

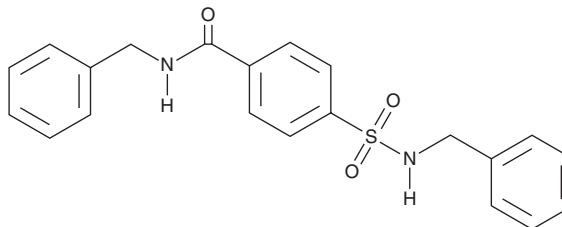
# PRODUCT INFORMATION



**TH263**

Item No. 37295

**CAS Registry No.:** 313520-94-4  
**Formal Name:** N-(phenylmethyl)-4-[[[(phenylmethyl)amino]sulfonyl]-benzamide  
**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>: 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

TH263 is supplied as a solid. A stock solution may be made by dissolving the TH263 in the solvent of choice, which should be purged with an inert gas. TH263 is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of TH263 in these solvents is approximately 10 mg/ml.

## Description

TH263 is an inactive analog of the inhibitor of LIM kinase 1 (LIMK1) and LIMK2 TH257 (Item No. 34236) that has been used as a negative control for the kinase activity of LIMK1 and LIMK2.<sup>1</sup>

## Reference

1. Hanke, T., Mathea, S., Woortman, J., *et al.* Development and characterization of type I, type II, and type III LIM-kinase chemical probes. *J. Med. Chem.* **65(19)**, 13264-13287 (2022).

### 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, 10/02/2023

## 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