

# PRODUCT INFORMATION

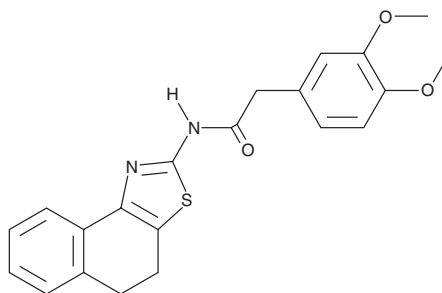


**ZINC00881524**

Item No. 30788

**CAS Registry No.:** 557782-81-7  
**Formal Name:** N-(4,5-dihydronaphtho[1,2-d]thiazol-2-yl)-3,4-dimethoxybenzeneacetamide

**MF:** C<sub>21</sub>H<sub>20</sub>N<sub>2</sub>O<sub>3</sub>S  
**FW:** 380.5  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 232, 282 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

ZINC00881524 is supplied as a crystalline solid. A stock solution may be made by dissolving the ZINC00881524 in the solvent of choice, which should be purged with an inert gas. ZINC00881524 is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of ZINC00881524 in these solvents is approximately 30 mg/ml.

ZINC00881524 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, ZINC00881524 should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. ZINC00881524 has a solubility of approximately 0.33 mg/ml in a 1:2 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

## Description

ZINC00881524 is a Rho-associated kinase (ROCK) inhibitor.<sup>1</sup> It decreases levels of ROCK1 in T47D and CAMA-1 breast cancer cells. ZINC00881524 decreases proliferation of T47D and CAMA-1 cells when used in combination with CNN1 knockdown.

## Reference

1. Wang, Z., Li, T.-E., Chen, M., *et al.* miR-106b-5p contributes to the lung metastasis of breast cancer via targeting CNN1 and regulating Rho/ROCK1 pathway. *Aging (Albany NY)* **12(2)**, 1867-1887 (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, 12/09/2022

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