

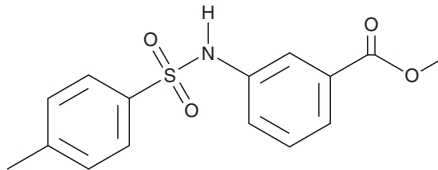
PRODUCT INFORMATION



MSAB

Item No. 36497

CAS Registry No.: 173436-66-3
Formal Name: 3-[[[4-methylphenyl)sulfonyl]amino]-
benzoic acid, methyl ester
MF: C₁₅H₁₅NO₄S
FW: 305.3
Purity: ≥98%
UV/Vis.: λ_{max}: 214 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

MSAB is supplied as a solid. A stock solution may be made by dissolving the MSAB in the solvent of choice, which should be purged with an inert gas. MSAB is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of MSAB in these solvents is approximately 10, 25, and 50 mg/ml, respectively.

Description

MSAB is an inhibitor of Wnt/β-catenin signaling.¹ It binds to β-catenin and promotes its proteasomal degradation in HCT116 cells. MSAB (2-10 μM) selectively inhibits the proliferation of Wnt-dependent cancer cells over non-cancerous cells and Wnt-independent cancer cells. It reduces Wnt3a-induced increases in β-catenin levels in a reporter assay using HEK293T cells. MSAB decreases the expression of the Wnt target genes *AXIN2*, *MYC*, *CCND1*, and *BMP4* in a concentration-dependent manner in HCT116 cells. It reduces tumor growth and increases intratumoral apoptosis in Wnt-dependent HCT116, HT115, and H23 mouse xenograft models when administered at doses of 10 and 20 mg/kg.

Reference

1. Hwang, S.-Y., Deng, X., Byun, S., *et al.* Direct targeting of β-catenin by a small molecule stimulates proteasomal degradation and suppresses oncogenic Wnt/β-catenin signaling. *Cell Rep.* **16(1)**, 28-36 (2016).

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, 05/31/2023

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