

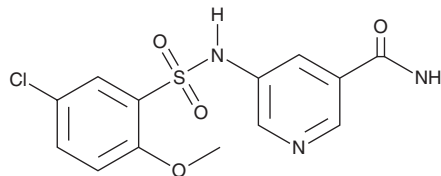
PRODUCT INFORMATION



SBI-425

Item No. 34626

CAS Registry No.: 1451272-71-1
Formal Name: 5-[[[(5-chloro-2-methoxyphenyl)sulfonyl]amino]-3-pyridinecarboxamide
MF: C₁₃H₁₂ClN₃O₄S
FW: 341.8
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

SBI-425 is supplied as a solid. A stock solution may be made by dissolving the SBI-425 in the solvent of choice, which should be purged with an inert gas. SBI-425 is soluble in DMSO.

Description

SBI-425 is an inhibitor of tissue-nonspecific alkaline phosphatase (TNAP; IC₅₀ = 0.016 μM).¹ It is selective for TNAP over a panel of 35 targets at 10 μM. SB-425 (10 and 30 mg/kg) increases survival and nearly completely inhibits the development of medial arterial calcification in a mouse model of chronic kidney disease-mineral bone disorder (CKD-MBD) induced by a high-adenine high-phosphate diet.²

References

1. Pinkerton, A.B., Sergienko, E., Bravo, Y., *et al.* Discovery of 5-((5-chloro-2-methoxyphenyl)sulfonamido) nicotinamide (SBI-425), a potent and orally bioavailable tissue-nonspecific alkaline phosphatase (TNAP) inhibitor. *Bioorg. Med. Chem. Lett.* **28(1)**, 31-34 (2018).
2. Tani, T., Fujiwara, M., Orimo, H., *et al.* Inhibition of tissue-nonspecific alkaline phosphatase protects against medial arterial calcification and improves survival probability in the CKD-MBD mouse model. *J. Pathol.* **250(1)**, 30-41 (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, 11/22/2022

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