

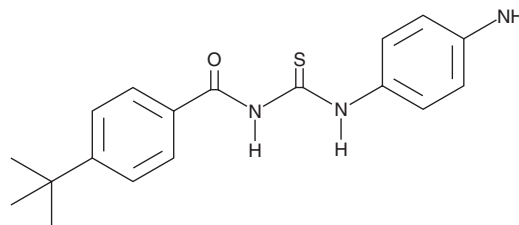
PRODUCT INFORMATION



Tenovin-3

Item No. 16288

CAS Registry No.: 1011301-27-1
Formal Name: N-[[[4-(aminophenyl)amino]thioxomethyl]-4-(1,1-dimethylethyl)-benzamide
MF: C₁₈H₂₁N₃OS
FW: 327.4
Purity: ≥98%
UV/Vis.: λ_{max}: 203, 253, 277 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-3 is supplied as a crystalline solid. A stock solution may be made by dissolving the tenovin-3 in the solvent of choice, which should be purged with an inert gas. Tenovin-3 is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of tenovin-3 in these solvents is approximately 0.25 and 0.5 mg/ml, respectively.

Description

Tenovin-3 is a small molecule activator of p53 transcriptional activity.¹ It increases p53 activity in MCF-7 cells when used at a concentration of 10 μM. It also increases the level of K40-acetylated α-tubulin in H1299 cells and inhibits SIRT2 protein deacetylase activity.

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, 10/26/2022

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