

PRODUCT INFORMATION

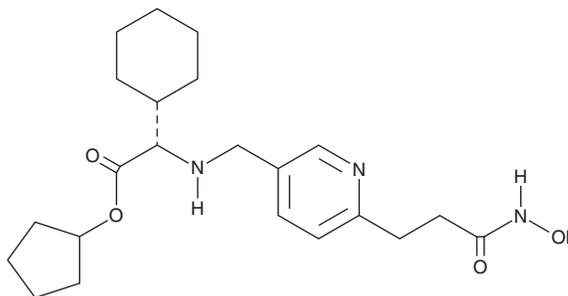


GSK3117391

Item No. 23846

CAS Registry No.: 1018673-42-1
Formal Name: αS-[[[6-[3-(hydroxyamino)-3-oxopropyl]-3-pyridinyl]methyl]amino]-cyclohexaneacetic acid, cyclopentyl ester

MF: C₂₂H₃₃N₃O₄
FW: 403.5
Purity: ≥98%
UV/Vis.: λ_{max}: 267 nm
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

GSK3117391 is supplied as a crystalline solid. A stock solution may be made by dissolving the GSK3117391 in the solvent of choice, which should be purged with an inert gas. GSK3117391 is soluble in the organic solvent DMSO.

Description

GSK3117391 is a histone deacetylase (HDAC) inhibitor (IC₅₀ = <100 nM in HeLa nuclear extracts).¹ It inhibits growth of U937 and HUT cancer cells (IC₅₀s = <330 nM).

Reference

1. Davidson, A.H., Moffat, D.F.C., Day, F.A., et al. HDAC Inhibitors *Chroma Therapeutics Ltd.* **WO 2008/040934 A1** (2008).

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, 12/17/2022

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