PRODUCT INFORMATION



4SC-202

Item No. 26175

CAS Registry No.: Formal Name:	1186222-89-8 (2E)-N-(2-aminophenyl)-3-[1-[[4- (1-methyl-1H-pyrazol-4-yl)phenyl] sulfonyl]-1H-pyrrol-3-yl]-2-propenamide,	N I I	H —
MF: FW: Purity:	$C_{23}H_{21}N_5O_3S \bullet C_7H_8O_3S$ 619.7 ≥98%		
Supplied as: Storage: Stability:	A solid -20°C ≥4 years	• CH ₃ C _e H ₄ SO ₃ H	

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

Laboratory Procedures

4SC-202 is supplied as a solid. A stock solution may be made by dissolving the 4SC-202 in the solvent of choice, which should be purged with an inert gas. 4SC-202 is soluble in the organic solvent DMSO.

Description

4SC-202 is an inhibitor of the class I histone deacetylases (HDACs) HDAC1-3 and the histone demethylase KDM1A.¹ It reduces cell viability in a panel of colorectal cancer (CRC) cell lines when used at concentrations ranging from 1 to 10 μ M and in patient-derived CRC cell lines when used at a concentration of 5 μ M.² 4SC-202 (5 μ M) induces cell cycle arrest at the G₂/M phase in HT-29 cells and primary human CRC cells and increases apoptosis in HT-29 cells in a concentration-dependent manner, an effect that is enhanced by the Akt inhibitors perifosine (Item No. 10008112) and MK-2206 (Item No. 11593). 4SC-202 (100 mg/kg, p.o., every other day) reduces tumor growth in an HT-29 mouse xenograft model when administered alone and to an enhanced degree when co-administered with oxaliplatin (Item No. 13106).

References

- 1. Maes, T., Carceller, E., Salas, J.A., et al. Advances in the development of histone lysine demethylase inhibitors. Curr. Opin. Pharmacol. 23, 52-60 (2015).
- 2. Zhijun, H., Shusheng, W., Han, M., et al. Pre-clinical characterization of 4SC-202, a novel class I HDAC inhibitor, against colorectal cancer cells. Tumour Biol. 37(8), 10257-10267 (2016).

WARNING THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFFTY 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.

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, 12/07/2022

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