

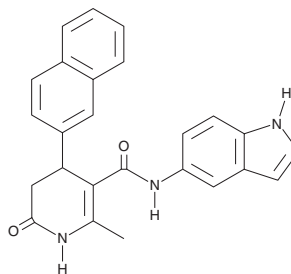
# PRODUCT INFORMATION



**GSK270822A**

Item No. 43717

**CAS Registry No.:** 864082-23-5  
**Formal Name:** 1,4,5,6-tetrahydro-N-1H-indazol-5-yl-2-methyl-4-(2-naphthalenyl)-6-oxo-3-pyridinecarboxamide  
**MF:** C<sub>24</sub>H<sub>20</sub>N<sub>4</sub>O<sub>2</sub>  
**FW:** 396.5  
**Purity:** ≥98%  
**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

GSK270822A is supplied as a solid. A stock solution may be made by dissolving the GSK270822A in the solvent of choice, which should be purged with an inert gas. GSK270822A is soluble in the organic solvent DMSO at a concentration of approximately 10 mg/ml.

## Description

GSK270822A is an inhibitor of G protein-coupled receptor kinase 2 (GRK2; IC<sub>50</sub> = 12.6 μM).<sup>1</sup> It is selective for GRK2 over GRK1, GRK5, and PKA (IC<sub>50</sub>s = 794, 63.1, and 50.1 μM, respectively).

## Reference

1. Homan, K.T., Larimore, K.M., Elkins, J.M., *et al.* Identification and structure-function analysis of subfamily selective G protein-coupled receptor kinase inhibitors. *ACS Chem. Biol.* **10**(1), 310-319 (2015).

### 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, 09/30/2025

## 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](http://WWW.CAYMANCHEM.COM)