

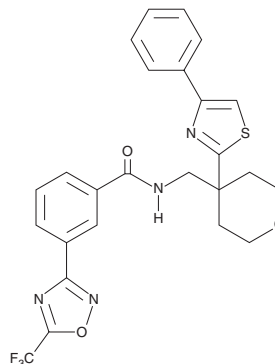
PRODUCT INFORMATION



TMP269

Item No. 17738

CAS Registry No.: 1314890-29-3
Formal Name: N-[[tetrahydro-4-(4-phenyl-2-thiazolyl)-2H-pyran-4-yl]methyl]-3-[5-(trifluoromethyl)-1,2,4-oxadiazol-3-yl]-benzamide
MF: C₂₅H₂₁F₃N₄O₃S
FW: 514.5
Purity: ≥95%
Supplied as: A crystalline 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

TMP269 is supplied as a crystalline solid. A stock solution may be made by dissolving the TMP269 in the solvent of choice, which should be purged with an inert gas. TMP269 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of TMP269 in these solvents is approximately 0.25, 10 and 25 mg/ml, respectively.

TMP269 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, TMP269 should first be dissolved in DMF and then diluted with the aqueous buffer of choice. TMP269 has a solubility of approximately 0.25 mg/ml in a 1:3 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

While class I HDACs are localized predominantly within the nucleus, class II HDACs shuttle into and out of the nucleus in response to intracellular signaling.¹ Class IIa HDACs, which include HDAC4, 5, 7, and 9, commonly act as corepressors and play diverse roles in cell biology.² TMP269 is a cell-permeable inhibitor of class IIa HDACs (IC₅₀s = 126, 80, 36, and 19 nM for HDAC4, 5, 7, and 9, respectively).³ It contains a metal-binding group, trifluoromethyloxadiazole that confers selectivity for class IIa HDACs.³

References

1. de Ruijter, A.J., van Gennip, A.H., Caron, H.N., *et al.* Histone deacetylases (HDACs): Characterization of the classical HDAC family. *Biochem. J.* **370**(Pt 3), 737-749 (2003).
2. Yang, X.-J. and Grégoire, S. Class II histone deacetylases: From sequence to function, regulation, and clinical implication. *Mol. Cell. Biol.* **25**(8), 2873-2884 (2005).
3. Lobera, M., Madauss, K.P., Pohlhaus, D.T., *et al.* Selective class IIa histone deacetylase inhibition via a nonchelating zinc-binding group. *Nat. Chem. Biol.* **9**, 319-325 (2013).

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