

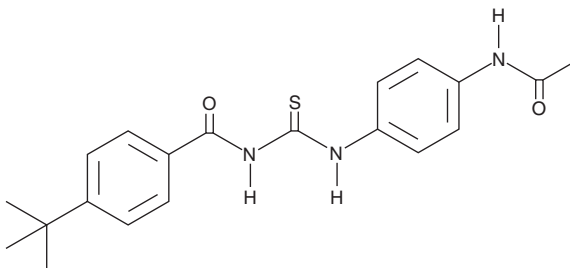
PRODUCT INFORMATION



Tenovin-1

Item No. 13085

CAS Registry No.: 380315-80-0
Formal Name: N-[[[4-(acetylamino)phenyl]amino]thioxomethyl-4-(1,1-dimethylethyl)]-benzamide
MF: C₂₀H₂₃N₃O₂S
FW: 369.5
Purity: ≥98%
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-1 is supplied as a crystalline solid. A stock solution may be made by dissolving the tenovin-1 in the solvent of choice, which should be purged with an inert gas. Tenovin-1 is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of tenovin-1 in these solvents is approximately 0.25 and 0.50 mg/ml, respectively.

Description

Tenovin-1 is a small molecule activator of p53 transcriptional activity. At 10 μM, it elevates p53 expression in MCF-7 cells within two hours of treatment and longer-term exposure significantly decreases the growth of BL2 Burkitt's lymphoma and ARN8 melanoma cells.¹ Functioning upstream of p53, tenovin-1 acts by inhibiting the deacetylase activity of purified human SIRT1 and SIRT2, members of NAD⁺-dependent class III histone deacetylases that belong to the sirtuin family.¹ While tenovin-1 demonstrates low genotoxicity, the compound has poor water solubility, which limits its uses *in vivo*.

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, 02/28/2024

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