

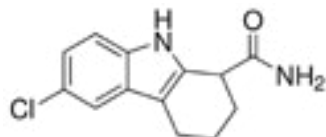
# EX-527

Catalog Number P005-5MG

Catalog Number P005-25MG

## FEATURES

- Selective SIRT1 inhibitor
- Does not inhibit HDACs
- Inhibitor of SIRT1 deacetylation of p53



ARBOR  
ASSAYS

## INTRODUCTION

EX-527 is a cell-permeable, selective inhibitor of SIRT1 (IC<sub>50</sub> = 98 nM *in vivo*, 38nM *in vitro*) that does not inhibit histone deacetylase (HDAC) or other sirtuin deacetylase family members (IC<sub>50</sub> values are 20 μM, 49 μM, > 100 μM and > 100 μM for SIRT2, SIRT3, HDAC and NADase respectively). EX-527 has been used to investigate the relationship between SIRT1-mediated deacetylation of p53, p53 activity, and cell survival following DNA damage.

<b>FORM:</b>	Light yellow powder
<b>MOLECULAR WEIGHT:</b>	284.71
<b>STORAGE:</b>	4°C, desiccated
<b>FORMULA:</b>	C <sub>13</sub> H <sub>13</sub> ClN <sub>2</sub> O
<b>CAS NUMBER:</b>	49843-98-3
<b>OTHER NAMES:</b>	6-Chloro-2,3,4,9-tetrahydro-1H-carbazole-1-carboxamide, racemate
<b>USES:</b>	Soluble to 18 mg/mL in DMSO and 10 mg/mL in Ethanol

## REFERENCES:

Napper, AD, et al. Discovery of indoles as potent and selective inhibitors of the deacetylase SIRT1. J.Med.Chem. 48:8045. (2005)

Solomon, JM, et al. Inhibition of SIRT1 catalytic activity increases p53 acetylation but does not alter cell survival following DNA damage. Mol.Cell.Biol. 26:28. (2006)

Nayagam, V.M., Wang, X., Tan, Y.C., et al. SIRT1 modulating compounds from high-throughput screening as anti-inflammatory and insulin-sensitizing agents. J Biomol Screen 11:959-967 (2006)

**FOR RESEARCH USE ONLY**

info@ArborAssays.com • www.ArborAssays.com

10Aug12