## SirtAct



Catalog No: 14080, 14081

Format: 25 mg, 5 mg

**Chemical Properties:** 

MW = 367.4 $C_{20}H_{25}N_5O_2$ 

CAS 839699-72-8

Physical Properties: Olive green powder

Names: SirtAct, Sirt1 activator 3, CAY10591, 2-Amino-N-cyclopentyl-1-(3-

methoxypropyl)-1H-pyrrolo[2,3-b]quinoxaline-3-carboxamide

**Pharmacology:** Sirt1 activator. It suppressed TNFα release from THP-1 cells and stimulated adipogenesis in 3T3L1 cells (ref 1). SIRT1 stimulation by SirtAct leads to decreased tissue factor mediated thrombogenicity in mice (ref 2). It inhibited NFκB activation and reduced inflammatory cytokine production in lamina propria mononuclear cells from IBD patients (ref 3). It prevented and cured experimental Colitis in a mouse model (ref 3).

Solubilization: May be dissolved in DMSO (50 mg/ml); Ethanol (20 mg/ml,

warm)

Fluorescent Properties: N/A

**Quality Control:** 

>98% (TLC); NMR (Conforms)

References:

1. VM Nayagam et al. J. Biomol. Screen. 2006, 11:959

2. SS Barbieri et al. Circulation 2012, 126:1373

3. R Caruso et al. Mucosal Immunol. 2014, May 21 Epub

**Storage and Guarantee:** Store desiccated as supplied at -20°C for up to 2 years. Store solutions at -20°C for up to 3 months. This product is guaranteed for 6 months from date of arrival.