

# Tenovin 1

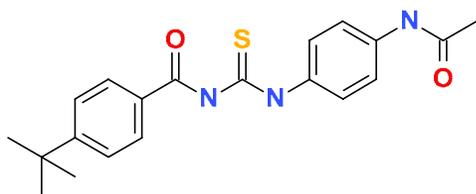
Cat. # SI9865

**Background:** Tenovin 1 inhibits the deacetylase activity of SirT1 and SirT2. One of the substrates for SirT1 is p53 which is stabilized by acetylation of Lys-382. The ability of p53 to bind DNA is also enhanced by acetylation of Lys-382. Hence, inhibition of SirT1 by tenovin 1 leads to increased stability and activity of p53 (Lain, Hollick et al. 2008).

**Application:** Studying the role of sirtuins in regulating cell growth

## Product Information

<b>CAS No.:</b>	380315-80-0
<b>Purity:</b>	99%; NMR consistent with structure
<b>Molecular Weight:</b>	369.48
<b>Physical State:</b>	Powder
<b>Quantity:</b>	10mg
<b>Solubility:</b>	200mM in DMSO
<b>Storage:</b>	Store dessicated as supplied at -20°C for 2 years
<b>Formula:</b>	C <sub>20</sub> H <sub>23</sub> N <sub>3</sub> O <sub>2</sub> S
<b>SMILES String:</b>	CC(=O)Nc1ccc(cc1)NC(=S)NC(=O)c2ccc(cc2)C(C)(C)C



## References

Lain, S., J. J. Hollick, et al. (2008). "Discovery, in vivo activity, and mechanism of action of a small-molecule p53 activator." *Cancer Cell* **13**(5): 454-463.

All products are for research use only • not intended for human or animal diagnostic or therapeutic uses  
Copyright © 2013 LifeSensors, Inc. All Rights Reserved