

Salermide

Catalog No: 14124, 14125

Format: 25 mg, 5 mg

Chemical Properties:

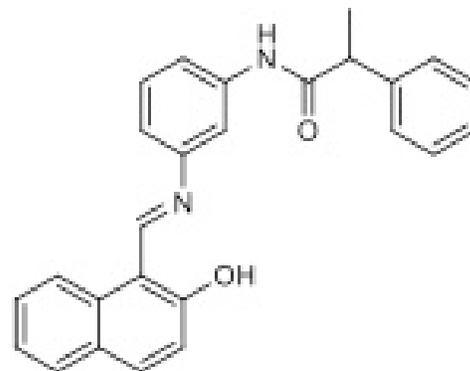
MW = 394.5

C₂₆H₂₂N₂O₂

CAS 1105698-15-4

Physical Properties: Yellow powder; MP = 175-177°C

Names: Salermide, N-[3-[[[2-Hydroxy-1-naphthalenyl)methylene]amino]phenyl]-α-methylbenzeneacetamide



Pharmacology: SIRT1 and SIRT2 inhibitor. Exhibits a stronger inhibitory effect on SIRT2 than on SIRT1 *in vitro*. Induces the reactivation of proapoptotic genes repressed by SIRT1 and causes massive apoptosis in cancer cells within 24 hours (ref 1). Displays protective effect in nematodes expressing muscular dystrophy protein (ref 2). Promotes neuronal apoptosis confirming the protective effect of SIRT1 (ref 3).

Solubilization: May be dissolved in DMSO (40 mg/ml); ethanol (10 mg/ml; warm); or DMF (30 mg/ml)

Fluorescent Properties: N/A

Quality Control:

>98% (TLC); NMR (Conforms)

References:

1. E Lara *et al. Oncogene* 2009, 28:781
2. MY Pasco *et al. J. Med. Chem.* 2010, 53:1407
3. Y Zhao *et al. Exp. Neurol.* 2012, 237:489

Storage and Guarantee: Store desiccated as supplied at -20° for up to 2 years. Store solutions at -20°C for up to 1 month. This product is for research use only and is not for use in diagnostic procedures. This product is guaranteed for 6 months from date of arrival.