

PRODUCT INFORMATION



VER-246608

Item No. 19182

CAS Registry No.: 1684386-71-7
Formal Name: N-[4-(2-chloro-5-methyl-4-pyrimidinyl)phenyl]-N-[[4-[[[(2,2-difluoroacetyl)amino]methyl]phenyl]methyl]-2,4-dihydroxy-benzamide

MF: C₂₈H₂₃ClF₂N₄O₄

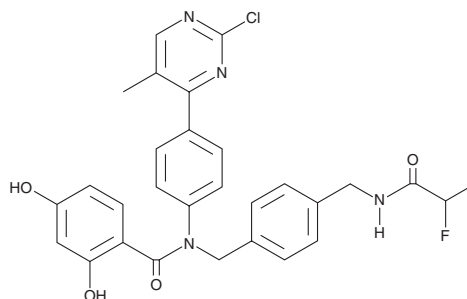
FW: 553.0

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

VER-246608 is supplied as a solid. A stock solution may be made by dissolving the VER-246608 in the solvent of choice, which should be purged with an inert gas. VER-246608 is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of VER-246608 in these solvents is approximately 2 mg/ml.

VER-246608 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, VER-246608 should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. VER-246608 has a solubility of approximately 0.5 mg/ml in a 1:1 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

VER-246608 is a pan-inhibitor of pyruvate dehydrogenase kinase (PDHK; IC₅₀s = 35, 84, 40, and 91 nM for PDHK1, -2, -3, and -4, respectively).¹ It is selective for PDHKs over heat shock protein 90 (Hsp90; IC₅₀ = >100 μM), as well as a panel of 97 kinases at 10 μM. VER-246608 (10 and 20 μM) decreases the production of L-lactate, a marker of glycolytic activity, in PC3 cells cultured in D-glucose- and L-glutamine-depleted media. It reduces the growth of, and induces cell cycle arrest at the G₁ phase in, serum-starved PC3 cells when used at a concentration of 20 μM.

Reference

1. Moore, J.D., Staniszewska, A., Shaw, T., et al. VER-246608, a novel pan-isoform ATP competitive inhibitor of pyruvate dehydrogenase kinase, disrupts Warburg metabolism and induces context-dependent cytostasis in cancer cells. *Oncotarget* **5(24)**, 12862-12876 (2014).

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, 12/22/2022

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