

PRODUCT INFORMATION



TCS HDAC6 20b

Item No. 14977

CAS Registry No.: 956154-63-5

Formal Name: 2-methyl-propanethioic acid, S-[(6S)-6-[[[(1,1-dimethylethoxy)carbonyl]amino]-7-oxo-7-(tricyclo[3.3.1.1^{3,7}]dec-1-ylamino)heptyl] ester

MF: C₂₆H₄₄N₂O₄S

FW: 480.7

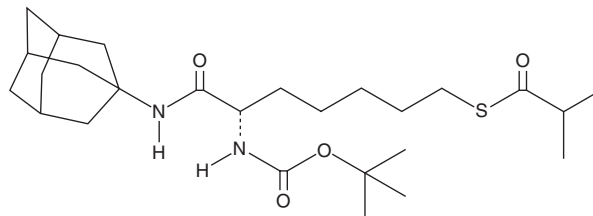
Purity: ≥95%

UV/Vis.: λ_{max}: 234 nm

Supplied as: A solution in ethanol

Storage: -20°C

Stability: ≥2 years



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

Laboratory Procedures

TCS HDAC6 20b is supplied as a solution in ethanol. To change the solvent, simply evaporate the ethanol under a gentle stream of nitrogen and immediately add the solvent of choice. Solvents such as DMSO purged with an inert gas can be used. The solubility of TCS HDAC6 20b in DMSO is approximately 20 mg/ml.

Description

TCS HDAC6 20b is an inhibitor of histone deacetylase 6 (HDAC6) that inhibits the growth of HCT116 cells synergistically when used at a concentration of 5 μM in combination with paclitaxel (Item No. 10461), but it has no activity on cell growth in HCT116 cells when used alone.¹ It inhibits cell growth of α-positive MCF-7 breast cancer cells treated with estrogen but not untreated cells.

Reference

1. Itoh, Y., Suzuki, T., Kouketsu, A., *et al.* Design, synthesis, structure--Selectivity relationship, and effect on human cancer cells of a novel series of histone deacetylase 6-selective inhibitors. *J. Med. Chem.* **50**(22), 5425-5438 (2007).

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, 11/04/2022

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