

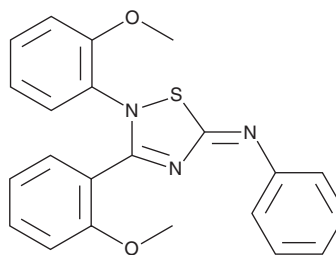
# PRODUCT INFORMATION



**JNJ-10229570**

Item No. 37768

**CAS Registry No.:** 524923-88-4  
**Formal Name:** N-[2,3-bis(2-methoxyphenyl)-1,2,4-thiadiazol-5(2H)-ylidene]-benzenamine  
**MF:** C<sub>22</sub>H<sub>19</sub>N<sub>3</sub>O<sub>2</sub>S  
**FW:** 389.5  
**Purity:** ≥95%  
**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

JNJ-10229570 is supplied as a solid. A stock solution may be made by dissolving the JNJ-10229570 in the solvent of choice, which should be purged with an inert gas. JNJ-10229570 is soluble in the organic solvent DMSO.

## Description

JNJ-10229570 is an antagonist of melanocortin receptor 1 (MC1R), MC4R, and MC5R (IC<sub>50</sub>s = 270, 240, and 200 nM, respectively, in radioligand binding assays).<sup>1</sup> It prevents bovine pituitary extract-induced increases in lipid droplets in primary human sebocytes when used at a concentration of 0.05 μM. JNJ-10229570 (1 μM) inhibits cholera toxin-induced increases in squalene and cholesterol esters, but not triglycerides, in primary human sebocytes. Topical application of JNJ-10229570 (0.05%) inhibits sebum secretion from, and sebaceous gland differentiation in, transplanted human skin in a mouse xenograft model.

## Reference

1. Eisinger, M., Li, W.-H., Anthonavage, M., *et al.* A melanocortin receptor 1 and 5 antagonist inhibits sebaceous gland differentiation and the production of sebum-specific lipids. *J. Dermatol. Sci.* **63**(1), 23-32 (2011).

### 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, 04/27/2023

## 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