

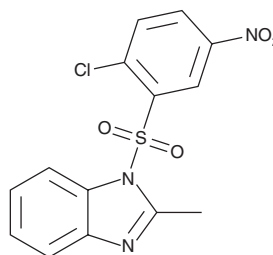
PRODUCT INFORMATION



BIM5078

Item No. 46433

CAS Registry No.: 337506-43-1
Formal Name: 1-[(2-chloro-5-nitrophenyl)sulfonyl]-2-methyl-1H-benzimidazole
MF: $C_{14}H_{10}ClN_3O_4S$
FW: 351.8
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

BIM5078 is supplied as a solid. A stock solution may be made by dissolving the BIM5078 in the solvent of choice, which should be purged with an inert gas. BIM5078 is slightly soluble (0.1-1 mg/ml) in a 1:1 solution of acetonitrile:water.

Description

BIM5078 is a hepatocyte nuclear factor 4 α (HNF4 α) antagonist, a peroxisome proliferator-activated receptor γ (PPAR γ) agonist, and a *meta*-substituted positional isomer of CAY10802 (Item No. 12031).¹ It inhibits the expression of endogenous insulin mRNA in T6PNE pancreatic β -cells (IC_{50} = 930 nM) and induces reporter gene expression in HeLa cells expressing a peroxisome proliferator-response element (PPRE).

Reference

1. Kiselyuk, A., Lee, S.H., Farber-Katz, S., *et al.* HNF4 α antagonists discovered by a high-throughput screen for modulators of the human insulin promoter. *Chem. Biol.* **19**(7), 806-818 (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, 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