

PRODUCT INFORMATION



BC-1382

Item No. 43143

CAS Registry No.: 1013753-99-5

Formal Name: N-[(1S)-2-[[[(2-methoxyphenyl)methyl]amino]-1-methyl-2-oxoethyl]-1-(phenylsulfonyl)-4-piperidinecarboxamide

MF: C₂₃H₂₉N₃O₅S

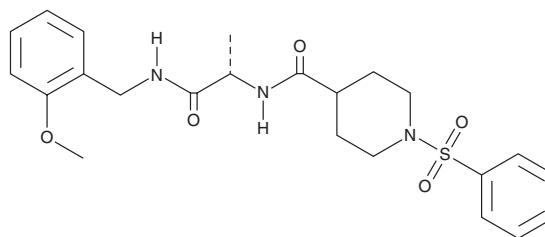
FW: 459.6

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

BC-1382 is supplied as a solid. A stock solution may be made by dissolving the BC-1382 in the solvent of choice, which should be purged with an inert gas. BC-1382 is sparingly soluble (1-10 mg/ml) in ethanol and DMSO.

Description

BC-1382 is an inhibitor of the protein-protein interaction between E3 ubiquitin-protein ligase HECTD2 and E3 SUMO-protein ligase PIAS1.¹ It inhibits the HECTD2-PIAS1 protein-protein interaction (IC₅₀ = ~5 nM in a pull-down assay) and increases the stability and levels of PIAS1. BC-1382 (10 mg/kg) decreases Tnf and Il-6 levels in bronchoalveolar lavage fluid (BALF) in mouse models of *P. aeruginosa*- or LPS-induced pneumonia.

Reference

1. Coon, T.A., McKelvey, A.C., Lear, T., *et al.* The proinflammatory role of HECTD2 in innate immunity and experimental lung injury. *Sci. Transl. Med.* **7**(295), 295ra109 (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, 04/09/2025

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