

Catalog No: 14028, 14029

Format: 25 mg, 5 mg

Chemical Properties:

MW = 248.7 $C_{13}H_{13}CIN_2O$

CAS 49843-98-3

Physical Properties: Light yellow powder

Names: EX-527, 6-Chloro-2,3,4,9-tetrahydro-1H-carbazole-1-carboxamide

(racemic)

Pharmacology: Selective SIRT1 inhibitor (IC₅₀ = 98 nM). Does not inhibit other HDACs or SIRT family members. Increases p53 acetylation following DNA damage (ref 1). Causes cell cycle arrest at G1 (ref 2). Inhibits differentiation and facilitates expansion of hematopoietic progenitor cells (ref 3). An extremely useful tool for determining the involvement of SIRT1 in cellular physiology (ref 4). Cell permeable.

Solubilization: May be dissolved in DMSO (18 mg/ml); Ethanol (10 mg/ml)

Fluorescent Properties: N/A

Quality Control:

>98% (TLC); NMR (Conforms)

References:

1. JM Solomon et al. Mol. Cell Biol. 2006, 26:28

2. B Peck et al. Mol. Cancer Ther. 2010, 9:844

3. T Peled et al. Exp. Hematol.. 2012, 40:342

4. JL Andersen et al. Mol. Cell 2011, 43:834

Storage and Guarantee: Store desiccated as supplied at 4°C for up to 2 years. Store solutions at -20°C for up to 2 months. This product is guaranteed for 6 months from date of arrival.

