

PRODUCT INFORMATION



Tenovin-6 (hydrochloride)

Item No. 13086

CAS Registry No.: 1011301-29-3
Formal Name: N-[[[4-[[5-(dimethylamino)-1-oxopentyl]amino]phenyl]amino]thioxomethyl]-4-(1,1-dimethylethyl)-benzamide, monohydrochloride

MF: C₂₅H₃₄N₄O₂S • HCl

FW: 491.1

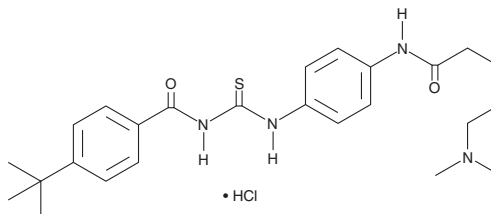
Purity: ≥95%

UV/Vis.: λ_{max}: 274 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

Tenovin-6 (hydrochloride) is supplied as a crystalline solid. A stock solution may be made by dissolving the tenovin-6 (hydrochloride) in the solvent of choice, which should be purged with an inert gas. Tenovin-6 (hydrochloride) is soluble in organic solvents such as ethanol and DMSO. The solubility of tenovin-6 (hydrochloride) in these solvents is approximately 1 and 20 mg/ml, respectively.

Description

The tumor suppressor function of p53 hinges on its ability to function as a transcription factor. Tenovin-6 is an analog of tenovin-1, a small molecule activator of p53 transcriptional activity. At 10 μM, this compound is slightly more effective than tenovin-1 at elevating p53 activity in MCF-7 cells and reduces growth of ARN8 melanoma xenograft tumors in SCID mice at a dose of 50 mg/kg.¹ Tenovin-6 inhibits the protein deacetylase activities of purified human SIRT1, SIRT2, and SIRT3 *in vitro* with IC₅₀ values of 21, 10, and 67 μM, respectively.¹

Reference

1. Lain, S., Hollick, J.J., Campbell, J., *et al.* Discovery, *in vivo* activity, and mechanism of action of a small-molecule p53 activator. *Cancer Cell* **13**(5), 454-463 (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, 11/08/2022

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