

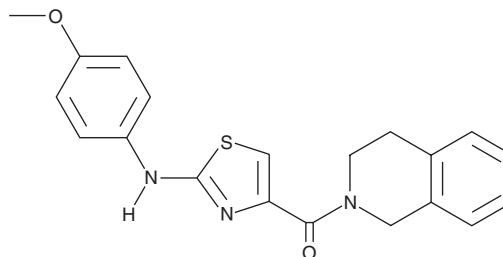
# PRODUCT INFORMATION



## GSK 417651A

Item No. 15268

**CAS Registry No.:** 736945-96-3  
**Formal Name:** (3,4-dihydro-2(1H)-isoquinolinyloxy)  
[2-[(4-methoxyphenyl)amino]-4-thiazolyl]-methanone  
**MF:** C<sub>20</sub>H<sub>19</sub>N<sub>3</sub>O<sub>2</sub>S  
**FW:** 365.5  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 271 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

GSK 417651A is supplied as a solid. A stock solution may be made by dissolving the GSK 417651A in the solvent of choice, which should be purged with an inert gas. GSK 417651A is soluble in DMSO.

### Description

GSK 417651A is an inhibitor of transient receptor potential canonical 3 (TRPC3) and TRPC6 (IC<sub>50</sub> = ~0.4 μM for both).<sup>1</sup> It inhibits migration of isolated human fibroblastic valve interstitial cells into the wounded area in a scratch test assay when used at a concentration of 2 μM.<sup>2</sup>

### References

1. Xu, X., Lozinskaya, I., Costell, M., *et al.* Characterization of small molecule TRPC3 and TRPC6 agonist and antagonists. *Biophys. J.* **104**(2), 454A (2013).
2. Al-Shammari, H., Latif, N., Sarathchandra, P., *et al.* Expression and function of mechanosensitive ion channels in human valve interstitial cells. *PLoS One* **15**(10), e0240532 (2020).

#### 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, 11/28/2022

#### 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