

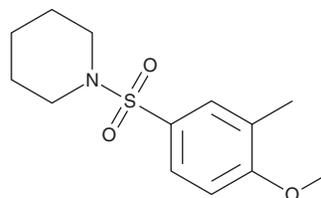
PRODUCT INFORMATION



AA92593

Item No. 35836

CAS Registry No.: 457961-34-1
Formal Name: 1-[(4-methoxy-3-methylphenyl)sulfonyl]-piperidine
MF: C₁₃H₁₉NO₃S
FW: 269.4
Purity: ≥98%
UV/Vis.: λ_{max}: 242 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

AA92593 is supplied as a solid. A stock solution may be made by dissolving the AA92593 in the solvent of choice, which should be purged with an inert gas. AA92593 is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of AA92593 in these solvents is approximately 10 and 16 mg/ml, respectively. AA92593 is also slightly soluble in ethanol.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of AA92593 can be prepared by directly dissolving the solid in aqueous buffers. The solubility of AA92593 in PBS (pH 7.2) is approximately 0.2 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

AA92593 is an opsin 4 (OPN4) receptor antagonist ($K_i = 16$ nM).¹ It inhibits light-induced calcium transients in CHO cells expressing human OPN4 ($IC_{50} = 665$ nM). AA92593 (10 μM) reverses light-induced calcium influx in rat intrinsically photosensitive retinal ganglion cells (ipRGCs). *In vivo*, AA92593 (30 mg/kg) reduces light-induced pupil constriction in mice.

Reference

1. Jones, K.A., Hatori, M., Mure, L.S., *et al.* Small-molecule antagonists of melanopsin-mediated phototransduction. *Nat. Chem. Biol.* **9**(10), 630-635 (2013).

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, 10/06/2022

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