

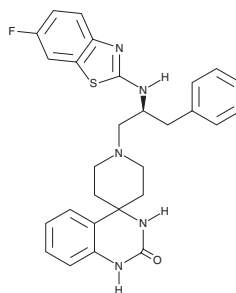
PRODUCT INFORMATION



NVS-ZP7-4

Item No. 33773

CAS Registry No.: 2349367-89-9
Formal Name: 1-[(2S)-2-[(6-fluoro-2-benzothiazolyl)amino]-3-phenylpropyl]-spiro[piperidine-4,4'-(1'H)-quinazolin]-2'-(3'H)-one
MF: C₂₈H₂₈FN₅OS
FW: 501.6
Purity: ≥98%
UV/Vis.: λ_{max}: 227, 256 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

NVS-ZP7-4 is supplied as a solid. A stock solution may be made by dissolving the NVS-ZP7-4 in the solvent of choice, which should be purged with an inert gas. NVS-ZP7-4 is soluble in ethanol and DMSO.

Description

NVS-ZP7-4 is an inhibitor of the zinc transporter SLC39A7 (ZIP7).¹ It induces apoptosis in TALL-1 human T cell acute lymphoblastic leukemia cells, an effect that can be reduced by expression of the ZIP7 point mutant ZIP7^{V430E}. NVS-ZP7-4 (20 μM) increases endoplasmic reticulum, but not cytosolic, zinc levels in U2OS cells. It inhibits ferroptosis induced by erastin (Item No. 17754) in MDA-MB-231 and RCC4 cells when used at a concentration of 10 μM.²

References

1. Nolin, E., Gans, S., Llamas, L., *et al.* Discovery of a ZIP7 inhibitor from a Notch pathway screen. *Nat. Chem. Biol.* **15**(2), 179-188 (2019).
2. Chen, P.-H., Wu, J., Xu, Y., *et al.* Zinc transporter ZIP7 is a novel determinant of ferroptosis. *Cell Death Dis.* **12**(2), 198 (2021).

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, 12/08/2022

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