

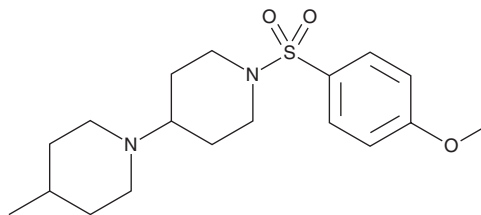
PRODUCT INFORMATION



TASIN-1

Item No. 21155

CAS Registry No.: 792927-06-1
Formal Name: 1'-[(4-methoxyphenyl)sulfonyl]-4-methyl-1,4'-bipiperidine
MF: C₁₈H₂₈N₂O₃S
FW: 352.5
Purity: ≥98%
UV/Vis.: λ_{max}: 241 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

TASIN-1 is supplied as a crystalline solid. A stock solution may be made by dissolving the TASIN-1 in the solvent of choice, which should be purged with an inert gas. TASIN-1 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of TASIN-1 in these solvents is approximately 25, 15, and 30 mg/ml, respectively.

TASIN-1 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, TASIN-1 should first be dissolved in DMF and then diluted with the aqueous buffer of choice. TASIN-1 has a solubility of approximately 0.2 mg/ml in a 1:4 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

TASIN-1 is a small molecule inhibitor that specifically kills cells with adenomatous polyposis coli (APC) truncations (IC₅₀ = 70 nM) but spares cells with wild-type APC.¹ TASIN-1 is effective *in vivo*, inhibiting tumor growth of colorectal cancer (CRC) cells with truncated APC, but not wild-type CRC cells, in a mouse xenograft model.¹ TASIN-1 decreases cholesterol biosynthesis in APC-truncated cells but not in wild-type APC cells.¹

Reference

1. Zhang, L., Theodoropoulos, P.C., Eskiocak, U., *et al.* Selective targeting of mutant adenomatous polyposis coli (APC) in colorectal cancer. *Sci. Transl. Med.* **8(361)**, 361ra140 (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, 12/01/2022

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