

PRODUCT INFORMATION



TH257

Item No. 34236

CAS Registry No.: 2244678-29-1

Formal Name: N-butyl-4-[(phenylamino)sulfonyl]-
N-(phenylmethyl)-benzamide

MF: C₂₄H₂₆N₂O₃S

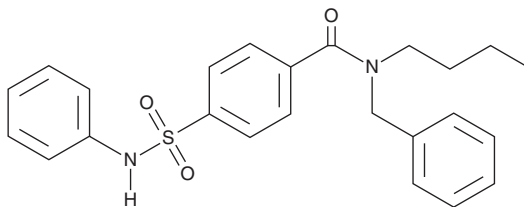
FW: 422.5

Purity: ≥98%

Supplied as: A solution in ethanol

Storage: -20°C

Stability: ≥2 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

TH257 is supplied as a solution in ethanol. To change the solvent, simply evaporate the ethanol under a gentle stream of nitrogen and immediately add the solvent of choice. TH257 is soluble in DMSO.

Description

TH257 is an inhibitor of LIM kinase 1 (LIMK1) and LIMK2 (IC₅₀s = 84 and 39 nM, respectively).¹

Reference

1. Manetti, F. Recent advances in the rational design and development of LIM kinase inhibitors are not enough to enter clinical trials. *Eur. J. Med. Chem.* **155**, 445-458 (2018).

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, 11/18/2022

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