid. A stock solution may be made by dissolving the BI-6015 in the solvent of choice, which should be purged with an inert gas. BI-6015 is soluble (≥ 10 mg/ml) in DMSO.

Description

BI-6015 is a hepatocyte nuclear factor 4 α (HNF4 α) antagonist and a *meta*-substituted positional isomer of CAY10803 (Item No. 12032).¹ It reduces the expression of HNF4 α in T6PNE and MIN6 cells at 2.5 and 5 μ M. It also reduces the expression of a reporter driven by the HNF4 α -responsive ornithine transcarbamylase (OTC) promoter in reporter assays using HepG2 and CV-1 cells. BI-6015 reduces HNF4 α binding to DNA in HepG2 cells at 10 and 20 μ M and induces hepatic steatosis in primary mouse hepatocytes at 5 μ M. It reduces hepatic HNF4 α levels and induces steatosis in the liver, but not the intestine or kidney, in mice. BI-6015 (20 mg/kg per day) also prevents increases in serum aspartate aminotransferase (AST), alanine transaminase (ALT), bilirubin, and bile acid levels in a mouse model of parenteral nutrition-associated cholestasis (PNAC).² It induces cytotoxicity in Hep3B hepatocellular carcinoma (HCC) cells and apoptosis in tumors in an orthotopic Hep3B mouse xenograft model.¹

References

1. Chao, S.H., Fujinaga, K., Marion, J.E., *et al.* Flavopiridol inhibits P-TEFb and blocks HIV-1 replication. *J. Biol. Chem.* **275**(37), 28345-28348 (2000).
2. Ghosh, S., Devereaux, M.W., Orlicky, D.J., *et al.* Pharmacologic inhibition of HNF4 α prevents parenteral nutrition associated cholestasis in mice. *Sci. Rep.* **13**(1), 7752 (2023).

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, 08/18/2026

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