

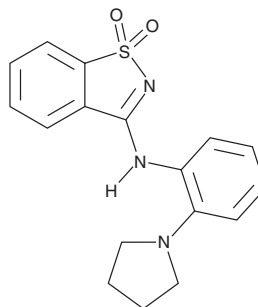
PRODUCT INFORMATION



M1001

Item No. 37393

CAS Registry No.: 874590-32-6
Formal Name: N-[2-(1-pyrrolidinyl)phenyl]-
1,2-benzisothiazol-3-amine,
1,1-dioxide
MF: C₁₇H₁₇N₃O₂S
FW: 327.4
Purity: ≥98%
UV/Vis.: λ_{max}: 210, 255 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

M1001 is supplied as a solid. A stock solution may be made by dissolving the M1001 in the solvent of choice, which should be purged with an inert gas. M1001 is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of M1001 in these solvents is approximately 10 and 15 mg/ml, respectively.

Description

M1001 is a modulator of hypoxia-inducible factor-2α (HIF-2α).¹ It stabilizes the aryl hydrocarbon receptor nuclear translocator (ARNT) and HIF-2α protein-protein interaction and promotes the gene expression of HIF-2α targets VEGF-A, CCND1, SLC2A1, and NDRG1, in 786-O cells when used at a concentration of 10 μM.

Reference

1. Wu, D., Su, X., Lu, J., *et al.* Bidirectional modulation of HIF-2 activity through chemical ligands. *Nat. Chem. Biol.* **15**(4), 367-376 (2019).

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, 02/20/2023

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